

# AlphaScreen & AlphaLISA

化学放大型 亲和发光  
均相检测

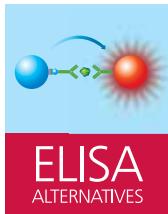
AlphaLISA INSTRUCTIONS  
FOR BEST RESULTS,  
DO NOT WASH

- REDUCE ASSAY TIME  
TO TWO HOURS
- WILL NOT WASH AWAY  
WEAK INTERACTIONS
- LEAVE OVERNIGHT  
IF NEEDED
- USE CONFIDENTLY RIGHT  
OUT OF THE BOX



从小分子到活细胞  
适于各种样品的  
检 测 系 统

PerkinElmer®



[www.perkinelmer.com](http://www.perkinelmer.com)

PerkinElmer®  
*For the Better*

# Alpha 技术：仅需混合操作的简单检测

Alpha (Amplified: 放大、Luminescence: 化学发光、Proximity: 接近、Homogeneous: 均相、Assay: 检测) 是基于微珠的高灵敏度均相 (Mix & Measure) 检测技术。

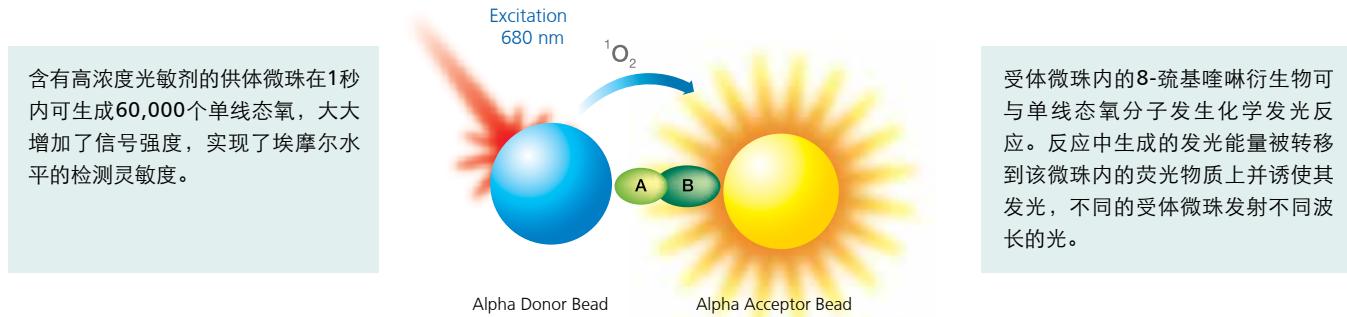
Alpha 技术可快速检测 (无需洗涤操作) 种类多样的各生物学分子间的相互作用。Alpha 是 Dade Behring 公司 (现为 Siemens AG 公司) 基于 LOCI (Luminescent Oxygen Channeling Immunoassay) 技术\*开发而成。PerkinElmer 在获得 Alpha 技术的独家销售权的情况下，针对生命科学研究领域生产和销售微珠和相关试剂。

\*Ullman EF, et al. "Luminescent oxygen channeling Immunoassay: measurement of particle binding kinetics by chemiluminescence." Proc Natl Acad Sci USA. 1994 Jun 7;91(12):5426-30.

## 原理

Alpha 技术使用供体和受体两种微珠。结合于供体微珠上的分子 (分子 A) 与结合于受体微珠上的分子 (分子 B) 发生相互作用，当两个微珠达到足够接近的状态时，能够检测出发光信号。

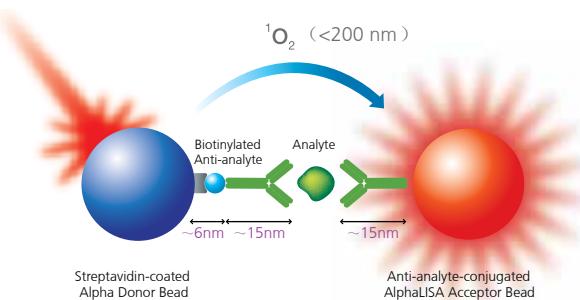
供体微珠表面的光敏剂经激光照射而受到激发，将周围的氧转变为激发态的单线态氧。单线态氧向供体微珠周围扩散，到达附近的受体微珠后可在微珠内引发化学发光反应，并最终发射出光。如果各微珠上的分子 A 与分子 B 不发生相互作用，则供体微珠产生的单线态氧不会到达受体微珠，从而不会引发化学发光反应。



## 特点

**均相 (Mix & Measure)**..... 无需分离和洗涤步骤，操作简便。

**非常适合生物分子的相互作用** ..... 不同于 FRET 检测，单线态氧的扩散距离可达到 200 nm，最大程度减少了供体/受体之间的距离和方向限制。因此可用于全长蛋白检测和使用抗体的夹心免疫测定等应用 (下图)。



**具有较宽范围的亲和性**..... 可检测亚 nM ~ μM (低亲和性) 范围的相互作用。

**低背景、高灵敏度**..... 通过激发光照射进行检测的方法会受到样品和微孔板生成的自发荧光的影响，产生较高的背景。Alpha 技术中，与单线态氧发生反应的受体微珠的发光时间较长 (半衰期: 300 毫秒)，因此可在激发光照射完一段时间后 (20 毫秒后) 再开始检测，从而有效避免自发荧光的影响。此外，检测波长比激发波长更短，而自发荧光的波长比激发波长更长，因此自发荧光的影响也会进一步降低。

**可实现自动化**..... 由于微珠的体积非常小 (直径 250 nm) 且不易发生沉淀，因此不会堵塞分液头，易于实现自动化分液操作，非常适合 HTS。

**可检测微量样品**..... 可检测低至 5 μL 的样品。不会改变试剂浓度或造成灵敏度下降，适用于常规的 96、384 和 1536 孔规格。贴上 TopSeal-A (粘附密封膜) 的情况下进行检测可有效避免样品蒸发。

## Alpha 微珠 (供体: 1 种; 受体: 4 种)

Alpha 微珠的直径很小, 约为 250 nm。这种直径远远小于 FMAT ( $6 \sim 20 \mu\text{m}$ ) 等微珠检测中所用微珠的直径。因此 Alpha 微珠可在水溶液中保持稳定并均匀分散成悬浊液, 经过数小时也不会发生沉淀。微珠采用聚苯乙烯制成, 并涂覆有葡聚糖聚合物以避免非特异性结合。各类生物分子通过葡聚糖上的醛基结合到微珠表面。

\*\* 1 个微珠表面涂覆有约 3,000 个链霉亲合素分子或 300 个 IgG 分子。

### 检测波长

受体微珠有 4 种含有不同荧光染料的类型可供选择。各种类型的检测波长不同, 因此可用于多靶点检测或根据样品种类来选择使用。专为使用血清/血浆的检测而开发的 AlphaLISA 受体微珠所在的波长区域可有效避免血清/血浆样品的荧光背景干扰 (300-600 nm) 以及氧合血红蛋白的吸收干扰 (<600 nm)。此外, 该款微珠的检测波长区域比 AlphaScreen 受体微珠更为狭窄, 可获得更强的光信号。

#### • AlphaScreen (红荧烯; 发射波长: 520-620 nm) :

适用于缓冲液、细胞培养上清液和细胞裂解液样品。

#### • AlphaLISA (铕; 发射波长: 615-623 nm) :

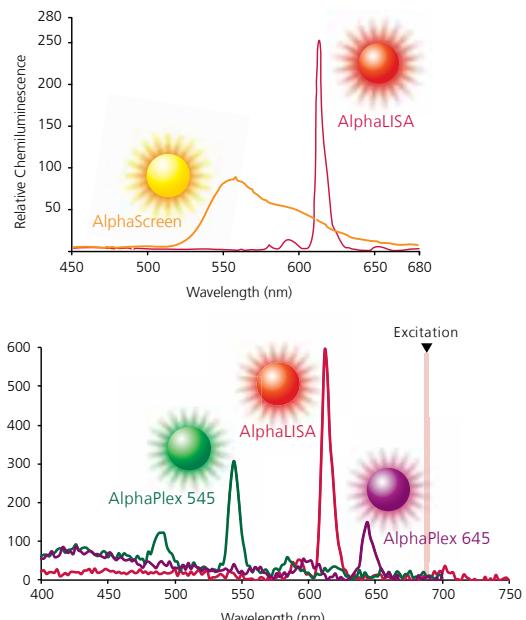
适用于缓冲液、细胞培养上清液和细胞裂解液样品。

#### • AlphaPlex 545 (铽; 发射波长: 515-555 nm) :

适用于缓冲液、细胞培养上清液和细胞裂解液样品。

#### • AlphaPlex 645 (钐; 发射波长: 645 nm) : 新品

适用于缓冲液、细胞培养上清液和细胞裂解液样品。



## 多靶点检测

通过联用 AlphaLISA 微珠和 AlphaPlex 微珠, 用户可同时检测多个目标, 即进行多靶点检测。

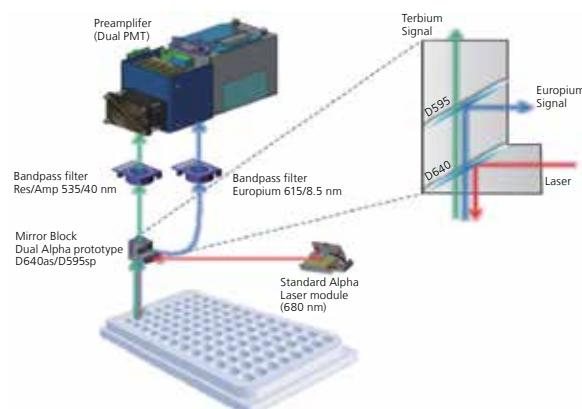
### 多靶点检测中的必要设备和附件

附件名称	产品货号
设备主机	EnVision Xcite / HTS 敬请询问
选件*1	Alpha 标准选件 2102-8340
	现场升级包 2108-8330
软件*2	EnVision 软件升级 2101-8020
	EnVision PC 升级包 2104-8040
连续检测多个波长 (各发光波长依次被检测)	
滤光器	Europium 2100-5090
	AlphaPlex Tb 2100-5930
	AlphaPlex Sm 2100-5940
单镜	AlphaScreen D640as (Eu-Tb) *3 2101-4010
	AlphaPlex D670as (Eu-Tb-Sm) 2102-5910
同时检测 2 个波长 (仅配备 2PMT 时)	
滤光器	见上述 —
双镜	Dual Alpha Eu/Tb 2102-5900
	Dual Alpha Eu/Sm or Tb/Sm 2102-5920

### 通过 Alpha 双镜模块实现同时检测

为 EnVision 配备第二检测器和 Alpha 双镜模块后, 就可以同时检测 AlphaLISA 微珠 (615 nm) 和 AlphaPlex 545 微珠 (545 nm) 的发射光信号。

没有配备第二检测器的情况下, 两种发射光信号将顺次被检出。



\*1 Alpha HTS 选件不能用于多靶点检测。

\*2 必须安装 EnVision 软件 1.13 版。同时需要升级 PC。

\*3 Alpha Standard 选件中的标配。

# Alpha 技术：仅需混合操作的简单检测

## Alpha 常见问答

**问题：** Alpha 检测采用何种设备进行检测？

我可以使用现有的荧光微孔板检测仪或光度计进行检测吗？

**回答：** Alpha 检测需要使用专用激光来发射 680 nm 的激发光。基于氙闪光灯的荧光微孔板检测仪或光度计不能用于Alpha检测。

AlphaScreen与AlphaLISA可使用PerkinElmer的EnVision (Alpha)、EnSpire (Alpha)、EnSight (Alpha)系统来检测。由于Victor未配备内置激光，因此不能用于Alpha检测。Alpha检测最适合使用 PerkinElmer 的检测设备。

AlphaPlex只能使用EnVision (Alpha选件) 来检测，并且还需要其他滤光器和光学模块。(关于AlphaPlex检测中所需附件的详细信息，请参见第 3 页)。



EnVision



EnSpire



EnSight

**问题：** 需要使用专用微孔板吗？

**回答：** 推荐使用颜色介于白色微孔板和灰色微孔板之间的灰白色 AlphaPlate (第 38 页)。Alpha 技术中使用发射光检测器来检测信号。检测发射光时，大多数情况使用白色微孔板，但使用 AlphaPlate 可抑制交叉干扰并提高检测精度。

**问题：** 听说 Alpha 不能承受较强的光，这是真的吗？

**回答：** 受体微珠对光不敏感，但链霉亲合素供体微珠对光较为敏感。在阳光直射或强光照射下，供体微珠受到激发并生成单线态氧，这类单线态氧会与供体微珠上结合的链霉亲合素发生反应，降低其效率。如果不注意的情况下使用过量光照射，就会造成信号强度降低，但大多数情况下，该检测都能获得良好的结果。最好将供体微珠相关步骤放在100Lux以下（右表）、限制了光强度的实验室中进行。例如，可关闭实验室中一半的照明灯，在关灯的一侧进行实验操作等。也可将绿色滤光器贴在照明灯上。Alpha检测中进行孵育时，请用不透明的微孔板或铝箔覆盖检测用微孔板，以达到遮光目的。

Lux 水平	
700- Lux	窗边
320-500 Lux	室内 (照明灯正下方)
30-70 Lux	室内 (照明灯非正下方/背光处)
10 Lux	桌下

**问题：** 缓冲液和化合物等会对 Alpha 造成影响吗？

**回答：** 单线态氧的清除剂 NaN3 (0.005% 以上)、μM 级的过渡金属 ( $\text{Fe}^{2+}$ 、 $\text{Fe}^{3+}$ 、 $\text{Cu}^{2+}$ 、 $\text{Ni}^{2+}$ 、 $\text{Al}^{2+}$ )、抗坏血酸以及基本结构为噻吩的化合物 (100 μM ~ 10 mM) 会对检测造成影响。虽然超过上述浓度时会造成信号降低，但大多数情况下仍可获得可接受的结果。一般使用的 HEPES、PBS、Tris 等缓冲液的浓度和 pH 值基本上不会影响检测。盐浓度在 300 mM 以下以及非离子表面活性剂的用量在 3% 以下时不会影响检测。此外，最高可耐受 5% DMSO、0.02% SDS 或 100 mM EDTA。

**问题：** 微珠能够与我手上现有的抗体或带标签的蛋白结合吗？

**回答：** 可以。我们提供未标记的微珠、标记有用于捕获二抗或表达标记的谷胱甘肽或镍螯合物等物质的供体微珠和受体微珠 (第 31 ~ 41 页)。并提供标记定制服务。

**问题：** 关于微珠的分液操作，有什么特别的技巧吗？

**回答：** 缓冲液中经常含有表面活性剂或 BSA 等物质，因此推荐使用反向移液法。由于是从顶部进行检测，因此如果有气泡的话会影响检测信号。采用移液操作的话不需要混合或搅拌。

# AlphaLISA：只需混合无需洗涤的ELISA

## AlphaLISA 无需洗涤的ELISA

AlphaLISA技术帮助您进行血清，血浆及细胞裂解液等复杂样品中靶蛋白的检测。无论小分子激素蛋白或大分子抗体蛋白均能通过简单加样无需洗涤的Alpha技术来检测

### AlphaLISA 免疫检测试剂盒

适用于细胞裂解液，血清及血浆样本中靶蛋白的检测。您只需准备样品，即刻就能进行检测。

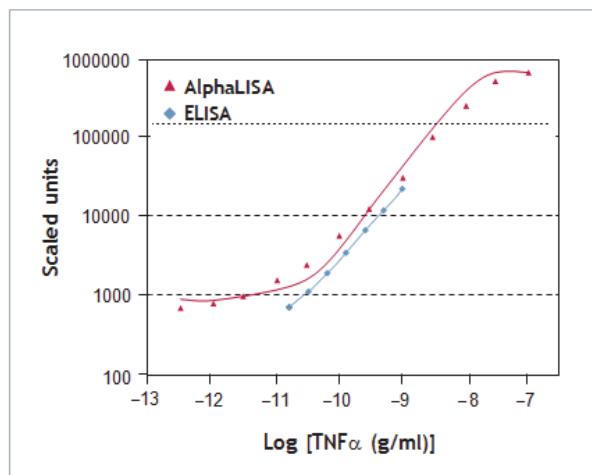
#### AlphaLISA vs ELISA

AlphaLISA 因其均相免洗的特性，可以采用比ELISA更小体积的检测体系。

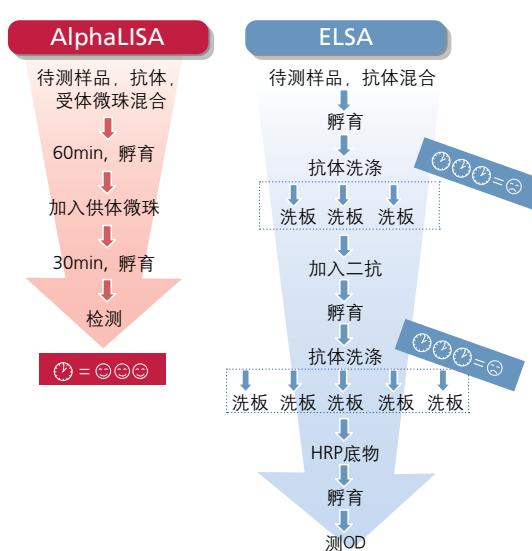
如果与自动加样系统或其他自动化系统整合，可以进行高通量筛选实验。

	AlphaLISA	ELISA
洗板	不要	必要
实验步骤	3 ~ 4 步	> 5 步
微孔板型号	96, 384, 1536-well	96-well
自动化程度	★★★	★
实验通量	高	低
样品需要量	少 (5 μL 以下)	多
灵敏度	★★★★	★★★
动态检测范围	2.5-5 logs	2 logs
性价比	★★★	★

#### 动态检测范围比



#### 实验步骤比



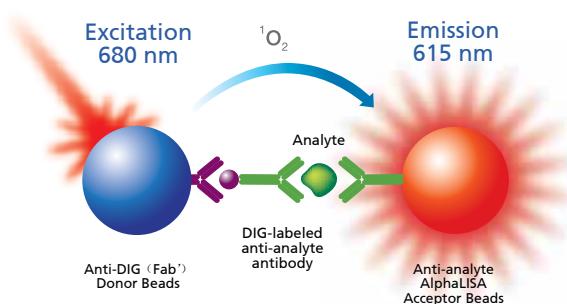
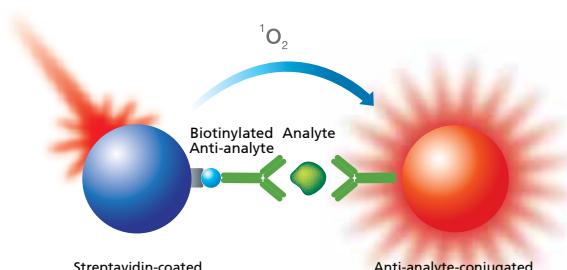
# AlphaLISA：只需混合无需洗涤的ELISA

## 免疫分析试剂盒原理

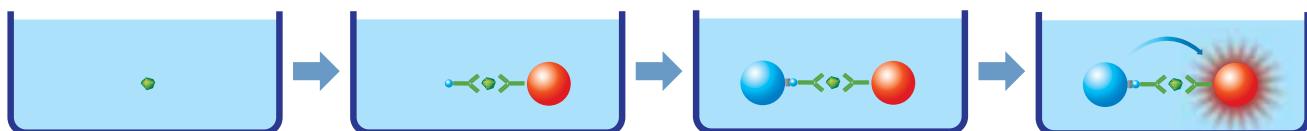
细胞裂解液、血清，或血浆样品中的待测物、分别被生物素化抗体和受体微珠偶联的抗体识别，形成三明治形式。之后加入链霉亲和素包被的供体微珠结合生物素化抗体。

## Biotin-free (DIG-based) 免疫分析试剂盒 新品

与常规使用的免疫分析试剂盒不同的是，前者用链霉亲和素包被的供体微珠识别生物素化抗体，该试剂盒用抗体高辛抗体包被的供体微珠识别地高辛标记的抗体。该试剂盒更适用于RPMI培养液的细胞学实验，能够避免培养液中的生物素干扰，得到更高的灵敏度和S/B值。



## 实验步骤



将5 $\mu\text{L}$ 样品（待测物）加入微孔板。

加入生物素化抗体，及与受体微珠偶联的抗体，共计20  $\mu\text{L}$ ，室温孵育60分钟。

加入25 $\mu\text{L}$ 链霉亲和素包被的供体微珠，室温孵育30分钟（总体积为50  $\mu\text{L}$ ）。

激光激发供体微珠，产生的单线态氧扩散至受体微珠附近，受体微珠接受单线态氧后发光。

The screenshot shows a search interface for Certificates of Analysis (COAs). It includes fields for 'Product ID' and 'Search term', and a 'Search' button. Below the search bar, there's a list of results with columns for 'Product ID', 'Product Name', and 'COA Number'. One result is highlighted: 'N08000 - AlphaLISA® HRP-CP Kit' with COA number 'COA-N08000-001'. The page also includes links for 'Product of Interest' and 'Certifications'.



### 试剂盒组分

- 抗体偶联的受体微珠
- 链霉亲和素包被的供体微珠
- 生物素化抗体
- 反应溶液, 10X
- 标准品冻干粉

出于环保考虑，产品未附带纸质版使用说明，您可以进入以下网址  
<http://www.perkinelmer.com/tools/COASearch>

输入产品货号或批号进行查询下载



产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
免疫炎性靶点							
Human CRP	500	AL233C	7.3pg/mL	1,000,000pg/mL	✓	X	X
	5000	AL233F					
	3μg	AL233S					
Human CXCL1/GRO-α	100	AL349HV					
	500	AL349C	0.5pg/mL	30,000pg/mL	✓	X	X
	5000	AL349F					
	0.1μg	AL349S					
Human CXCL10/IP-10	500	AL259C	8.6pg/mL	100,000pg/mL	✓	X	X
	5000	AL259F					
	0.3μg	AL259S					
	500	AL280C	1.6pg/mL	100,000pg/mL	✓	X	
Human CXCL9/MIG	5000	AL280F					
	0.3μg	AL280S					
	500	AL213C	14pg/mL	300,000pg/mL	✓	X	
Human G-CSF	5000	AL213F					
	1μg	AL213S					
	500	AL216C	1.6pg/mL	10,000pg/mL	✓	X	X
Human GM-CSF	5000	AL216F					
	0.03μg	AL216S					
	500	AL517C	2pg/mL	10,000pg/mL	X	✓	X
Mouse GM-CSF	5000	AL517F					
	10ng	AL517S					
	500	AL297C	8.2pg/mL	100,000pg/mL	✓	X	X
Human IFN-α	5000	AL297F					
	0.3μg	AL297S					
	500	AL217C	11pg/mL	100,000pg/mL	✓	X	X
Human IFN-γ	5000	AL217F					
	0.3μg	AL217S					
	100	AL327HV					
Human IFN-γ biotin-free	500	AL327C	2.4pg/mL	10,000pg/mL	✓	X	X
	5000	AL327F					
	0.3μg	AL327S					
Mouse IFN-γ	500	AL501C	1.1pg/mL	100,000pg/mL	X	✓	X
	5000	AL501F					
	0.1μg	AL501S					
Human IL-1α	500	AL238C	1.3pg/mL	30,000pg/mL	✓	X	X
	5000	AL238F					
	0.1μg	AL238S					
Human IL1β	500	AL220C	0.6pg/mL	30,000pg/mL	✓	X	X
	5000	AL220F					
	0.1μg	AL220S					
Mouse IL1β	500	AL503C	11.5pg/mL	100,000pg/mL	X	✓	X
	5000	AL503F					
	0.1μg	AL503S					
Rat IL1β	500	AL518C	6.2pg/mL	100,000pg/mL	X	X	✓
	5000	AL518F					
	0.3μg	AL518S					
Human IL1RA	100	AL358HV					
	500	AL358C	0.43pg/mL	10,000pg/mL	✓		
	5000	AL358F					
Human Vimentin	0.1μg	AL358S					
	100	AL372HV	Coming soon				
	500	AL372C					
Human COX2	5000	AL372F					
	1μg	AL372S					
	100	AL376HV	Coming soon				
Human Claudin 7	500	AL376C					
	5000	AL376F					
	1μg	AL376S					
Human BCL2	100	AL381HV	Coming soon				
	500	AL381C					
	5000	AL381F					
	1μg	AL381S					
其他							
Human IL2	500	AL221C					
	5000	AL221F					
	0.3μg	AL221S					
Human IL2	100	AL333HV					
	500	AL333C	2.6pg/mL	30,000pg/mL	✓	X	X
	5000	AL333F					
Mouse IL2	0.3μg	AL333S					
	500	AL511C	1.8pg/mL	30,000pg/mL	X	✓	X
	5000	AL511F					
Human IL3	0.1μg	AL511S					
	500	AL222C	11pg/mL	300,000pg/mL	✓	X	X
	5000	AL222F					
Human IL5	0.3μg	AL222S					
	500	AL267C	4.5pg/mL	100,000pg/mL	✓	X	X
	5000	AL267F					
Human IL6	0.3μg	AL267S					
	500	AL223C	1.3pg/mL	30,000pg/mL	✓	X	X
	5000	AL223F					
Mouse IL6	0.1μg	AL223S					
	500	AL504C	1.4pg/mL	100,000pg/mL	X	✓	✓
	5000	AL504F					
Human IL7	0.3μg	AL504S					
	500	AL268C	8.5pg/mL	100,000pg/mL	✓	X	X
	5000	AL268F					
Human IL8	0.3μg	AL268S					
	500	AL512C	24.8pg/mL	300,000pg/mL	X	✓	✓
	5000	AL512F					
Mouse IL7	0.3μg	AL512S					
	500	AL224C	1.1pg/mL	30,000pg/mL	✓		
	5000	AL224F					
Human IL8 Biotin-free	0.1μg	AL224S					
	100	AL328HV					
	500	AL328C	3.6pg/mL	30,000pg/mL	✓		
Human IL8	5000	AL328F					
	0.1μg	AL328S					
	500	AL218C	4.4pg/mL	30,000pg/mL	✓	X	X
Human IL10	5000	AL218F					
	0.3μg	AL218S					
	500	AL502C	13.5pg/mL	300,000pg/mL	X	✓	X
Mouse IL10	5000	AL502F					
	0.3μg	AL502S					
	500	AL293C	8.6pg/mL	300,000pg/mL	✓		
Human IL11	5000	AL293F					
	0.3μg	AL293S					
	500	AL239C	4.9pg/mL	300,000pg/mL	✓	X	X
Human IL12(P70)	5000	AL239F					
	0.3μg	AL239S					
	500	AL240C	3.9pg/mL	100,000pg/mL	✓	X	X
Human IL13	5000	AL240F					
	0.1μg	AL240S					
	100	AL382HV					
Human IL-12 p40/IL-23 p40	500	AL382C	0.45pg/mL	100,000pg/mL	✓		
	5000	AL382F					
	1μg	AL382S					
Alpha Tubulin	100	AL383HV					
	500	AL383C	77.1pg/mL	1,000,000pg/mL	✓		
	5000	AL383F					
Human BCL2	1μg	AL383S					
	100	AL384HV					
	500	AL384C	30pg/mL	1,000,000pg/mL	✓	X	
Human BCL2	5000	AL384F					
	1μg	AL384S					

# AlphaLISA：只需混合无需洗涤的ELISA

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
Human IL15	500	AL283C	1.2pg/mL	100,000pg/mL	✓	X	X
	5000	AL283F					
	0.1µg	AL283S					
Human IL15 Biotin-Free	100	AL360HV					
	500	AL360C	1.9pg/mL	100,000pg/mL	✓		
	5000	AL360F					
Mouse IL15	0.1µg	AL360S					
	500	AL513C	20.7pg/mL	300,000pg/mL	X	✓	✓
	5000	AL513F					
Human IL17	0.3µg	AL513S					
	500	AL219C	14pg/mL	300,000pg/mL	✓	X	X
	5000	AL219F					
Human IL-17A Biotin-Free	1µg	AL219S					
	100	AL346HV					
	500	AL346C	13.5pg/mL	30,000pg/mL	✓		
Mouse IL17A	5000	AL346F					
	0.3µg	AL346S					
	500	AL514C	4.1pg/mL	100,000pg/mL	X	✓	✓
Mouse Interleukin -17A Homodimer	5000	AL514F					
	0.1µg	AL514S					
	100	AL540HV					
Human IL18	500	AL540C	3.5pg/mL	300,000pg/mL	X	✓	X
	5000	AL540F					
	0.3µg	AL540S					
Human/Rat IL22	500	AL241C	5pg/mL	300,000pg/mL	✓	X	X
	5000	AL241F					
	0.3µg	AL241S					
Human IL28B	100	AL320HV					
	500	AL320C	18pg/mL	1,000,000pg/mL	✓	X	✓
	5000	AL320F					
Human IP10	1µg	AL320S					
	100	AL365HV					
	500	AL365C	16pg/mL	100,000pg/mL	✓	X	
Human TNFa	5000	AL365F					
	0.3µg	AL365S					
	100	AL326HV					
Human TNF-a Biotin-free	500	AL326C	3.3pg/mL	300,000pg/mL	✓	X	X
	5000	AL326F					
	0.3µg	AL326S					
Mouse TNFa	500	AL208C	2.2pg/mL	30,000pg/mL	✓	X	X
	5000	AL208F					
	0.1µg	AL208S					
Human TNF-a Biotin-free	100	AL325HV					
	500	AL325C	0.013ng/mL	30ng/mL	✓	X	X
	5000	AL325F					
Mouse TNF-a Biotin-Free	0.1µg	AL325S					
	500	AL505C	2pg/mL	30,000pg/mL	X	✓	✓
	5000	AL505F					
Anti-Trypsin (A1AT)	0.1µg	AL505S					
	100	AL541HV					
	500	AL541C	0.007ng/mL	70ng/mL	X	✓	✓
Human NGAL	5000	AL541F					
	0.1µg	AL541S					
	100	AL344HV	Coming soon				
Human NGAL	500	AL344C					
	5000	AL344F					
	0.3µg	AL344S					
Human NGAL	100	AL364HV	Coming soon				
	500	AL364C					
	5000	AL364F					
	0.3µg	AL364S					

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
<b>内分泌疾病相关</b>							
Human Adiponectin	500	AL209C					
	5000	AL209F	45pg/mL	1,000,000pg/mL	✓	X	X
	1µg	AL209S					
Mouse Adiponectin	500	AL515C					
	5000	AL515F	3.1pg/mL	100,000pg/mL	X	✓	X
	0.3µg	AL515S					
Human Albumin	500	AL294C					
	5000	AL294F	104.8pg/mL	1,000,000pg/mL	✓	X	X
	1µg	AL294S					
Human Albumin biotin free	100	AL363HV					
	500	AL363C	102pg/mL	1,000,000pg/mL	✓		
	5000	AL363F					
Mouse Albumin	1µg	AL363S					
	500	AL508C					
	5000	AL508F	716pg/mL	1,000,000pg/mL	X	✓	X
Human Factor VIII	3µg	AL508S					
	100	AL342HV					
	500	AL342C	87pg/mL	100,000pg/mL	✓		
Human Factor VIII Biotin-Free	5000	AL342F					
	0.3µg	AL342S					
	100	AL362HV					
Human Fibrinogen	500	AL362C	71.2pg/mL	300,000pg/mL	✓		
	5000	AL362F					
	0.3µg	AL362S					
Human GLP-1	100	AL352HV					
	500	AL352C	9.2pg/mL	1,000,000pg/mL	✓	X	X
	5000	AL352F					
Human Glucagon	1µg	AL352S					
	500	AL215C	27pg/mL	30,000pg/mL	✓	✓	✓
	5000	AL215F					
Human Growth Hormone	0.1µg	AL215S					
	100	AL359HV					
	500	AL359C	6pg/mL	100,000pg/mL	✓		
Human Hyaluronic Acid	5000	AL359F					
	0.1µg	AL359S					
	100	AL312HV					
Human IGF1	500	AL312C	20pg/mL	300,000pg/mL	✓		
	5000	AL312F					
	1vial	AL312S					
Human IGF2	500	AL214C					
	5000	AL214F	5pg/mL	100,000pg/mL	✓	X	X
	0.3µg	AL214S					
Human Insulin	100	AL354HV					
	500	AL354C	0.3ng/mL	3,000ng/mL	✓		
	5000	AL354F					
Human IGF2	3µg	AL354S					
	500	AL236C	150pg/mL	100,000pg/mL	✓	X	X
	5000	AL236F					
Human IGF2	1µg	AL236S					
	500	AL237C	128pg/mL	300,000pg/mL	✓	X	✓
	5000	AL237F					
Human Insulin	1µg	AL237S					
	500	AL204C	0.8µIU/mL	3,000µIU/mL	✓	✓	✓
	5000	AL204F					
Human Insulin	0.01IU	AL204S					

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
Human Insulin HS	100	AL350HV					
	500	AL350C	0.21μIU/mL	10,000μIU/mL	✓	✓	✓
	5000	AL350F					
	0.01IU	AL350S					
Human Leptin	500	AL225C					
	5000	AL225F	50pg/mL	100,000pg/mL	✓	X	X
	0.3μg	AL225S					
Mouse Leptin	500	AL521C					
	5000	AL521F	16pg/mL	300,000pg/mL	✓	✓	✓
	1μg	AL521S					
Human Prolactin	500	AL227C					
	5000	AL227F	125pg/mL	300,000pg/mL	✓	X	X
	1μg	AL227S					
Human C-peptide	500	AL299C					
	5000	AL299F	23.5pg/mL	30,000pg/mL	✓	X	X
	0.1μg	AL299S					
Mouse C-peptide	500	AL510C					
	5000	AL510F	11.1pg/mL	30,000pg/mL	✓	✓	✓
	0.1μg	AL510S					
Human BMP-7	100	AL369HV					
	500	AL369C					
	5000	AL369F	4pg/mL	300,000pg/mL	✓	✓	
	0.3μg	AL369S					
Human COL1A1	100	AL371HV					
	500	AL371C					
	5000	AL371F	6pg/mL	300,000pg/mL	✓		
Human Perlecan	0.3μg	AL371S					
	100	AL377HV					
	500	AL377C					
Human Laminin	5000	AL377F	4.4pg/mL	300,000pg/mL	✓		
	1μg	AL377S					
	100	AL368HV					
Human E-Cadherin	500	AL368C	0.1pg/mL	30,000pg/mL	✓	X	
	5000	AL368F					
	3μg	AL368S					
Human TGF-β1	100	AL370HV					
	500	AL370C					
	5000	AL370F	3.5pg/mL	300,000pg/mL	✓	✓	✓
	0.3μg	AL370S					
<b>血管新生靶点</b>							
Human TGF-β1 Biotin-Free	100	AL336HV					
	500	AL336C	9pg/mL	100,000pg/mL	✓	✓	
	5000	AL336F					
	0.3μg	AL336S					
Human VEGFA	100	AL361HV					
	500	AL361C					
	5000	AL361F	9.2pg/mL	300,000pg/mL	✓		
Human VEGFA	0.3μg	AL361S					
	500	AL201C					
	5000	AL201F	2.2pg/mL	100,000pg/mL	✓	X	X
Mouse VEGFA	0.3μg	AL201S					
	500	AL520C					
	5000	AL520F	1.5pg/mL	100,000pg/mL	✓	✓	✓
Human VEGFB	0.1μg	AL520S					
	500	AL251C					
	5000	AL251F	140pg/mL	1,000,000pg/mL	✓	✓	
	1μg	AL251S					

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
Mouse VEGFC	500	AL252C					
	5000	AL252F	6.2pg/mL	100,000pg/mL	✓	X	X
	0.3μg	AL252S					
	500	AL253C					
Human VEGFD	5000	AL253F	29.9pg/mL	300,000pg/mL	✓	✓	
	0.15μg	AL253S					
	<b>肿瘤标志物</b>						
Human AFP	500	AL230C					
	5000	AL230F	40.3pg/mL	300,000pg/mL	✓		
	0.3μg	AL230S					
Human β-NGF	500	AL246C					
	5000	AL246F	32pg/mL	300,000pg/mL	✓	✓	X
	1μg	AL246S					
Human CA125	500	AL256C					
	5000	AL256F	1.5U/mL	3,000U/mL	✓		
	1500U	AL256S					
Human E-Cadherin	100	AL370HV					
	500	AL370C	3.5pg/mL	300,000pg/mL	✓	✓	✓
	5000	AL370F					
Human Caspase-3 active	0.3μg	AL370S					
	500	AL278C					
	5000	AL278F	126pg/mL	300,000pg/mL	✓	X	X
Human CXCL11/I-TAC	0.3μg	AL278S					
	500	AL260C					
	5000	AL260F	1.4pg/mL	30,000pg/mL	✓	X	
Human EGF	0.1μg	AL260S					
	100	AL339HV					
	500	AL339C	0.24pg/mL	100,000pg/mL	✓		
Human EGFR	5000	AL339F					
	0.1μg	AL339S					
	100	AL340HV					
Human EPO	500	AL340C	1.46pg/mL	300,000pg/mL	✓	✓	
	5000	AL340F					
	0.3μg	AL340S					
Human ERBB2/HER2	500	AL206C					
	5000	AL206F	1mIU/mL	30,000mIU/mL	✓	X	X
	30IU	AL206S					
Human Fibronectin (h Plasma)	500	AL235C					
	5000	AL235F	54.1pg/mL	300,000pg/mL	✓		
	1μg	AL235S					
Human HGFR/c-MET	100	AL351HV					
	500	AL351C	148pg/mL	1,000,000pg/mL	✓	X	X
	5000	AL351F					
Human IFN-β	1μg	AL351S					
	500	AL281C					
	5000	AL281F	23.1pg/mL	1,000,000pg/mL	✓		
Human VEGFR	3μg	AL281S					
	500	AL265C					
	5000	AL265F	9.6pg/mL	100,000pg/mL	✓	X	X
Human VEGFR	0.3μg	AL265S					

# AlphaLISA：只需混合无需洗涤的ELISA

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
Human MMP1	500	AL242C	82.6pg/mL	300,000pg/mL	✓	X	
	5000	AL242F					
	0.3μg	AL242S					
Human MMP2	500	AL269C	260pg/mL	300,000pg/mL	✓	X	C
	5000	AL269F					
	0.3μg	AL269S					
Human MMP3	500	AL284C	25.8pg/mL	300,000pg/mL	✓	X	
	5000	AL284F					
	0.3μg	AL284S					
Human MMP7	100	AL378HV					
	500	AL378C	35pg/mL	1,000,000pg/mL	✓	X	
	5000	AL378F					
Human MMP8	1μg	AL378S					
	100	AL374HV					
	500	AL374C	10pg/mL	100,000pg/mL	✓		
Human MMP9	5000	AL374F					
	0.1μg	AL374S					
	500	AL243C	10.9pg/mL	300,000pg/mL	✓	X	
Human MMP9	5000	AL243F					
	0.3μg	AL243S					
Mouse MMP9	500	AL519C	19.9pg/mL	300,000pg/mL	X	✓	✓
	5000	AL519F					
	0.3μg	AL519S					
Human MMP12	100	AL375HV					
	500	AL375C	6.1pg/mL	10,000pg/mL	✓		
	5000	AL375F					
Human MMP13	0.1μg	AL375S					
	100	AL318HV					
	500	AL318C	6.5pg/mL	100,000pg/mL	✓		
Human Neprilysin	5000	AL318F					
	0.1μg	AL318S					
	100	AL337HV					
Human Nidogen-1	500	AL337C	8.9pg/mL	300,000pg/mL	✓	✓	
	5000	AL337F					
	1μg	AL337S					
Human Oncostatin M	100	AL367HV					
	500	AL367C	17pg/mL	300,000pg/mL	✓		
	5000	AL367F					
Human PIP Collagen	0.3μg	AL367S					
	100	AL341HV					
	500	AL341C	0.6pg/mL	100,000pg/mL	✓	X	
Human PSA	5000	AL341F					
	0.1μg	AL341S					
	100	AL353HV					
Human TFF3	500	AL353C	10pg/mL	300,000pg/mL	✓		
	5000	AL353F					
	0.3μg	AL353S					
Human TIMP1	500	AL228C					
	5000	AL228F	1.2ng/mL	1,000ng/mL	✓		
	3μg	AL228S					
Human VCAM-1	500	AL273C					
	5000	AL273F	6.1pg/mL	30,000pg/mL	✓		
	0.1μg	AL273S					
Human Fibrinogen	500	AL272C	520pg/mL	100,000pg/mL	✓	X	X
	5000	AL272F					
	0.1μg	AL272S					
Human VCAM-1	100	AL338HV					
	500	AL338C	0.92pg/mL	300,000pg/mL	✓		
	5000	AL338F					
Human Transferin	1μg	AL338S					
	100	AL352HV					
	500	AL352C	9.2pg/mL	1,000,000pg/mL	✓	X	X
	5000	AL352F					
	1μg	AL352S					

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
肿瘤免疫治疗							
Human PD-1	100	AL343HV					
	500	AL343C	9pg/mL	100,000pg/mL	✓	X	
	5000	AL343F					
Human PDL-1	0.3μg	AL343S					
	100	AL355HV					
	500	AL355C	2pg/mL	300,000pg/mL	✓	X	
Human PD-1 and PDL-1 Binding	5000	AL355F					
	0.3μg	AL355S					
	100	AL356HV					
Human D-dimer	500	AL356C	2nM				
	5000	AL356F					
	1.3μg/2.1μg	AL356S					
心血管疾病相关							
Human α2 macroglobulin	500	AL277C	575pg/mL	1,000,000pg/mL	✓	X	X
	5000	AL277F					
	3μg	AL277S					
Human Cardiac Troponin I	500	AL295C	4.7pg/mL	100,000pg/mL	✓	X	
	5000	AL295F					
	0.1μg	AL295S					
Human D-dimer	500	AL290C	41.3pg/mL	1,000,000pg/mL	✓		
	5000	AL290F					
	1μg	AL290S					
Human ICAM-1	500	AL282C	3.4pg/mL	100,000pg/mL	✓	X	X
	5000	AL282F					
	1μg	AL282S					
Human Myeloperoxidase	500	AL245C	67.1pg/mL	300,000pg/mL	✓	✓	X
	5000	AL245F					
	1μg	AL245S					
Human Myoglobin	500	AL285C	1pg/mL	100,000pg/mL	✓	X	X
	5000	AL285F					
	0.3μg	AL285S					
Human PAI-1	500	AL286C	39pg/mL	300,000pg/mL	✓	X	X
	5000	AL286F					
	1μg	AL286S					
Human PCSK9	500	AL270C	1056pg/mL	300,000pg/mL	✓	X	
	5000	AL270F					
	0.3μg	AL270S					
Human Plasminogen	500	AL248C	4.4ng/mL	10,000ng/mL	✓	X	
	5000	AL248F					
	10μg	AL248S					
Human Prorenin	500	AL249C	8.1pg/mL	300,000pg/mL	✓		
	5000	AL249F					
	0.3μg	AL249S					
Human tPA	500	AL250C	11.1pg/mL	300,000pg/mL	✓	X	X
	5000	AL250F					
	1μg	AL250S					
Human Transferin	100	AL311HV					
	500	AL311C	29pg/mL	1,000,000pg/mL			✓
	5000	AL311F					
	1vial	AL311S					

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
<b>CNS相关</b>							
Human Amyloid $\beta$ 1-15	500	AL274C	8.1pg/mL	10,000pg/mL	✓		
	5000	AL274F					
	0.1 $\mu$ g	AL274S					
Human Amyloid $\beta$ 1-40	500	AL202C	88pg/mL	100,000pg/mL	✓	✓	✓
	5000	AL202F					
	0.3 $\mu$ g	AL202S					
Human Amyloid beta 1-40 HS	100	AL373HV					
	500	AL373C	42pg/mL	10,000pg/mL	✓	X	X
	5000	AL373F					
Mouse Amyloid $\beta$ 1-40	500	AL506C	31.7pg/mL	30,000pg/mL	x	✓	✓
	5000	AL506F					
	0.3 $\mu$ g	AL506S					
Human Amyloid $\beta$ 1-42	500	AL203C					
	5000	AL203F	300pg/mL	300,000pg/mL	✓	✓	✓
	1 $\mu$ g	AL203S					
Human Amyloid $\beta$ 1-42(HS)	500	AL276C	20.8pg/mL	30,000pg/mL	✓	✓	
	5000	AL276F					
	0.1 $\mu$ g	AL276S					
Mouse Amyloid $\beta$ 1-42	500	AL507C	25.2pg/mL	100,000pg/mL	x	✓	✓
	5000	AL507F					
	0.3 $\mu$ g	AL507S					
Human Amyloid $\beta$ 1-x	500	AL288C	4.5pg/mL	30,000pg/mL	✓		
	5000	AL288F					
	5 $\mu$ g	AL288S					
Human Amyloid Oligomers HS	500	AL334C	42pg/mL	300,000pg/mL	✓		
	5000	AL334F					
	5 $\mu$ g	AL334S					
Human A $\beta$ 42	100	AL330HV					
	500	AL330C	14.7pg/mL	10,000pg/mL	✓		
	5000	AL330F					
Human BDNF	1 $\mu$ g	AL330S					
	100	AL317HV					
	500	AL317C	3.3pg/mL	300,000pg/mL	✓		
Human Tau	5000	AL317F					
	0.3 $\mu$ g	AL317S					
	500	AL271C					
Human CCL2	5000	AL271F	14.3pg/mL	300,000pg/mL	✓		
	0.3 $\mu$ g	AL271S					
	500	AL244C	3.8pg/mL	100,000pg/mL	✓	x	x
Mouse CCL2/MCP-1	5000	AL244F					
	0.3 $\mu$ g	AL244S					
	500	AL509C	6.9pg/mL	100,000pg/mL	x	✓	✓
Human CCL2/MCP-1 $\alpha$	5000	AL509F					
	0.3 $\mu$ g	AL509S					
	500	AL257C					
Human CCL2/MCP-1 $\beta$	5000	AL257F	0.8pg/mL	30,000pg/mL	✓	x	x
	0.3 $\mu$ g	AL257S					
	500	AL258C	2.8pg/mL	100,000pg/mL	✓	x	x
Human CCL5/RANTES	5000	AL258F					
	0.1 $\mu$ g	AL258S					
	500	AL289C	0.8pg/mL	100,000pg/mL	✓	x	x
	5000	AL289F					
	0.3 $\mu$ g	AL289S					

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠
<b>激素检测</b>							
Mouse CCL5/RANTES	500	AL516C	2.6pg/mL	30,000pg/mL	x	✓	✓
	5000	AL516F					
	0.3 $\mu$ g	AL516S					
Human Frataxin	100	AL322HV					
	500	AL322C	8pg/mL	300,000pg/mL	✓	✓	
	5000	AL322F					
Human GDNF	0.3 $\mu$ g	AL322S					
	100	AL319HV					
	500	AL319C	3.8pg/mL	100,000pg/mL	✓		
Human TDP-43	5000	AL319F					
	0.1 $\mu$ g	AL319S					
	100	AL357HV					
Cortisol	500	AL357C	14.2pg/mL	300,000pg/mL	✓		
	5000	AL357F					
	1 $\mu$ g	AL357S					
<b>自体吞噬</b>							
Human Progesterone	100	AL314HV					
	500	AL314C	177pg/mL	3,600,000pg/mL			
	5000	AL314F					
Human Somatostatin	100 $\mu$ M	AL314S					
	100	AL335HV					
	500	AL335C	0.59nM	1,000nM	✓		
Human testosterone	5000	AL335F					
	100 $\mu$ M	AL335S					
	100	AL329HV					
Human P62	500	AL329C	55pg/mL	1,637,000pg/mL	✓		
	5000	AL329F					
	2nmol	AL329S					
Human LC38	100	AL324HV					
	500	AL324C	0.7nM	875nM	✓		
	5000	AL324F					
	20 $\mu$ M	AL324S					
<b>病毒类检测</b>							
Human HIV p24	500	AL305C	929pg/mL	3,000,000pg/mL	✓		
	5000	AL305F					
	3 $\mu$ g	AL305S					
HIV P24 Biotin-Free	500	AL306C	7.9ng/mL	3,000ng/mL	✓	✓	C
	5000	AL306F					
	3 $\mu$ g	AL306S					

# AlphaLISA: 只需混合无需洗涤的ELISA

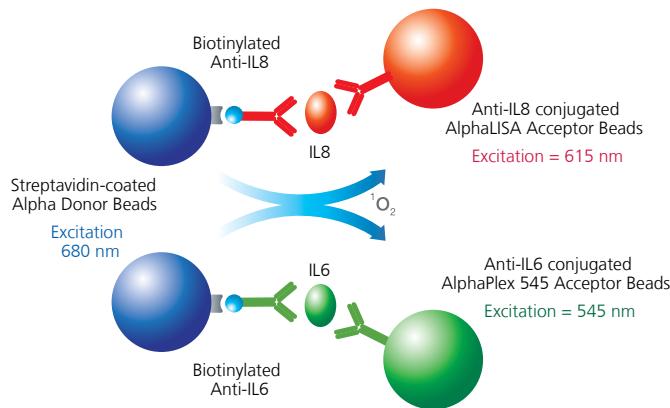
产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应						
					人	小鼠	大鼠				
抗体类检测											
抗体滴度											
Human total IgG	500	AL205C	0.24ng/mL	1,000ng/mL	✓	X	X				
	5000	AL205F									
	3µg	AL205S									
Mouse total IgG	100	AL529HV									
	500	AL529C	36pg/mL	3,000,000pg/mL		✓					
	5000	AL529F									
	0.1µg	AL529S									
Human FCGR3A/CD16a (176Phe/F158)	100	AL347HV									
	500	AL347C	22.6pM								
	5000	AL347F									
	1.5µg	AL347S									
Human FCGR3A/CD16a (176Val/V158)	100	AL348HV									
	500	AL348C	4.3pM								
	5000	AL348F									
	1µg	AL348S									
抗体分型											
Human IgG1 (isotyping)	100	AL307HV									
	500	AL307C	47pg/mL	1,000,000pg/mL	✓		X				
	5000	AL307F									
	1µg	AL307S									
Human IgG2 (isotyping)	100	AL308HV									
	500	AL308C	6.65ng/mL	100,000ng/mL	✓						
	5000	AL308F									
	100µg	AL308S									
Human IgG3 (isotyping)	100	AL309HV									
	500	AL309C	500pg/mL	3,000,000pg/mL	✓						
	5000	AL309F									
	3µg	AL309S									
Human IgG4 (isotyping)	100	AL310HV									
	500	AL310C	300pg/mL	500,000pg/mL	✓						
	5000	AL310F									
	10µg	AL310S									
Human IgA	500	AL262C	449pg/mL	3,000,000pg/mL	✓	X					
	5000	AL262F									
	10µg	AL262S									
Human IgM	500	AL263C	511pg/mL	3,000,000pg/mL	✓	X	X				
	5000	AL263F									
	3µg	AL263S									
Human IgG1 (isotyping)	100	AL522HV									
	500	AL522C	117pg/mL	3,000,000pg/mL	X	✓					
	5000	AL522F									
	3µg	AL522S									
Mouse IgG2a (isotyping)	100	AL523HV									
	500	AL523C	23pg/mL	1,000,000pg/mL		✓					
	5000	AL523F									
	3µg	AL523S									
Mouse IgG2b (isotyping)	100	AL524HV									
	500	AL524C	60pg/mL	1,000,000pg/mL		✓					
	5000	AL524F									
	1µg	AL524S									
Mouse IgG3 (isotyping)	100	AL525HV									
	500	AL525C	60pg/mL	1,000,000pg/mL		✓					
	5000	AL525F									
	1µg	AL525S									

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应						
					人	小鼠	大鼠				
PK/PD											
杂质检测											
Anti-PEG IgM	100	AL313HV									
	500	AL313C	6.65ng/mL	100,000ng/mL							
	5000	AL313F									
	0.1mg	AL313S									
安全性检测											
Mouse Albumin	500	AL508C	716pg/mL	1,000,000pg/mL	X	✓	X				
	5000	AL508F									
	3µg	AL508S									
Anti-PEG IgM	100	AL313HV									
	500	AL313C	6.65ng/mL	100,000ng/mL							
	5000	AL313F									
	100µg	AL313S									
Human Transferin	100	AL311HV									
	500	AL311C	29pg/mL	1,000,000pg/mL		✓					
	5000	AL311F									
	1 vial	AL311S									

产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应			产品名称	规格 检测点	货号	检出下限	检出上限	交叉反应		
					人	小鼠	大鼠						人	小鼠	大鼠
动物健康															
Bovine Immunoglobulin A	100	AL53HV						Cry1Ab/1Ac	100	AL401HV	Coming soon				
	500	AL536C	0.25ng/mL	3,000pg/mL	X				500	AL401C					
	5000	AL536F							5000	AL401F					
	3µg	AL536S							0.5µg	AL401S					
Bovine total IgG	100	AL530HV						Cry1F	100	AL402HV					
	500	AL530C	530pg/mL	3000,000pg/mL	X	X			500	AL402C	111pg/mL	300,000pg/mL			
	5000	AL530F							5000	AL402F					
	0.1µg	AL530S							0.5µg	AL402S					
Bovine IgG1	100	AL531HV						Cry2	100	AL403HV					
	500	AL531C	1,700pg/mL	3000,000pg/mL	X				500	AL403C	12pg/mL	300,000pg/mL			
	5000	AL531F							5000	AL403F					
	3µg	AL531S							0.5µg	AL403S					
Bovine IgG2	100	AL532HV						Chloramphenicol	100	AL385HV	Coming soon				
	500	AL532C	2,100pg/mL	3000,000pg/mL					500	AL385C					
	5000	AL532F							5000	AL385F					
	3µg	AL532S							0.5µg	AL385S					
Bovine IgM	100	AL533HV													
	500	AL533C	450pg/mL	10,000pg/mL											
	5000	AL533F													
	3µg	AL533S													
Bovine IL-1α	100	AL537HV													
	500	AL537C	2.58pg/mL	300,000pg/mL	X										
	5000	AL537F													
	0.3µg	AL537S													
Bovine IL-1β	100	AL539HV													
	500	AL539C	20pg/mL	300,000pg/mL	X										
	5000	AL539F													
	0.3µg	AL539S													
Bovine IL-2	100	AL543HV													
	500	AL543C	30pg/mL	1,000,000pg/mL											
	5000	AL543F													
	1µg	AL543S													
Bovine IL-4	100	AL542HV													
	500	AL542C	1.8pg/mL	100,000pg/mL											
	5000	AL542F													
	0.1µg	AL542S													
Bovine IL-6	100	AL538HV													
	500	AL538C	0.01ng/mL	37ng/mL											
	5000	AL538F													
	0.3µg	AL538S													
Bovine/Ovine IL-8	100	AL545HV	Coming soon												
	500	AL545C													
	5000	AL545F													
	1µg	AL545S													
Bovine/Ovine IL-10	100	AL546HV	Coming soon												
	500	AL546C													
	5000	AL546F													
	1µg	AL546S													
Bovine IL-17A	100	AL547HV													
	500	AL547C	Coming soon												
	5000	AL547F													
	1µg	AL547S													
Bovine TNF-α	100	AL534HV													
	500	AL534C	1pg/mL	300,000pg/mL											
	5000	AL534F													
	0.3µg	AL534S													
Bovine IFN-γ	100	AL535HV													
	500	AL535C	10pg/mL	300,000pg/mL											
	5000	AL535F													
	0.3µg	AL535S													

# AlphaPlex (Tb/Sm) 多标记免疫分析试剂

该免疫分析试剂盒使用AlphaPlex545受体微珠。可以与其他AlphaLISA免疫分析试剂盒联合使用。实现单孔样品检测两种靶点。



## AlphaPlex检测原理

生物素化的抗IL-6抗体和生物素化的抗IL-8抗体分别与链霉亲和素偶联的供体微珠结合。另一个位点的抗IL-6抗体和抗IL-8抗体分别偶联在AlphaLISA 和 AlphaPlex545受体微珠上。抗体与分析物结合后拉近了受体和供体微珠。在680nm激光激发下供体微珠激发出活性单线态氧，受体微珠接收后分别产生615nm(AlphaLISA)和545nm(AlphaPlex)。可以使用Envision的Alpha模块进行信号检测。

产品	规格(点数)	货号
免疫炎症相关		
AlphaPlex 545 Human IFN- $\gamma$ Kit	100	AP217TB-HV
	500	AP217TB-C
	5,000	AP217TB-F
AlphaPlex 545 Human IL-1 $\beta$ Kit	100	AP220TB-HV
	500	AP220TB-C
	5,000	AP220TB-F
AlphaPlex 545 Human IL2 Kit	100	AP221TB-HV
	500	AP221TB-C
	5,000	AP221TB-F
AlphaPlex 545 Human IL6 Kit	100	AP223TB-HV
	500	AP223TB-C
	5,000	AP223TB-F
AlphaPlex 545 Human IL17 Kit	100	AP219TB-HV
	500	AP219TB-C
	5,000	AP219TB-F
AlphaPlex 545 Human TNF- $\alpha$ Kit	100	AP208TB-HV
	500	AP208TB-C
	5,000	AP208TB-F
AlphaPlex 545 Human/rat IL22 Kit	100	AP320TB-HV
	500	AP320TB-C
	5,000	AP320TB-F
AlphaPlex 545 Human IP10 Kit	100	AP326TB-HV
	500	AP326TB-C
	5,000	AP326TB-F
内分泌疾病相关		
AlphaPlex 545 Human Insulin Kit	100	AP204TB-HV
	500	AP204TB-C
	5,000	AP204TB-F
心血管疾病相关		
AlphaPlex 545 Human ICAM-1 Kit	100	AP282TB-HV
	500	AP282TB-C
	5,000	AP282TB-F

产品	规格(点数)	货号
免疫炎症相关		
AlphaPlex 645 Human IFN- $\gamma$ Kit	100	AP217SM-HV
	500	AP217SM-C
	5,000	AP217SM-F
AlphaPlex 645 Human IL-1 $\beta$ Kit	100	AP220SM-HV
	500	AP220SM-C
	5,000	AP220SM-F
AlphaPlex 645 Human IL2 Kit	100	AP221SM-HV
	500	AP221SM-C
	5,000	AP221SM-F
AlphaPlex 645 Human IL17 Kit	100	AP219SM-HV
	500	AP219SM-C
	5,000	AP219SM-F
AlphaPlex 645 Human TNF- $\alpha$ Kit	100	AP208SM-HV
	500	AP208SM-C
	5,000	AP208SM-F
AlphaPlex 645 Human IP10 Kit	100	AP326SM-HV
	500	AP326SM-C
	5,000	AP326SM-F
内分泌疾病相关		
AlphaPlex 645 Human Insulin Kit	100	AP204SM-HV
	500	AP204SM-C
	5,000	AP204SM-F
心血管疾病相关		
AlphaPlex 645 Human ICAM-1 Kit	100	AP282SM-HV
	500	AP282SM-C
	5,000	AP282SM-F

试剂盒成分：生物素化抗体，抗体偶联的AlphaPlex受体微珠，链霉亲和素偶联的供体微珠，标准品，反应溶液

试剂盒规格：100点试剂盒每孔反应体系为100 $\mu$ L, 500/5000点试剂盒每孔反应体系为50 $\mu$ L

保存条件：4°C

## AlphaLISA 结合检测试剂盒 新品

检测配体与受体结合行为的检测试剂盒，可用于抑制剂研究等。

### 特点

- 高信噪比

使用 AlphaLISA 进行检测可获得非常高的信噪比。

- 可随时使用

试剂盒中包含有检测所需的试剂和方案，可立即开始实验。

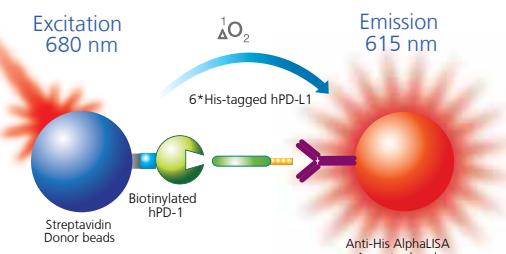
- 高通量

检测中仅需进行混合操作，非常适合筛选和大量样品处理。

### 免疫检查点

检测人 PD-1 与 PD-L1 结合行为的试剂盒。

我们还提供检测人 PD-1 与 PD-L1 的免疫测定试剂盒。（请参见第 10 页免疫检查点相关信息。）



产品名称	数量 (检测点)	产品货号
AlphaLISA	100	AL356HV
PD-1 and PD-L1 Binding kit	500	AL356C
	10,000	AL356F

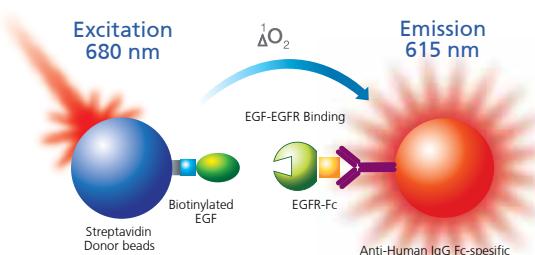
成分：抗 His AlphaLISA 受体微珠、链霉亲合素供体微珠、生物素化人 PD-1、His-PD-L1、AlphaLISA 免疫测定缓冲液 (10x)。

数量：1 检测点相当于按试剂盒推荐量 (AL356HV 为 100  $\mu\text{L}$ 、AL356C/F 为 40  $\mu\text{L}$ ) 使用时，1/2 AreaPlate 96 孔板中 1 个孔的用量。

保存：+ 4°C

### 受体 - 配体结合

检测人 EGF 和 EGFR 结合行为的试剂盒。

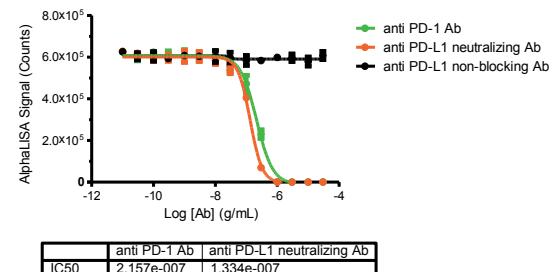


产品名称	数量 (检测点)	产品货号
AlphaLISA	100	AL366HV
EGF and EGFR Binding kit	500	AL366C
	10,000	AL366F



S/N	1600 (存在 5 nM PD-1 时)
Kd (app)	2.5 nM (存在 5 nM PD-1 时)

#### PD-1/PD-L1 结合抑制试验

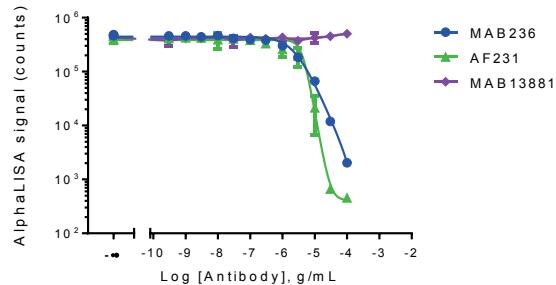


存在人 PD-1 (5 nM) 和人 PD-L1 (5 nM) 时，加入抗 PD-1 抗体（绿色）和抗 PD-L1 中和抗体（橙色）进行反应，观察到 PD-1/PD-L1 的结合受到抑制，信号下降。

#### 试剂盒性能

S/N	947 (存在 3.125 nM EGFR 时)
Kd (app)	2.5 nM (存在 3.125 nM EGFR 时)

#### EGF/EGFR 结合抑制试验



存在 EGF (3.5 nM) 和 EGFR (3.5 nM) 时，加入抗人 EGF 抗体（绿色：AF231）和抗人 EGFR 抗体（蓝色：MAB235）进行反应。观察到 EGF/EGFR 的结合受到抑制，信号下降，IC50 分别为 3.9  $\mu\text{g}/\text{mL}$  和 2.2  $\mu\text{g}/\text{mL}$ 。

## 细胞磷酸化检测 Alpha SureFire

*Surefire* 是一种用“Mix Only”的方式来检测细胞内特异性被激活的磷酸化蛋白的方法。

### Alpha SureFire / SureFire Ultra 新品

在传统的AlphaScreen *Surefire* 检测方法中，用于调控细胞的抗体会与Protein A竞争结合，从而抑制AlphaScreen *Surefire* 的信号。新的AlphaLISA *Surefire Ultra*采用成熟的CaptSure™技术代替Protein A，不仅实现了试剂盒抗体捕获的特异性，而且具有更低的检测背景。新的试剂盒利用AlphaLISA受体微珠代替AlphaScreen受体微珠，对血红蛋白淬灭干扰具有更加出色的光谱特性。

### 检测特点：

- 只需混合( Mix & Measure)

细胞样本经过刺激、裂解后，只需要将试剂组分加入裂解液，无需任何多余的洗涤步骤。

- 高灵敏度

检测灵敏度大于等于Western Blot方法；可以检测过表达蛋白靶点及内源水平表达的蛋白；可以使用稳定表达细胞系或原代细胞及干细胞进行试验。

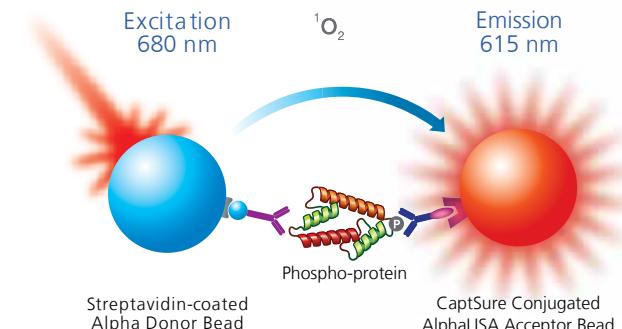
- 满意的Z'值

可以兼容96、384或1536孔板的不同检测模式，均可得到很高的Z'值。

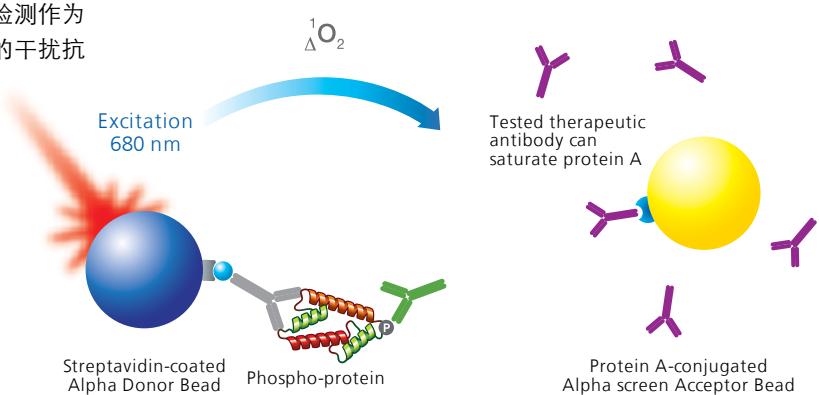
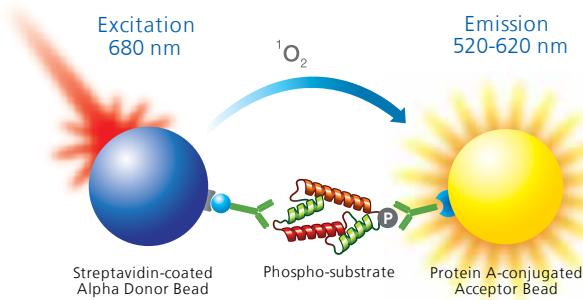
- 适用于抗体药物的开发使用，适应组织裂解液样品 (*Surefire Ultra*)

在药物研发过程中，我们进行细胞学水平磷酸化蛋白检测作为功能学研究方法。当待测药物为抗体分子时，样品中的干扰抗体与Protein A进行非特异性结合从而干扰检测信号（如右下图）。*Surefire Ultra*方法使用新的Captsure标签，有效的避免了这种干扰。而且，与传统的*Surefire*检测相比，*Surefire Ultra*也更加适合应用于肝脏这类器官、组织样品的检测（当操作组织样品时，请进行匀浆及生物素剔除等预处理）。

### AlphaLISA Surefire Ultra 新品



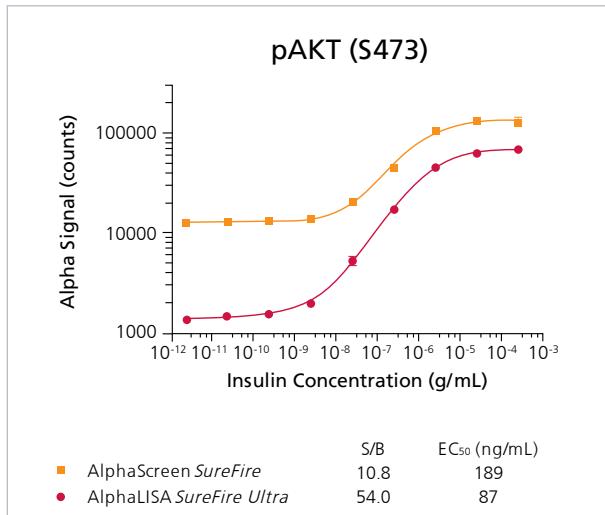
### AlphaScreen Surefire



## 检测步骤



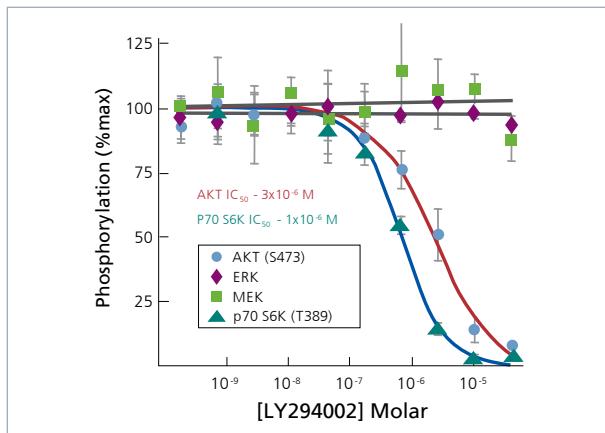
## Alpha Surefire及Surefire Ultra 检测结果对比



### 结果对比

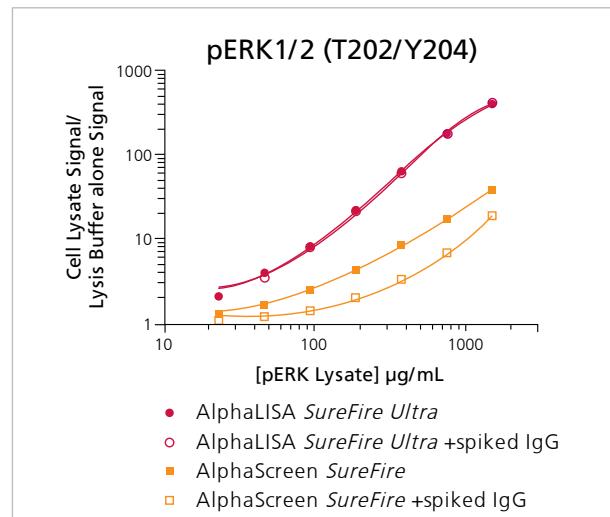
MCF7细胞(40,000细胞/孔)培养过夜，2小时无血清培养。用insulin刺激30分钟后裂解细胞。同时用AKT(S473) Surefire及 Surefire Ultra试剂盒检测。Surefire Ultra试剂盒得到的S/B ratio更高，检测结果表现出更高的灵敏度。

## AlphaScreen 激酶检测示例



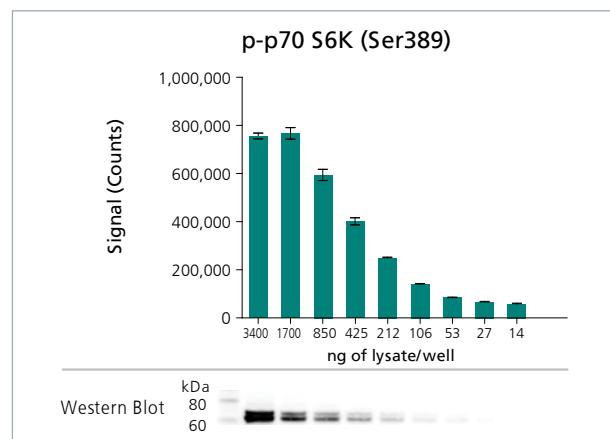
### PI3K 通路抑制剂特异性检测

NIH-3T3细胞经过 PI3K 通路抑制剂 LY294002处理后再用PDGF刺激。分别用AlphaLISA Surefire Ultra试剂盒检测磷酸化AKT (Ser473)、P70S6K (T389) , ERK (T202/Y204) , MEK1 (S217/S221) 蛋白。ERK通路蛋白（磷酸化ERK及MEK）没有受到抑制剂的作用，但观察到PI3K通路蛋白（磷酸化AKT及p70S6K）受到特异性的磷酸化表达抑制。



### 与IgG混合后检测结果对比

EGF刺激A431细胞后，在裂解液中加入rabbit IgG (10 µg/mL)，分别用Surefire和Surefire Ultra磷酸化ERK(T202/204)检测试剂盒检测梯度稀释的细胞裂解液，Surefire试剂盒的检测信号在IgG抗体的影响下有所降低，但是IgG抗体的加入对Surefire Ultra的检测信号不产生任何影响。



### AlphaScreen SureFire vs 高灵敏 Western Blot

用IGF1刺激MCF7细胞后裂解，分别用western lightening及AlphaLISA Surefire Ultra试剂盒检测细胞内源磷酸化p70S6K (Ser389) 蛋白的结果对比。

# AlphaLISA SureFire Ultra 试剂盒

## AlphaLISA SureFire Ultra 检测试剂盒 新品

AKT 通路		
产品名称	检测点	货号
AlphaLISA Surefire Ultra p-AKT1/2/3 (T308) kit	100	ALSU-PAKT-A-HV
	500	ALSU-PAKT-A500
	10,000	ALSU-PAKT-A10K
	50,000	ALSU-PAKT-A50K
AlphaLISA Surefire Ultra p-AKT 1/2/3 (S473) Kit	100	ALSU-PAKT-B-HV
	500	ALSU-PAKT-B500
	10,000	ALSU-PAKT-B10K
	50,000	ALSU-PAKT-B50K
AlphaLISA Surefire Ultra Total AKT1 Kit	100	ALSU-TAKT-B-HV
	500	ALSU-TAKT-B500
	10,000	ALSU-TAKT-B10K
	50,000	ALSU-TAKT-B50K

NFkB 通路		
产品名称	检测点	货号
AlphaLISA Surefire Ultra p-IKK $\alpha$ (Ser176/180) Kit	100	ALSU-PIKKA-A-HV
	500	ALSU-PIKKA-A500
	10,000	ALSU-PIKKA-A10K
	50,000	ALSU-PIKKA-A50K
AlphaLISA Surefire Ultra p-NFkB (S536) Kit	100	ALSU-PNFKB-A-HV
	500	ALSU-PNFKB-A500
	10,000	ALSU-PNFKB-A10K
	50,000	ALSU-PNFKB-A50K

TGF $\beta$ /SMAD 通路		
产品名称	检测点	货号
AlphaLISA Surefire Ultra SMAD1 (Ser463/465) Kit	100	ALSU-PSM1-A-HV
	500	ALSU-PSM1-A500
	10,000	ALSU-PSM1-A10K
	50,000	ALSU-PSM1-A50K
AlphaLISA Surefire Ultra SMAD3 (Ser423/425) Kit	100	ALSU-PSM3-A-HV
	500	ALSU-PSM3-A500
	10,000	ALSU-PSM3-A10K
	50,000	ALSU-PSM3-A50K

MAPK 通路		
产品名称	检测点	货号
AlphaLISA Surefire Ultra p-ERK (T202/Y204) Kit	100	ALSU-PERK-A-HV
	500	ALSU-PERK-A500
	10,000	ALSU-PERK-A10K
	50,000	ALSU-PERK-A50K
AlphaLISA Surefire Ultra Total ERK1/2 kit	100	ALSU-TERK-A-HV
	500	ALSU-TERK-A500
	10000	ALSU-TERK-A10K
	50000	ALSU-TERK-A50K
AlphaLISA Surefire Ultra p-p38 MAPK (T180/Y182) Kit	100	ALSU-PP38-A-HV
	500	ALSU-PP38-A500
	10000	ALSU-PP38-A10K
	50000	ALSU-PP38-A50K
AlphaLISA Surefire Ultra Total p38 MAPK Kit	100	ALSU-TP38-A-HV
	500	ALSU-TP38-A500
	10000	ALSU-TP38-A10K
	50000	ALSU-TP38-A50K
AlphaLISA Surefire Ultra p-JNK1/2/3 (Thr183/Tyr185)	100	ALSU-PJNK-A-HV
	500	ALSU-PJNK-A500
	10000	ALSU-PJNK-A10K
	50000	ALSU-PJNK-A50K
AlphaLISA Surefire Ultra p-MEK1(Ser218/222)	100	ALSU-PMEK1-A-HV
	500	ALSU-PMEK1-A500
	10000	ALSU-PMEK1-A10K
	50000	ALSU-PMEK1-A50K

mTOR 信号/转录调控试剂盒		
产品名称	检测点	货号
AlphaLISA Surefire Ultra 4E-BP1(Thr37/46)	100	ALSU-P4EBP-A-HV
	500	ALSU-P4EBP-A500
	10,000	ALSU-P4EBP-A10K
	50,000	ALSU-P4EBP-A50K
AlphaLISA Surefire Ultra p70 S6K (T389) Kit	100	ALSU-PP70-A-HV
	500	ALSU-PP70-A500
	10,000	ALSU-PP70-A10K
	50,000	ALSU-PP70-A50K
AlphaLISA Surefire Ultra p-eIF2 $\alpha$ (S51) Kit	100	ALSU-PEIF2-A-HV
	500	ALSU-PEIF2-A500
	10,000	ALSU-PEIF2-A10K
	50,000	ALSU-PEIF2-A50K
AlphaLISA Surefire Ultra p-eIF4E (S209) Kit	100	ALSU-PEIF4-A-HV
	500	ALSU-PEIF4-A500
	10,000	ALSU-PEIF4-A10K
	50,000	ALSU-PEIF4-A50K
AlphaLISA Surefire Ultra p-mTOR(Ser2448) Kit	100	ALSU-PMTOR-A-HV
	500	ALSU-PMTOR-A500
	10,000	ALSU-PMTOR-A10K
	50,000	ALSU-PMTOR-A50K
AlphaLISA Surefire Ultra p-RPS6(Ser240/244) Kit	100	ALSU-PS6R-A-HV
	500	ALSU-PS6R-A500
	10,000	ALSU-PS6R-A10K
	50,000	ALSU-PS6R-A50K

试剂盒组分：Lysis buffer, Activation Buffer, Dilution Buffer, Reaction Buffer (含抗体), Streptavidin Donor beads, CaptSure Acceptor beads, Positive control lysate.

检测点数：100检测点的每孔检测量是以 1/2 AreaPlate 96孔板中每孔检测总体积为60ul的体系计算；500, 10,000及50,000检测点的每孔检测量是以384孔板中每孔检测总体积为20ul的体系计算得出。

储存条件：+ 4°C (Control lysate -20°C)

## AlphaLISA SureFire Ultra 检测试剂盒

受体酪氨酸激酶		
产品名称	检测点	货号
AlphaLISA Surefire Ultra p-EGFR (Tyr1068) Kit	100	ALSU-PEGFR-A-HV
	500	ALSU-PEGFR-A500
	10,000	ALSU-PEGFR-A10K
	50,000	ALSU-PEGFR-A50K
AlphaLISA Surefire Ultra p-IGF1 Receptor b	100	ALSU-PIGFR-A-HV
	500	ALSU-PIGFR-A500
	10,000	ALSU-PIGFR-A10K
	50,000	ALSU-PIGFR-A50K
AlphaLISA Surefire Ultra p-Insuline Receptor b Kit	100	ALSU-PINR-A-HV
	500	ALSU-PINR-A500
	10,000	ALSU-PINR-A10K
	50,000	ALSU-PINR-A50K
AlphaLISA Surefire Ultra VEGR-Receptor 2 (Tyr1175) Kit	100	ALSU-PVGFR-A-HV
	500	ALSU-PVGFR-A500
	10,000	ALSU-PVGFR-A10K
	50,000	ALSU-PVGFR-A50K

MAPK 通路		
产品名称	检测点	货号
AlphaLISA Surefire Ultra p-ERK (T202/Y204) Kit	100	ALSU-PERK-A-HV
	500	ALSU-PERK-A500
	10,000	ALSU-PERK-A10K
	50,000	ALSU-PERK-A50K
AlphaLISA Surefire Ultra Total ERK1/2 kit	100	ALSU-TERK-A-HV
	500	ALSU-TERK-A500
	10000	ALSU-TERK-A10K
	50000	ALSU-TERK-A50K
AlphaLISA Surefire Ultra p-p38 MAPK (T180/Y182) Kit	100	ALSU-PP38-A-HV
	500	ALSU-PP38-A500
	10000	ALSU-PP38-A10K
	50000	ALSU-PP38-A50K
AlphaLISA Surefire Ultra Total p38 MAPK Kit	100	ALSU-TP38-A-HV
	500	ALSU-TP38-A500
	10000	ALSU-TP38-A10K
	50000	ALSU-TP38-A50K
AlphaLISA Surefire Ultra p-JNK1/2/3 (Thr183/Tyr185)	100	ALSU-PJNK-A-HV
	500	ALSU-PJNK-A500
	10000	ALSU-PJNK-A10K
	50000	ALSU-PJNK-A50K
AlphaLISA Surefire Ultra p-MEK1(Ser218/222)	100	ALSU-PMEK1-A-HV
	500	ALSU-PMEK1-A500
	10000	ALSU-PMEK1-A10K
	50000	ALSU-PMEK1-A50K

其他		
产品名称	检测点	货号
AlphaLISA Surefire Ultra p-CREB (Ser133) Kit	100	ALSU-PCREB-A-HV
	500	ALSU-PCREB-A500
	10,000	ALSU-PCREB-A10K
	50,000	ALSU-PCREB-A50K
AlphaLISA Surefire Ultra Total GAPDH Kit	100	ALSU-TGAPD-A-HV
	500	ALSU-TGAPD-A500
	10,000	ALSU-TGAPD-A10K
	50,000	ALSU-TGAPD-A50K
AlphaLISA Surefire Ultra p-SLP76(Ser376) Kit	100	ALSU-PSLP-A-HV
	500	ALSU-PSLP-A500
	10,000	ALSU-PSLP-A10K
	50,000	ALSU-PSLP-A50K

## AlphaLISA SureFire Ultra MultiPlex (总蛋白及磷酸化蛋白检测)

产品名称	检测点	货号
AKT 通路		
Phospho-ERK1/2(Eu) + Total ERK 545(Tb)	100	MPSU-PTERK-K-HV
	500	MPSU-PTERK-K500
	10,000	MPSU-PTERK-K10K
	50,000	MPSU-PTERK-K50K
MAPK 通路		
p-ERK1/2 (Eu) + Total ERK 545 (Tb) Kit	100	MPSU-PTERK-K-HV
	500	MPSU-PTERK-K-500
	10,000	MPSU-PTERK-K-10K
	50,000	MPSU-PTERK-K-50K
p-ERK (Eu) + p-AKT(Ser473) (Tb) Kit	100	coming soon
	500	coming soon
	10,000	coming soon
	50,000	coming soon
p-ERK (Eu) (directly coupled) *	100	coming soon
	500	coming soon
	10,000	coming soon
	50,000	coming soon

试剂盒组分: Lysis buffer, Activation Buffer, Dilution Buffer, Reaction Buffer (含抗体), Streptavidin Donor beads, CaptSure Acceptor beads, Positive control lysate.

检测点数: 100检测点的每孔检测量是以 1/2 AreaPlate 96孔板中每孔检测总体积为60ul的体系计算; 500、10,000及50,000检测点的每孔检测量是以384孔板中每孔检测总体积为20ul的体系计算得出。

储存条件: + 4°C (Control lysate -20°C)。

## AlphaLISA SureFire Ultra Buffer及Beads

产品名称	检测点	货号
AlphaLISA Surefire Ultra Activation Buffer	10ml	ALSU-AB-10mL
	100ml	ALSU-AB-100mL
AlphaLISA Surefire Ultra Dilution Buffer	10ml	ALSU-DB-10mL
	100ml	ALSU-DB-100mL
AlphaLISA Surefire Ultra Lysis Buffer	10ml	ALSU-LB-10mL
	100ml	ALSU-LB-100mL
AlphaScreen Streptavidin Donor Beads-2mg/mL	60ul	ALSU-ASDB-0.06mL
	1.2ml	ALSU-ASDB-1.2mL
	6ml	ALSU-ASDB-6mL
AlphaLISA CaptSure Acceptor Beads-2mg/mL	60ul	ALSU-ACAB-0.06mL
	1.2ml	ALSU-ACAB-1.2mL
	6ml	ALSU-ACAB-6mL

储存条件: +4°C。

# AlphaLISA SureFire Ultra 试剂盒

## AlphaScreen SureFire (384-well) 及AlphaScreen SureFire HV (96-well) 检测试剂盒

AKT 通路		
产品名称	检测点	货号
AKT1 (Thr308) Kit	500	TGRA2S500
	10,000	TGRA2S10K
	50,000	TGRA2S50K
AKT1/2/3 (Thr308) Kit	500	TGRA3S500
	10,000	TGRA3S10K
	50,000	TGRA3S50K
AKT1/2/3 (Ser473) Kit	100	TGRA4SHV100
	500	TGRA4S500
	10,000	TGRA4S10K
	50,000	TGRA4S50K
AKT1 (Ser473) Kit	500	TGRA5S500
	10,000	TGRA5S10K
	50,000	TGRA5S50K
AKT1/2 (Ser473) Kit	500	TGRAS500
	10,000	TGRAS10K
	50,000	TGRAS50K
Total AKT1 Kit	500	TGRA1S500
	10,000	TGRA1S10K
	50,000	TGRAS50K
Total AKT1/2/3 Kit	100	TGRTAPSHV100
	500	TGRTAPS500
	10,000	TGRTAPS10K
	50,000	TGRTAPS50K
BAD (Ser136) Kit	500	TGRB2S500
	10,000	TGRB2S10K
	50,000	TGRB2S50K
BAD (Ser112) Kit	500	TGRBS500
	10,000	TGRBS10K
	50,000	TGRBS50K
Caspase9 (Ser196) Kit	500	TGRC9S500
	10,000	TGRC9S10K
	50,000	TGRC9S50K
GSK3a (Ser21) Kit	500	TGRGAS500
	10,000	TGRGAS10K
	50,000	TGRGAS50K
GSK3b (Ser9) Kit	500	TGRGBS500
	10,000	TGRGBS10K
	50,000	TGRGBS50K
mTOR (Ser2481) Kit	500	TGRMT2S500
	10,000	TGRMT2S10K
	50,000	TGRMT2S50K
PDK1 (Ser241) Kit	500	TGRPS500
	10,000	TGRPS10K
	50,000	TGRPS50K

MAPK 通路		
产品名称	检测点	货号
ATF1 (Ser63) Kit	500	TGRAF1S500
	10,000	TGRAF1S10K
	50,000	TGRAF1S50K
ATF2 (Thr71) Kit	500	TGRAF2S500
	10,000	TGRAF2S10K
	50,000	TGRAF2S50K
c-Jun (Ser73) kit	500	TGRCJ2S500
	10,000	TGRCJ2S10K
	50,000	TGRCJ2S50K
c-Jun (Ser63) kit	500	TGRCJS500
	10,000	TGRCJS10K
	50,000	TGRCJS50K
ELK1 (Ser383) Kit	500	TGRELK500
	10,000	TGRELK10K
	50,000	TGRELK50K
ERK1/2 (Thr202/Tyr204) kit	100	TGRESBH100
	500	TGRESB500
	10,000	TGRESB10K
	50,000	TGRESB50K
Total ERK kit	100	TGRTESBH100
	500	TGRTESB500
	10,000	TGRTESB10K
	50,000	TGRTESB50K
HSP27 (Ser82) Kit	500	TGRHS27S500
	10,000	TGRHS27S10K
	50,000	TGRHS27S50K
JNK (pan) (Thr183/Tyr185) Kit	500	TGRJPS500
	10,000	TGRJPS10K
	50,000	TGRJPS50K
JNK1/3 (Thr183/Tyr185) Kit	500	TGRJS500
	10,000	TGRJS10K
	50,000	TGRJS50K
Total JNK Kit	500	TGRTJS500
	10,000	TGRTJS10K
	50,000	TGRTJS50K
MKK3/6 (Ser189/207) Kit	500	TGRMK3S500
	10,000	TGRMK3S10K
	50,000	TGRMK3S50K
MKK4 (Ser257/Thr261) Kit	500	TGRMK4S500
	10,000	TGRMK4S10K
	50,000	TGRMK4S50K
MEK1 (Ser217/221) Kit	500	TGRMS500
	10,000	TGRMS10K
	50,000	TGRMS50K
p38 MAPK Kit	500	TGR38S500
	10,000	TGR38S10K
	50,000	TGR38S50K
Total p38 MAPK Kit	500	TGRT38S500
	10,000	TGRT38S10K
	50,000	TGRT38S50K

mTOR 信号/转录调控试剂盒			NFkB 通路		
产品名称	检测点	货号	产品名称	检测点	货号
4E-BP1 (Thr70)	500	TGR4E2S500	NFkB p65 (Ser536) Kit	100	TGRNFSHV100
	10,000	TGR4E2S10K		500	TGRNFS500
	50,000	TGR4E2S50K		10,000	TGRNFS10K
4E-BP1 (Thr37/46)	500	TGR4E5S500		50,000	TGRNFS50K
	10,000	TGR4E5S10K		100	TGRTNFSHV100
	50,000	TGR4E5S50K	Total NFkB p65 Kit	500	TGRTNFS500
p70 S6K (Thr421/424) Kit	500	TGR702S500		10,000	TGRTNFS10K
	10,000	TGR702S10K		50,000	TGRTNFS50K
	50,000	TGR702S50K	IkBa (Ser32/36) Kit	500	TGRIKS500
p70 S6K (Thr229) Kit	500	TGR703S500		10,000	TGRIKS10K
	10,000	TGR703S10K		50,000	TGRIKS50K
	50,000	TGR703S50K	Total IkBa Kit	500	TGRTIKS500
p70 S6K (Thr389) Kit	500	TGR705S500		10,000	TGRTIKS10K
	10,000	TGR705S10K		50,000	TGRTIKS50K
	50,000	TGR705S50K	IKKa (Ser176/180) Kit	500	TGRKAS500
Total p70 S6K Kit	500	TGRT705S500		10,000	TGRKAS10K
	10,000	TGRT705S10K		50,000	TGRKAS50K
	50,000	TGRT705S50K	IKKb (Ser177/181) Kit	500	TGRKBS500
eIF4E (Ser209) Kit	500	TGREIF4S500		10,000	TGRKBS10K
	10,000	TGREIF4S10K		50,000	TGRKBS50K
	50,000	TGREIF4S50K	mTOR (Ser2448) Kit	500	TGRMTS500
mTOR (Ser2448) Kit	500	TGRMTS500		10,000	TGRMTS10K
	10,000	TGRMTS10K		50,000	TGRMTS50K
	50,000	TGRMTS50K	S6 RP (Ser240/244) Kit	500	TGRS6P2S500
S6 RP (Ser240/244) Kit	500	TGRS6P2S500		10,000	TGRS6P2S10K
	10,000	TGRS6P2S10K		50,000	TGRS6P2S50K
	50,000	TGRS6P2S50K	S6 RP (Ser235/236) Kit	500	TGRS6PS500
S6 RP (Ser235/236) Kit	500	TGRS6PS500		10,000	TGRS6PS10K
	10,000	TGRS6PS10K		50,000	TGRS6PS50K
	50,000	TGRS6PS50K	TGFb/SMAD 通路		
产品名称	检测点	货号	产品名称	检测点	货号
SMAD1 (Ser463/465) Kit	500	TGR4SM1S500	STAT1 (Tyr701) Kit	500	TGRS1S500
	10,000	TGR4SM1S10K		10,000	TGRS1S10K
	50,000	TGR4SM1S50K		50,000	TGRS1S50K
SMAD2 (Ser465/467) Kit	500	TGR4SM2S500	STAT3 (Tyr705) Kit	100	TGRS3SHV100
	10,000	TGR4SM2S10K		500	TGRS3S500
	50,000	TGR4SM2S50K		10,000	TGRS3S10K
SMAD3 (Ser423/425) Kit	500	TGR4SM1S500		50,000	TGRS3S50K
	10,000	TGR4SM1S10K	Total STAT3	100	TGRTS3SHV100
	50,000	TGR4SM1S50K		500	TGRTS3S500
STAT5 (Tyr694/699) Kit	500	TGR4SM2S500		10,000	TGRTS3S10K
	10,000	TGR4SM2S10K		50,000	TGRTS3S50K
	50,000	TGR4SM2S50K	STAT5 (Tyr694/699) Kit	500	TGRS5S500
STAT5 (Tyr694/699) Kit	500	TGR4SM1S500		10,000	TGRS5S10K
	10,000	TGR4SM1S10K		50,000	TGRS5S50K
	50,000	TGR4SM1S50K		50,000	TGRS5S50K

# AlphaLISA SureFire Ultra 试剂盒

受体酪氨酸激酶		
产品名称	检测点	货号
ALK (Tyr1604) Kit	500	TGRAL2S500
	10,000	TGRAL2S10K
	50,000	TGRAL2S50K
ALK (Tyr1586) Kit	500	TGRALS500
	10,000	TGRALS10K
	50,000	TGRALS50K
Total ALK Kit	500	TGRTALS500
	10,000	TGRTALS10K
	50,000	TGRTALS50K
c-Met (Tyr1234/1235) Kit	500	TGRCMS500
	10,000	TGRCMS10K
	50,000	TGRCMS50K
EGFR (Tyr1068) Kit	500	TGRERS500
	10,000	TGRERS10K
	50,000	TGRERS50K
ErbB2 (Tyr1221/1222) Kit	500	TGREB2S500
	10,000	TGREB2S10K
	50,000	TGREB2S50K
IGF1 Receptor (Tyr1135/1136) Kit	500	TGRIGFS500
	10,000	TGRIGFS10K
	50,000	TGRIGFS50K
Insuline Receptor (Tyr1150/1151) Kit	500	TGRIRS500
	10,000	TGRIRS10K
	50,000	TGRIRS50K
VEGR-Receptor2 (Tyr1175) Kit	500	TGRVRS500
	10,000	TGRVRS10K
	50,000	TGRVRS50K

其他		
产品名称	检测点	货号
b-catenin (Ser45) Kit	500	TGRBCS500
	10,000	TGRBCS10K
	50,000	TGRBCS50K
CHK1 (Ser345) Kit	500	TGRCHK1S500
	10,000	TGRCHK1S10K
	50,000	TGRCHK1S50K
CHK2 (Thr68) Kit	500	TGRCHK2S500
	10,000	TGRCHK2S10K
	50,000	TGRCHK2S50K
Cofilin (Ser3) Kit	500	TGRCOFS500
	10,000	TGRCOFS10K
	50,000	TGRCOFS50K
CREB (Ser133) Kit	500	TGRCBS500
	10,000	TGRCBS10K
	50,000	TGRCBS50K
GAPDH Kit	500	TGRGDS500
	10,000	TGRGDS10K
	50,000	TGRGDS50K
Histone H3 (Ser10) Kit	500	TGRH3S500
	10,000	TGRH3S10K
	50,000	TGRH3S50K
p53 (Ser15) Kit	100	TGRP53SHV100
	500	TGRP53S500
	10,000	TGRP53S10K
p53 (Ser392) Kit	50,000	TGRP53S50K
	500	TGRP532S500
	10,000	TGRP532S10K
Total p53 Kit	50,000	TGRP532S50K
	100	TGRTP53SHV100
	500	TGRTP53S500
Topoisomerase II a (Thr1343) Kit	10,000	TGRT2AS10K
	50,000	TGRT2AS50K
	500	TGRT2AS500
Topoisomerase II a (Ser1106) Kit	10,000	TGRT2AS10K
	50,000	TGRT2AS50K
	500	TGRT2A2S500
Total Topoisomerase II a Kit	10,000	TGRTT2AS10K
	50,000	TGRTT2AS50K
	500	TGRTT2AS500

试剂盒组分： Lysis buffer, Activation Buffer, Reaction Buffer (含抗体) ,Positive control lysate, Negative control lysate.

检测点数： 100检测点的每孔检测量是以 1/2 AreaPlate 96孔板中每孔检测总体积为50ul的体系计算； 500、10,000及50,000检测点的每孔检测量是以384孔板中每孔检测总体积为11ul的体系计算得出。

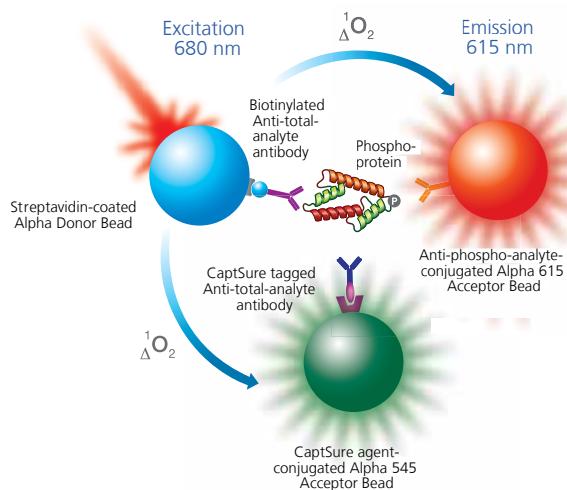
储存条件： + 4°C (Control lysate -20°C)

## AlphaLISA SureFire Ultra 多靶点检测试剂盒

新品

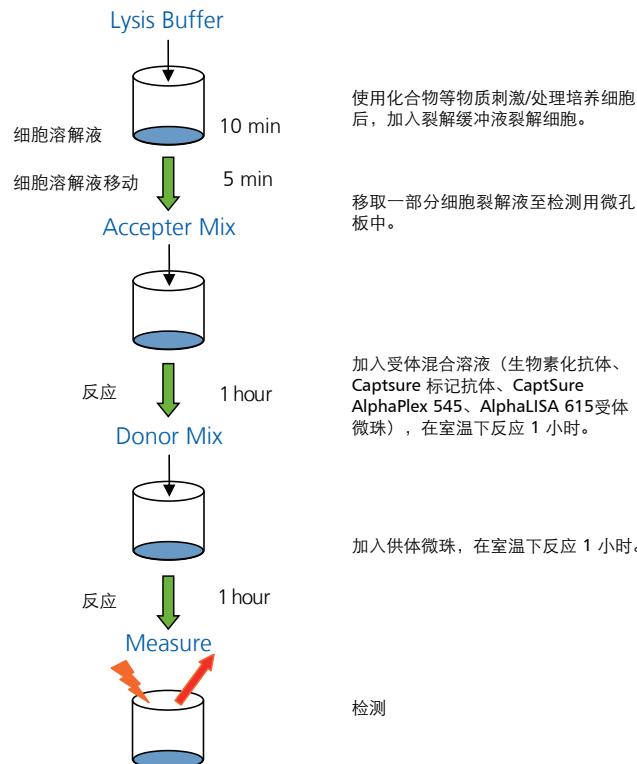
Alpha SureFire Ultra 多靶点检测试剂盒利用 2 种不同波长的 Alpha 受体微珠、AlphaLISA 615(Eu) 和 AlphaPlex 545(Tb)，可同时检测细胞裂解液中的内源性磷酸化蛋白质和总蛋白质。

### SureFire Ultra 多靶点检测的原理

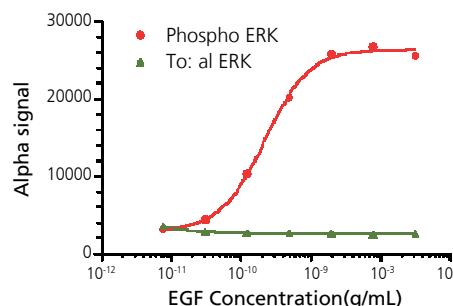


细胞裂解液中的目标蛋白质结合了 CaptSure 标记抗体和 AlphaPlex 545 受体，其磷酸化部位还结合了抗磷酸化抗体 AlphaLISA 受体微珠。此外，目标蛋白质还结合了生物素化抗体与链霉亲合素供体微珠。

采用激发光(680nm)照射后，AlphaLISA受体微珠(615nm)和 AlphaPlex 545 受体微珠 (545nm) 同时发射光信号以供检测。



### SureFire Ultra 多靶点检测示例



### 同时检测磷酸化 ERK 和总 ERK

将A431细胞接种到微孔板（40,000个细胞/孔，培养基为MEM+10%FCS）上，孵育过夜。进行2小时的血清饥饿处理后，用EGF刺激10分钟。加入SureFire Ultra裂解缓冲液（100 μL/孔）裂解细胞，取细胞裂解液10μL通过SureFire Ultra多靶点检测方法同时检测 p-ERK 和总 ERK。

产品名称	数量 (检测点)	产品货号
<b>AKT 信号通路</b>		
Alpha SureFire® Ultra™ Multiplex	100	MPSU-PTAKT-K-HV
p-Akt 1/2/3 (S473) (Eu) + Total Akt1 (Tb) Kit	500	MPSU-PTAKT-K-500
	10,000	MPSU-PTAKT-K-10K
	50,000	MPSU-PTAKT-K-50K
<b>MAPK 信号通路</b>		
Alpha SureFire® Ultra™ Multiplex	100	MPSU-PTERK-K-HV
p-ERK 1/2 (Eu) + Total ERK 545 (Tb) Kit	500	MPSU-PTERK-K-500
	10,000	MPSU-PTERK-K-10K
	50,000	MPSU-PTERK-K-50K

成分：Alpha多靶点反应缓冲液 1 (含生物素化抗体)、Alpha多靶点反应缓冲液 2 (含CaptSure标记抗体)、Alpha 615抗磷酸化抗体受体微珠、Alpha 545 CaptSure 受体微珠、链霉亲合素供体微珠、激活缓冲液、稀释缓冲液、裂解缓冲液、阳性对照裂解液。

数量：100 检测点试剂盒中的 1 检测点相当于按试剂盒推荐量 (100 μL) 使用时，96 孔微孔板中 1 个孔的用量，500/10,000 检测点试剂盒中的 1 检测点相当于按试剂盒推荐量 (40 μL) 使用时，1/2 AreaPlate-96 孔板中 1 个孔的用量。

产品名称	数量	检测点 (96/384 微孔板)
SureFire Ultra	60 μL	200/600
Terbium Captsure	1.2 mL	4,000/12,000
Acceptor Beads,2mg/mL	6 mL	20,000/60,000

成分：Alpha 545 CaptSure 受体微珠。

数量：60 μL 的 Alpha 545 Captsure 供体微珠相当于使用 96 孔微孔板检测时 200 个孔的用量或使用 384 孔微孔板检测时 600 个检测点的用量。

### 产品价格

产品名称	数量 (检测点)	产品货号
Alpha SureFire Ultra MultiPlex kit	100	120,000
	500	400,000
	10,000	on request
	50,000	on request
SureFire Ultra Terbium CaptSure Acceptor Beads	60 μL	39,000
	1.2 mL	390,000
	6 mL	on request

# AlphaScreen PhosphoSensor 及 Tyrosine Kinase

## 无需磷酸化抗体的检测方法：AlphaScreen PhosphoSensor

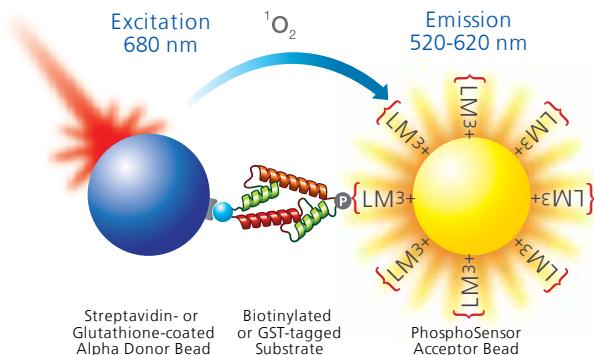
AlphaScreen PhosphoSensor 受体微珠上包被有 Lewis 金属离子用来捕获磷酸化基团 ( $LM3^+$ )。当生物素化底物被激酶磷酸化后，PhosphoSensor 受体微珠结合磷酸根基团，同时链霉亲和素 (SA) 包被的供体微珠结合生物素，此时可以检测到有效信号。如果您没有特异性的磷酸化抗体，此试剂盒是很好的选择。

产品名称	检测点	货号
AlphaScreen PhosphoSensor Assay kit	1,000	6760307D
	10,000	6760307M

试剂盒组分：phosphosensor受体微珠，链霉亲和素供体微珠，生物素标记肽，缓冲液。

检测点数：以384孔板25ul总反应体系的反应为一个检测点的计算量。

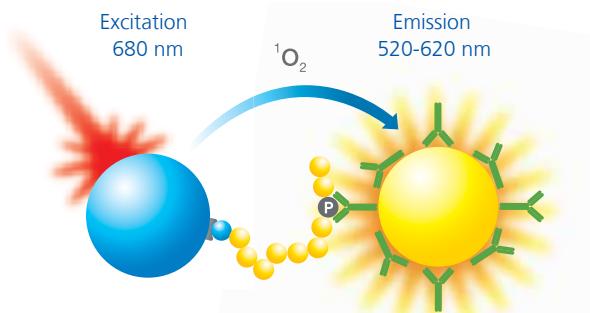
储存条件：+ 4°C。



## 需要使用抗体的检测方法

### 酪氨酸激酶磷酸化检测试剂盒

AlphaScreen 磷酸化酪胺检测试剂盒 (PY20, PT66, P-Tyr-100) 在酪氨酸激酶的激活下可以检测蛋白或多肽底物的磷酸化水平。



产品名称	检测点	货号
Phosphotyrosine (PY20) Assay kit	500	6760601C
	10,000	6760601M
Phosphotyrosine (PT66) Assay kit	500	6760602C
	10,000	6760602M
Phosphotyrosine (P-Tyr-100) Assay kit	500	6760620C
	10,000	6760620M

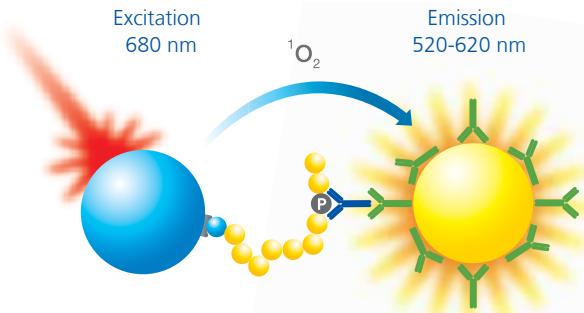
试剂盒组分：链霉亲和素包被供体微珠，抗磷酸化酪胺抗体包被的受体微珠，生物素化抗磷酸化酪胺多肽，对照缓冲液。

检测点数：以384孔板25ul总反应体系的反应为一个检测点的计算量。

储存条件：+ 4°C。

### 丝氨酸/苏氨酸激酶磷酸化检测试剂盒

丝氨酸/苏氨酸激酶活性的检测可以使我们市售的以下试剂：抗丝氨酸/苏氨酸特异性抗体及 AlphaScreen IgG 检测试剂盒 (Protein A或种属特异性二抗)。



产品名称	检测点	货号
Mouse IgG detection kit	500	6760606M
	10,000	6760606R
Rabbit IgG detection kit	500	6760607M
	10,000	6760607R
Protein A IgG detection kit	500	6760617M
	10,000	6760617R

试剂盒组分：链霉亲和素包被供体微珠，二抗包被或Protein A包被的受体微珠，生物素化抗体，对照缓冲液。

检测点数：以384孔板25ul总反应体系的反应为一个检测点的计算量。

储存条件：+ 4°C。

## cAMP 检测

### cAMP 检测试剂盒

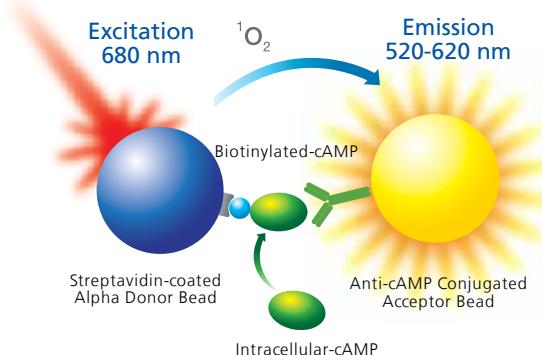
AlphaScreen cAMP检测是基于细胞产生的游离cAMP及生物素化cAMP探针竞争的检测原理。生物素化的cAMP会将供体微珠和受体微珠之间的距离拉近从而产生一个特异性信号。当细胞内源性cAMP含量增加时，细胞内的cAMP会替代生物素化cAMP探针与供体微珠相结合，产生的Alpha信号会随之降低，信号降低的量与细胞内cAMP的量成一定比例。

产品名称	检测点	货号
AlphaScreen cAMP assay kit	1,000	6760625D
	10,000	6760625M
	50,000	6760625R
cAMP biotin supplement	10,000	6760301M
	50,000	6760301R
	100	AL315HV
AlphaLISA cAMP assay kit	500	AL315C
	5,000	AL315F

试剂盒组分：链霉亲和素供体微珠，抗cAMP抗体标记受体微珠，生物素标记cAMP，cAMP标准品，缓冲液。

检测点数：以384孔板25 ul总反应体系的反应为一个检测点的计算量。

储存条件：+ 4°C。



## 磷酸化AKT、ERK检测

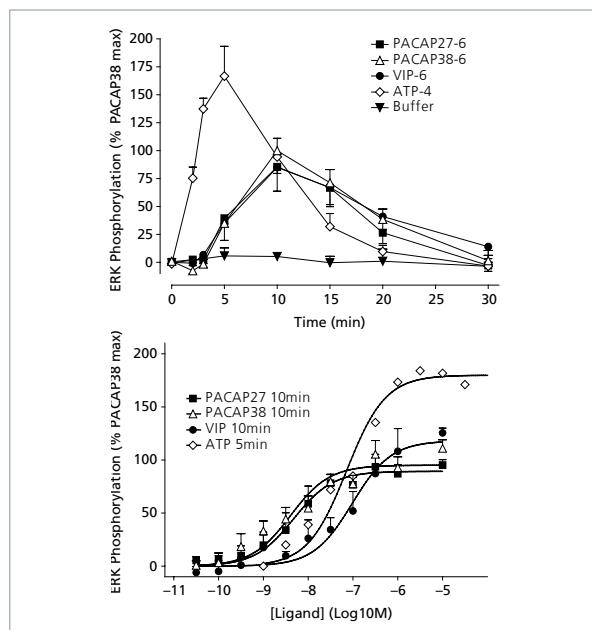
### AlphaScreen Surefire AKT、ERK 检测试剂盒

Gq、Gi协同作用的GPCR靶点通常很难用cAMP或钙流的方法检测。对于这类靶点，我们会选择ERK盒AKT来进行检测。具体检测原理及产品信息，请参见p16~p22页。

### 针对小鼠神经干细胞(mNSC) 进行AlphaScreen Surefire 检测示例 (右图)

利用AlphaScreen Surefire ERK试剂盒检测PAC (Pituitary adenylate cyclase activating peptide【PACAP】receptor1) 信号通路。右侧上图展示了在不同配体作用下随着时间推移ERK磷酸化水平的变化。右侧下图展示了随着不同配体浓度的变化ERK磷酸化水平的变化。将不同配体加入mNSC细胞培养液中，培养基移除后，加入Surefire裂解液，随后遵照试剂盒说明书进行磷酸化ERK检测。

Ref; Lodge AP, Langmead CJ, Daniel G, Anderson GW, Werry TD., Performance of mouse neural stem cells as a screening reagent: characterization of PAC1 activity in medium-throughput functional assays. J Biomol Screen. 2010 Feb;15(2):159-68. Neurosciences CED, GlaxoSmithKline Research & Development, Harlow, Essex, UK.



## cGMP 检测

### cGMP 检测试剂盒

此试剂盒是用来检测样品中cGMP含量的。它可以检测治疗性靶点，如磷酸二酯酶（phosphodiesterase）、昔酸环化酶（guanylate cyclase）、一氧化氮合成等等。检测是以竞争法检测为基础：生物素化cGMP与样品中游离的cGMP竞争特异性抗cGMP抗体位点。使用此试剂需要同时购买Protein A (6760617) 检测试剂盒。

产品名称	检测点	货号
Biotinylated cGMP Supplement with Antibody kit	10,000	6760308M
	50,000	6760308R
	500	6760617C
AlphaScreen Protein A kit	10,000	6760617M
	50,000	6760617R
AlphaLISA cGMP assay kit	100	AL316HV
	500	AL316C
	5,000	AL316F

试剂盒组分：Biotinylated cGMP Supplement with Antibody kit: 生物素化cGMP，抗cGMP抗体。

Protein A kit: 链霉亲和素供体微珠，Protein A包被受体微珠。

检测点数：以384孔板25ul总反应体系的反应为一个检测点的计算量。

储存条件：+ 4°C。

# Alpha Biologics: 生物药物相关应用

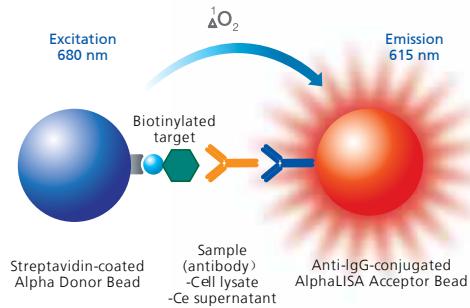
## Biologics

Alpha技术——为抗体药物研发提供可靠性和重现性更好的数据和结果

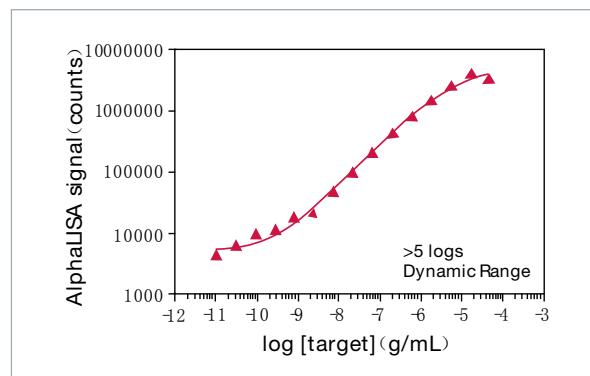
### 基于细胞的筛选

Alpha技术具有宽泛的线性范围，可满足对样品进一步稀释的更高要求。

#### 杂交瘤筛选 (vs ELISA)



Vainshtein I, Kurose S, Vickroy J, Russell Grove J, Liang M. Fast and efficient soluble antigen-antibody screening using homogeneous AlphaScreen assay. Poster Presentation. SBS 11th Annual Conference. September 2004, Orlando (USA)



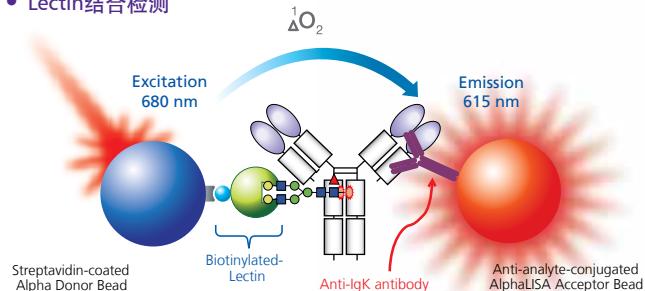
#### CHO细胞产生的抗体量检测

##### • AlphaLISA免疫检测试剂盒

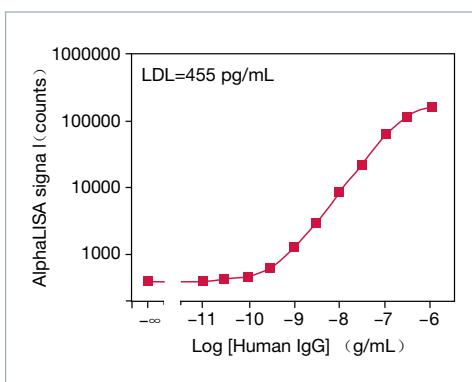
描述	数量 (well)	货号	检测下限	
			in Buffer	in Serum
Human IgG Immunoassay kit	500	AL205C	0.24ng/mL	N/A
	5,000	AL205F		

产品特点等方面的信息，请参阅11页。

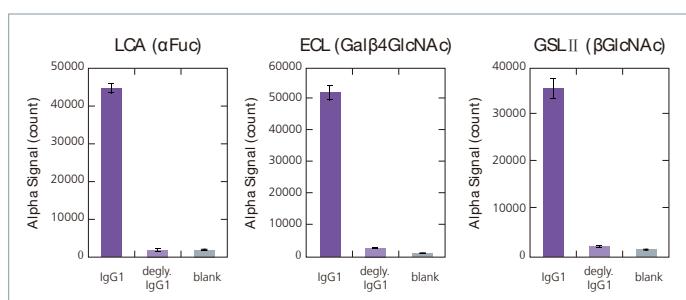
##### • Lectin结合检测



M Onitsuka, T Omasa. Rapid evaluation of N-glycosylation status of antibodies with chemiluminescent lectin-binding assay. Journal of Bioscience and Bioengineering. 2014 Dec.



Human IgG免疫检测试剂盒标准曲线



使用生物素化凝集素及完整的IgG1来检测去糖基化的IgG1。

## 同型筛选

### AlphaLISA 免疫检测试剂盒

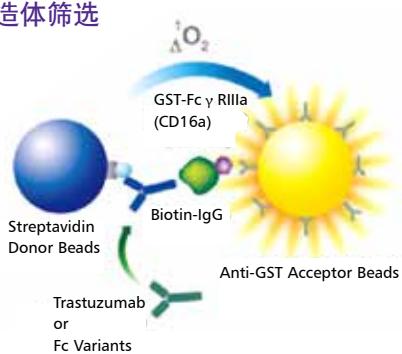
产品名称	数量 (well)	货号	检测下限	检测上限
Human IgG1 (isotyping) immunoassay kit	100	AL307HV	47 pg/mL	1,000,000 pg/mL
	500	AL307C		
	5,000	AL307F		
Human IgG2 (isotyping) immunoassay kit	100	AL308HV	—	—
	500	AL308C		
	5,000	AL308F		
Human IgG3 (isotyping) immunoassay kit	100	AL309HV	500 pg/mL	3,000,000 pg/mL
	500	AL309C		
	5,000	AL309F		
Human IgG4 (isotyping) Immunoassay kit	100	AL310HV	300 pg/mL	500,000 pg/mL
	500	AL310C		
	5,000	AL310F		
Human IgA Immunoassay kit	500	AL262C	449 pg/mL	3,000,000 pg/mL
	5,000	AL262F		
Human IgE Immunoassay kit	500	AL292C	6.4 pg/mL	3,000,000 pg/mL
	5,000	AL292F		
Human IgM Immunoassay kit	500	AL263C	511 pg/mL	3,000,000 pg/mL
	5,000	AL263F		
Mouse IgG1 (isotyping) Immunoassay kit	100	AL522HV	117 pg/mL	3,000,000 pg/mL
	500	AL522C		
	5,000	AL522F		

产品名称	数量 (well)	货号	检测下限	检测上限
Mouse IgG2a (isotyping) Immunoassay kit	100	AL523HV	23 pg/mL	1,000,000 pg/mL
	500	AL523C		
	5,000	AL523F		
Mouse IgG2b (isotyping) immunoassay kit	100	AL524HV	60 pg/mL	1,000,000 pg/mL
	500	AL524C		
	5,000	AL524F		
Mouse IgG3 (isotyping) immunoassay kit	100	AL525HV	60 pg/mL	1,000,000 pg/mL
	500	AL525C		
	5,000	AL525F		
Mouse IgA (isotyping) Immunoassay kit	100	AL528HV	00 pg/mL	1,000,000 pg/mL
	500	AL528C		
	5,000	AL528F		
Mouse IgE (isotyping) Immunoassay kit	100	AL526HV	40 pg/mL	1,000,000 pg/mL
	500	AL526C		
	5,000	AL526F		
Mouse IgM (isotyping) Immunoassay kit	100	AL527HV	104 pg/mL	1,000,000 pg/mL
	500	AL527C		
	5,000	AL527F		

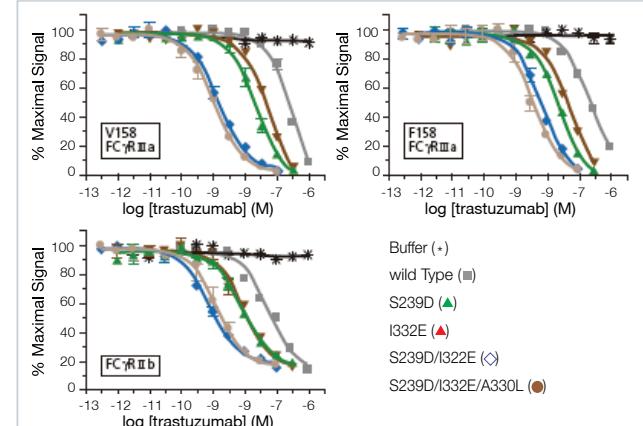
产品特点等方面的信息，请参阅5~13页。

## 抗体工程

### Fc区域改造体筛选



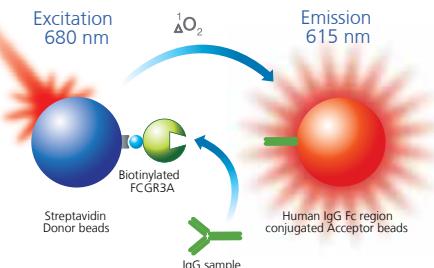
Lazar GA, Dang W, Karki S, Vafa O, Peng JS, Hyun L, Chan C, Chung HS, Eivazi A, Yoder SC, Vielmetter J, Carmichael DF, Hayes RJ, Dahiyat BI. Engineered antibody Fc variants with enhanced effector function. Proc Natl Acad Sci U S A. 2006 Mar 14; 103 (11) : 4005-1 Xencor-USA



曲妥珠单抗的Fc区域改造体亲和力结合试验：FcγRIIa (V158) 的变种 (左上)，FcγRIIa (F158) (右上)，FcγRIIa (V158) (下)。

### Fcγ受体结合检测试剂盒

该试剂盒基于人IgG Fc区域与Fc受体之间的结合，并在此基础上采用竞争法检测来鉴定抗体与Fc受体的亲和力，适用于诸如高ADCC活性的抗体筛选。



生物素化的Fc受体 (FCGR3A) 与链霉亲和素包被的供体微珠相结合，其与偶联了血清衍生人IgG (Fc区) 的受体微珠结合并产生发光信号。当用于评价IgG样品亲和力时，若其亲和力更强，便会与IgG (Fc区) 偶联受体微珠竞争与Fc受体 (FCGR3A) 结合并抑制信号，信号将随着IgG样品的量增加而降低。

### • AlphaLISA 免疫检测试剂盒

产品名称	数量 (well)	货号
AlphaLISA FCGR3A (176Val) Binding Kit	100	AL347HV
	500	AL347C
	5,000	AL347F
Biotinylated Human FCGR3A (176Phe)	1.5 µg	AL347S
	100	AL348HV
	500	AL348C
AlphaLISA FCGR3A (176Val) Binding Kit	5,000	AL348F
	1.5 µg	AL348S

数量：1个检测点相当于按试剂盒推荐量 (40 µL) 使用时，384孔板或96孔板中1个孔的用量。

产品特点等方面的信息，请参阅5~13页。

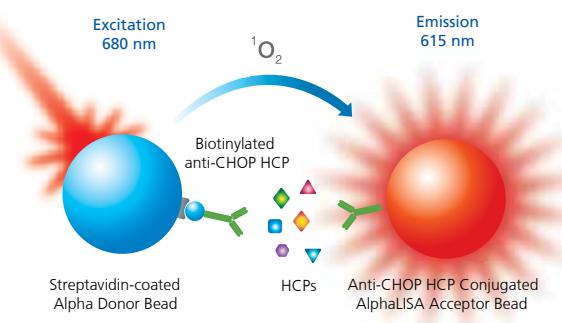
# Alpha Biologics: 生物药物相关应用

## 工艺开发

### AlphaLISA免疫检测试剂盒

产品名称	检测点	货号	检测下限	检测上限
CHO cell protein (CHO-P)	500 5,000	AL210C AL210F	181 pg/mL 130 pg/mL	300,000 pg/mL 1,000,000 pg/mL
Immunoassay kit (Broad Reactivity)	500 5,000	AL301C AL301F		
PER.C6 HCP Immunoassay kit	500 5,000	AL302C AL302F	1486 pg/mL 10.8 pg/mL	1,000,000 pg/mL 300,000 pg/mL
Residual Protein A Immunoassay kit	500 5,000	AL287C AL287F		
Host cell residual DNA Immunoassay kit	500 5,000	AL331C AL331F	0.06 ng/mL	300,000 pg/mL

### CHO Host Cell Protein Detection Assay



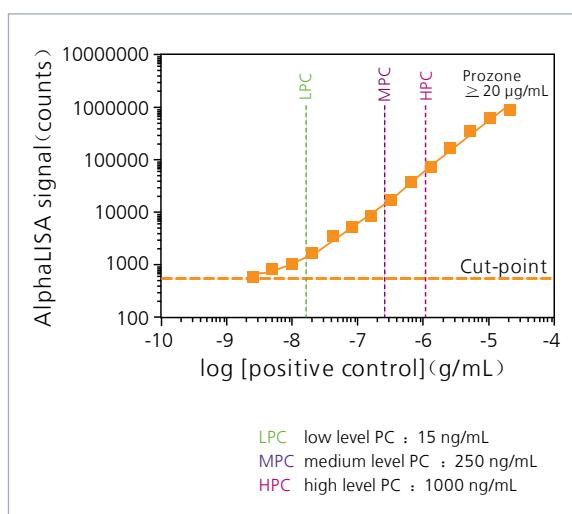
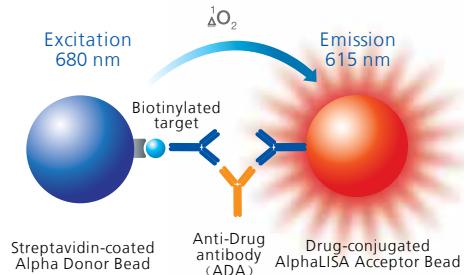
## PK / PD 免疫原性

### AlphaLISA免疫检测试剂盒

产品名称	数量 (well)	货号	檢出下限 in Buffer	檢出下限 in Serum	檢出上限	人	小鼠	大鼠	牛	猪	狗	猩猩	猴
Human IgG4 (PK/PD study) Immunoassay	500 5000	AL304C AL304F	18 pg/mL		35 pg/mL	300,000 pg/mL	0	×	×				
Anti-PEG IgM Immunoassay kit	100 500 5000	AL313HV AL313C AL313F	6.7 ng/mL			100,000 pg/mL							

### ADA检测

- 高灵敏度符合FDA的指导原则 (250-500 ng/mL)
- 卓越的重现性 ( $\leq 10\% \text{CV}$ )
- 最小的基质效应 (血清成分的干扰)
- +酸处理样品可耐受药物抗体达100  $\mu\text{g/mL}$ ,  
-酸处理为10  $\mu\text{g/mL}$
- 宽泛动态范围

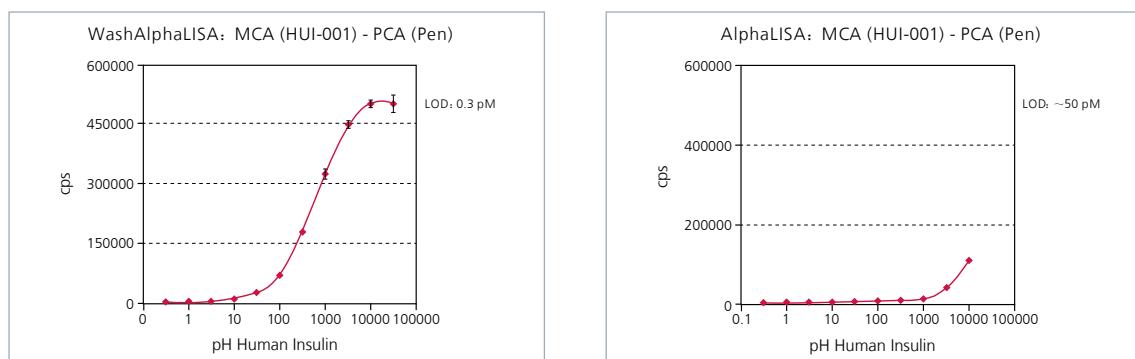
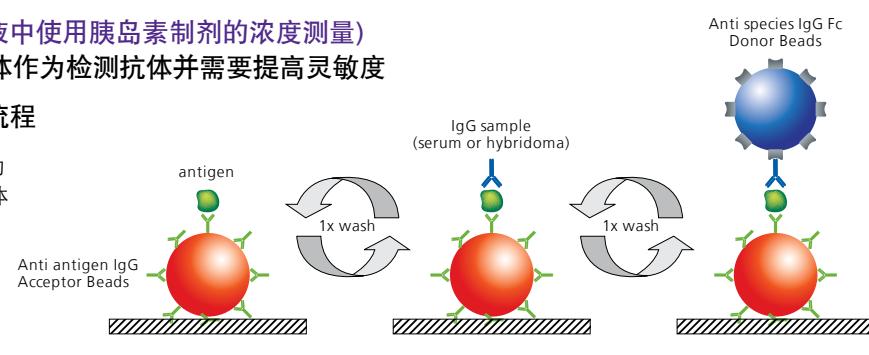


**Tips: Semi-Wash Alpha (血液中使用胰岛素制剂的浓度测量)**  
当使用低纯度（特异性）抗体作为检测抗体并需要提高灵敏度

● 受体微珠包被的制备流程

使用high binding高结合力  
微孔板预涂受体微珠受体

↓  
离心, 20度放置1晚  
↓  
缓冲液洗涤1次



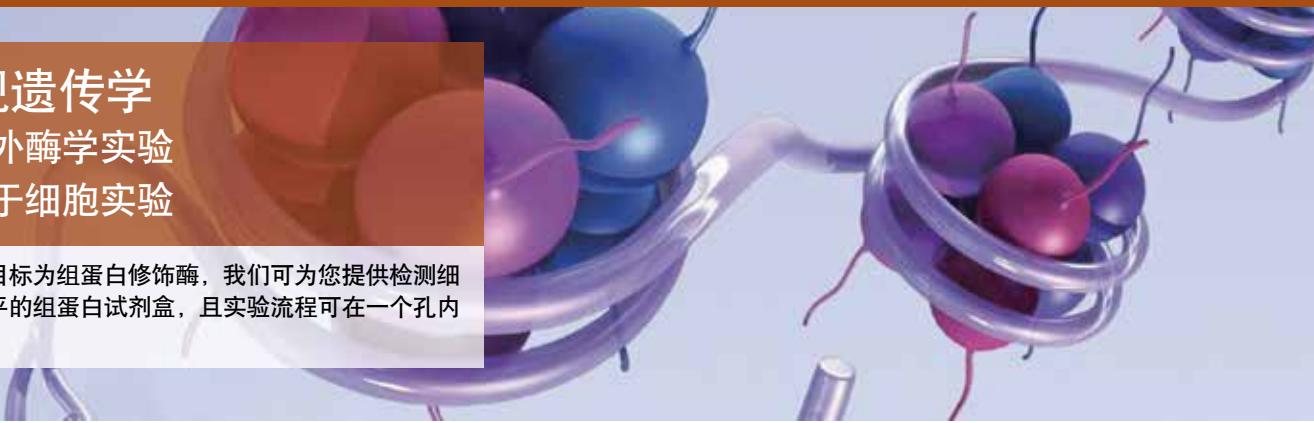
结论：使用未纯化多克隆抗体检测胰岛素（PCA-Pen）。左图为使用Alpha半洗涤步骤，右图为使用直接混合检测。结果显示当使用半洗涤步骤时，即便使用纯度低的多克隆抗体也能够高检测灵敏度。

Ref: Fritz Poulsen, Wash-LOCI-A Semi-Heterogeneous Version of the LOCI Technology Allowing Removal of Unbound Material After Each Assay Step., Chapter 16, Trends in Immunolabelled and Related Techniques.

## 表观遗传学

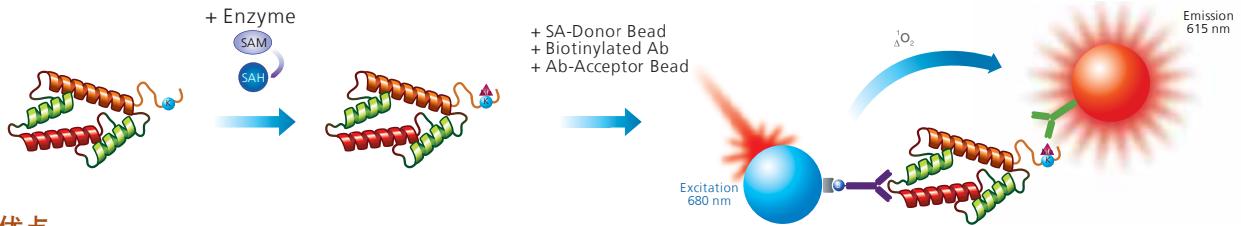
- 体外酶学实验
- 基于细胞实验

筛选的目标为组蛋白修饰酶，我们可为您提供检测细胞学水平的组蛋白试剂盒，且实验流程可在同一个孔内完成。



## 体外酶学实验

在被特定的酶修饰过的组蛋白H3或p53中，加入生物素化的抗体和链霉亲和素包被的供体微珠，受体微珠上的抗体识别修饰过的氨基酸区域，形成三明治夹心，组蛋白夹在中间，使供体微珠与受体微珠相互靠近，产生反应并检测出信号。也可使用生物素化多肽作为底物。还可通过使用已甲基化或乙酰化的底物，进行脱甲基酶或脱乙酰基酶活性的检测。



### 优点

- 简单操作 ..... 均相法测定，不需要洗板和分板。
- 高特异性 ..... 使用的抗体具有组蛋白修饰基团的高特异性。
- 可直接使用 ..... 提供以下酶的测定。

### Ready to Use

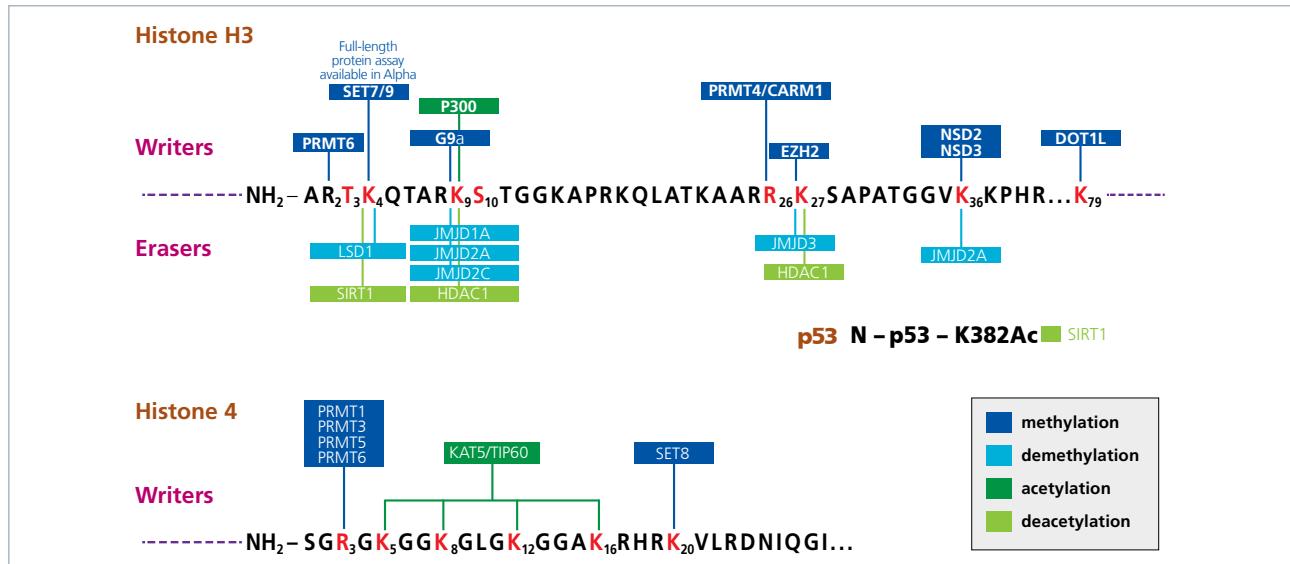
<http://www.perkinelmer.com/category/alpha-reagents/epigenetics/>

对于下列酶的测定，我们提供相应的技术说明。



! PerkinElmer 公司并不出售酶和底物。我们会提供酶以及一些必要试剂的供应商、抗体特异性数据等，详情请参阅技术说明和产品数据表。

产品数据表，可以从[www.perkinelmer.com/tools/coasearch](http://www.perkinelmer.com/tools/coasearch)下载。



## 受体微珠

甲基化修饰									
Mark	抗体特异性			Ready to Use		Anti-mark 受体微珠			
	me1	me2	me3	Methyl-transferase	Demethylase	货号	规格	检测点	
H3K4 unmodified	×	×	×	LSD1		AL119C	250 µg	500	
						AL119M	5 mg	10,000	
						AL119R	25 mg	50,000	
H3K4me1-2	○	○	×	SET7/9		AL116C	250 µg	500	
						AL116M	5 mg	10,000	
						AL116R	25 mg	50,000	
H3K9/K27 unmodified	×	×	×	JMJD1A		AL138C	250 µg	500	
						AL138M	5 mg	10,000	
						AL138R	25 mg	50,000	
H3K9me2	×	○	×	G9a	JMJD2A	AL117C	250 µg	500	
						AL117M	5 mg	10,000	
						AL117R	25 mg	50,000	
H3K27me2-1	○	○	×	EZH2	JMJD3	AL121C	250 µg	500	
						AL121M	5 mg	10,000	
						AL121R	25 mg	50,000	
H3K27me3	×	○	○			AL122C	250 µg	500	
						AL122M	5 mg	10,000	
						AL122R	25 mg	50,000	
H3K36me2	○	○	×	JMJD2A		AL123C	250 µg	500	
						AL123M	5 mg	10,000	
						AL123R	25 mg	50,000	
H3K36me2 (Nucleosome)	×	○	△	NSD2 NSD3		AL152C	250 µg	500	
						AL152M	5 mg	10,000	
						AL152R	25 mg	50,000	
H3R2me2	○	○	×	PRMT6		AL139C	250 µg	500	
						AL139M	5 mg	10,000	
						AL139R	25 mg	50,000	
H3 protein (C-ter)				DOT1L		AL147C	250 µg	500	
						AL147M	5 mg	10,000	
						AL147R	25 mg	50,000	
H3/H4R me generic	○	○	-	PRMT4/CARM1		AL151C	250 µg	500	
						AL151M	5 mg	10,000	
						AL151R	25 mg	50,000	
H4K20me1	○	×	×	SET8		AL145C	250 µg	500	
						AL145M	5 mg	10,000	
						AL145R	25 mg	50,000	
H4R3me		○ (2a) △ (2s)		PRMT1 PRMT3 PRMT5 PRMT6		AL150C	250 µg	500	
						AL150M	5 mg	10,000	
	△					AL150R	25 mg	50,000	
O-GlcNAc 修饰									
Mark	抗体特异性			Ready to Use		货号	规格	检测点	
S400-O-GlcNAc	× (S1011-O-GlcNAc)			OGA (N-acetylglucosaminidase)		AL144C	250 µg	500	
						AL144M	5 mg	10,000	
						AL144R	25 mg	50,000	

数量：在384孔板中进行测定（反应体系为25 µL）所得，250 µg的受体微珠可测500个点。

保存：+4°C。

# AlphaLISA Epigenetics

## 受体微珠

乙酰化修饰						
Mark	Ready to Use		Anti-mark Acceptor Beads			检测点
	Acetyl-transferase	Deacetylase	货号	规格		
H3K4 unmodified		SIRT1	AL119C	250 µg	500	
			AL119M	5 mg	10,000	
			AL119R	25 mg	50,000	
			AL138C	250 µg	500	
H3K9/K27 unmodified		HDAC1	AL138M	5 mg	10,000	
			AL138R	25 mg	50,000	
			AL114C	250 µg	500	
			AL114M	5 mg	10,000	
H3K9ac	p300		AL114R	25 mg	50,000	
			AL120C	250 µg	500	
			AL120M	5 mg	10,000	
			AL120R	25 mg	50,000	
H3K27ac		HDAC1	AL143C	250 µg	500	
			AL143M	5 mg	10,000	
			AL143R	25 mg	50,000	
			Non-Histone Target			
p53 K382ac		SIRT1	AL124C	250 µg	500	
			AL124M	5 mg	10,000	
			AL124R	25 mg	50,000	

数量：在384孔板中进行测定（反应体积为25 µL）所得，250 µg的受体微珠可测500个点。

保存：+ 4°C

## 生物素化的抗组蛋白抗体

甲基化修饰								
Mark	抗体特异性			Ready to Use		Anti-mark 受体微珠		
	me1	me2	me3	Methyl-transferase	Demethylase			
H3 protein (C-ter)				SET7/9		AL118C	2 µg	500
				NSD2		AL118M	40 µg	10,000
				NSD3		AL118R	200 µg	50,000
H4 protein (C-ter)				SET8		AL146C	2 µg	500
						AL146M	40 µg	10,000
						AL146R	200 µg	50,000
H3K79me2	△	○	○	DOT1L		AL148C	2 µg	500
						AL148M	40 µg	10,000
						AL148R	200 µg	50,000

数量：在384孔板中进行测定（反应体积为25 µL）所得，250 µg的受体微珠可测500个点。

保存：+ 4°C

## 供体微珠

产品名称	货号	规格	检测点
Streptavidin-coated Donor beads	6760002S	1 mg	2,000
	6760002	5 mg	10,000
	6760002B	50 mg	100,000

数量：在384孔板中进行测定（反应体系25 µL）所得，1 mg的微珠可测2,000个点。

保存：+ 4°C。

## 组蛋白检测试剂盒

产品名称	货号	检测点
Total Histone H4 detection kit	AL726HV	100
	AL726C	500
	AL726F	5,000

## 反应缓冲液

产品名称	货号	规格	检测点
AlphaLISA	AL008C	10 mL	20,000
5X Epigenetics Buffer 1	AL008F	100 mL	200,000

包含：AlphaLISA 5X Epigenetics Buffer1, AlphaLISA 30X Epigenetics Buffer Supplement。

保存：+ 4°C。

## AlphaPlex 545受体微珠

产品名称	货号	规格	检测点
Anti-unmodified H3K4	AP119TB-C	250 µg	500
AlphaPlex 545	AP119TB-M	5 mg	10,000
Acceptor Beads	AP119TB-R	25 mg	50,000

## 基于细胞的实验

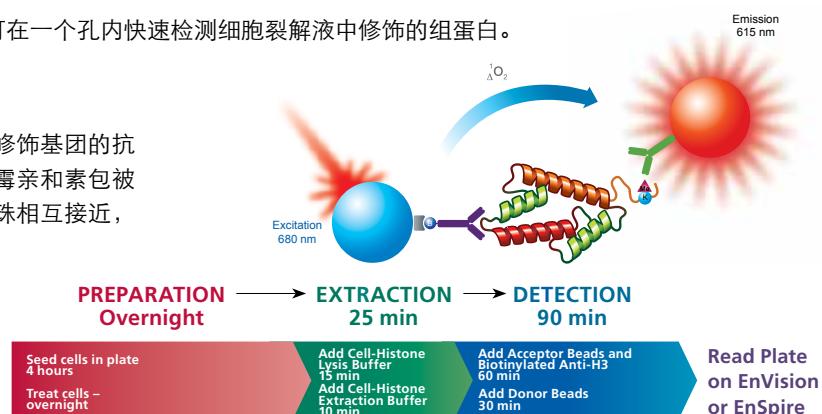
可检测细胞内组蛋白修饰的试剂盒。无需洗涤，可在同一个孔内快速检测细胞裂解液中修饰的组蛋白。

### 原理

细胞裂解液中含有被修饰的组蛋白，依次加入抗修饰基团的抗体标记的受体微珠，生物素化抗组蛋白抗体与链霉亲和素包被的供体微珠。修饰的组蛋白使供体微珠和受体微珠相互接近，产生信号并检测。

### 实验操作步骤

全部实验步骤可在同一个孔内完成。



PREPARATION  
Overnight → EXTRACTION  
25 min → DETECTION  
90 min



## 细胞学水平检测试剂盒

Mark	产品名称	检测点	货号
H3K9ac	AlphaLISA H3K9ac Cellular Detection kit	100	AL714HV
		500	AL714C
		5,000	AL714F
		100	AL716HV
H3K4me2	AlphaLISA H3K4me2 Cellular Detection kit	500	AL716C
		5,000	AL716F
		100	AL717HV
		500	AL717C
H3K9me2	AlphaLISA H3K9me2 Cellular Detection kit	5,000	AL717F
		100	AL727HV
		500	AL727C
		5,000	AL727F
Unmodified H3K4	AlphaLISA unmodified Cellular Detection kit	100	AL719HV
		500	AL719C
		5,000	AL719F
		100	AL720HV
H3K27ac	AlphaLISA H3K27ac Cellular Detection kit	500	AL720C
		5,000	AL720F
		100	AL721HV
		500	AL721C
H3K27me2-1	AlphaLISA H3K27me2-1 Cellular Detection kit	5,000	AL721F
		100	AL722HV
		500	AL722C
		5,000	AL722F
H3K36me	AlphaLISA H3K36me Cellular Detection kit	100	AL723HV
		500	AL723C
		5,000	AL723F
		100	AL748HV
H3K79me2	AlphaLISA H3K79me2 Cellular Detection kit	500	AL748C
		5,000	AL748F
		100	AL724HV
		500	AL724C
H2A-K119ub	AlphaLISA H2A-K119ub Cellular Detection kit	5,000	AL724F
		100	AL725HV
		500	AL725C
		5,000	AL725F
Total H2A	AlphaLISA Total H2A Cellular Detection kit	100	AL729C
		500	AL729F
Total H4	AlphaLISA Total Histone H4 Cellular Detection kit	500 新品	AL729F

数量：在1/2 AreaPlate-96孔板中使用推荐的体积(50 $\mu\text{L}$ )时，可使用的总量。

保存：+ 4°C

### 试剂盒组分

- 抗修饰基团的组蛋白H3的抗体标记的受体微珠
- 链霉亲和素包被的供体
- 生物素化抗组蛋白H3（C-末端）抗体
- Cell-Histone Lysis buffer
- Cell-Histone Extraction buffer
- Cell-Histone Detection (10X) buffer

- ① 其他实验中所需要的试剂以及实验步骤、抗体特异性等数据，请参考产品数据表。  
[www.perkinelmer.com/tools/coasearch](http://www.perkinelmer.com/tools/coasearch)



## AlphaLISA Cell-Histone Buffer Set

AlphaLISA细胞学水平的组蛋白检测试剂盒内已包含缓冲液，您也可以单独购买。

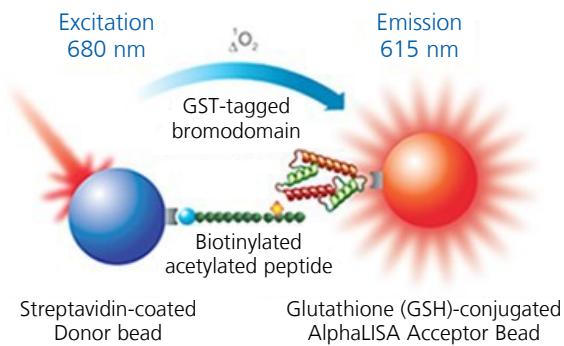
产品名称	检测点	货号
AlphaLISA	500	AL009C
Cell-Histone Buffer Set	5,000	AL009F

保存：+ 4°C

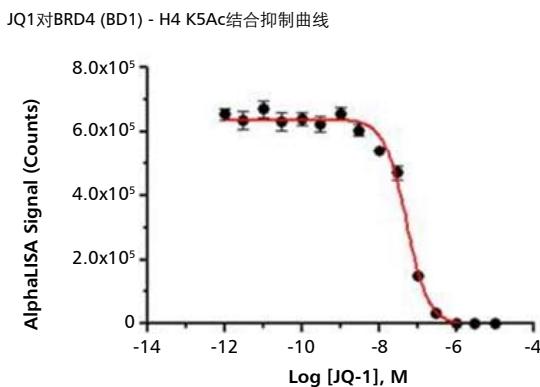
# AlphaLISA Epigenetics

## AlphaLISA Bromodomain蛋白结合抑制实验

Bromodomain (BRDs) 被认为是乙酰化组蛋白残基的“Reader”，与多种疾病相关。BRD4是BET家族（BD1,BD2）的一员，因其是重要的药物研发靶点而受到关注。我们的试剂盒可以用来检测BRD4与乙酰化组蛋白多肽的相互作用。AlphaLISA试剂盒可用于抑制实验的高通量筛选平台。



GST-tagged-bromodomain (GST-BRD) 与生物素化的乙酰化多肽 (biotin-H4) 反应后，再加入链霉亲和素包被的供体微珠和谷胱甘肽标记的受体微珠。GST-BRD与 biotin-H4相互作用从而使得供体微珠与受体微珠相互接近，产生615nm的信号。



BRD4 (BD1) 和 H4K5Ac 多肽的终浓度为 25 nM, Alpha试剂盒检测出浓度范围为 3 pM—10 μM 的 JQ1 抑制曲线。

产品名称	Mark (H4 ac)				检测点	货号
	K5	K8	K12	K16		
AlphaLISA					500	AL601C
BRD4_BD1 / H4 K5Ac Binding Kit	○				5,000	AL601F
AlphaLISA					500	AL602C
BRD4_BD2 / H4 K5Ac Binding Kit	○				5,000	AL602F
AlphaLISA					500	AL603C
BRD4_BD1 / H4 K8Ac Binding Kit	○				5,000	AL603F
AlphaLISA					500	AL604C
BRD4_BD2 / H4 K8Ac Binding Kit	○				5,000	AL604F
AlphaLISA					500	AL605C
BRD4_BD1 / H4 K12Ac Binding Kit	○				5,000	AL605F
AlphaLISA					500	AL606C
BRD4_BD2 / H4 K12Ac Binding Kit	○				5,000	AL606F
AlphaLISA					500	AL609C
BRD4_BD1 / H4 K5,8,12,16Ac Binding Kit	○	○	○	○	5,000	AL609F
AlphaLISA					500	AL610C
BRD4_BD2 / H4 K5,8,12,16Ac Binding Kit	○	○	○	○	5,000	AL610F
AlphaLISA					500	AL611C
BRD4_BD1 / H4 K5,8,12Ac Binding Kit	○	○	○		5,000	AL611F
AlphaLISA					500	AL612C
BRD4_BD2 / H4 K5,8,12Ac Binding Kit	○	○	○		5,000	AL612F
AlphaLISA					500	AL613C
BRD4_BD1 / H4 K5,12Ac Binding Kit	○	○			5,000	AL613F
AlphaLISA					500	AL614C
BRD4_BD2 / H4 K5,12Ac Binding Kit	○	○			5,000	AL614F
AlphaLISA					500	AL615C
BRD4_BD1 / H4 K8,16Ac Binding Kit	○		○		5,000	AL615F
AlphaLISA					500	AL616C
BRD4_BD2 / H4 K8,16Ac Binding Kit	○		○		5,000	AL616F

产品名称	检测点	货号
AlphaLISA Total Histone H4 Cell Detection Kit	500 5,000	AL729C AL729F

### 试剂盒组分

- Glutathione AlphaLISA Acceptor beads
- Streptavidin-coated Donor beads
- Biotinylated H4 peptide, lyophilized
- GST-tagged BRD4
- AlphaLISA 5X Epigenetics Buffer 1
- AlphaLISA 30X Epigenetics Buffer Supple- ment



\*\* EpiCypher Inc.是一家专业研究表观遗传学和染色质领域的公司，提供真正高质量的产品。

该试剂盒是与EpiCypher公司合作研发。

# Protein-Protein Interaction

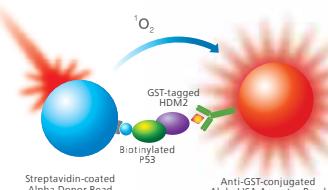
利用供体与受体微珠结合，可建立检测蛋白与蛋白相互作用的目标化合物筛选系统。



通过选择组合Alpha工具盒已有的供体微珠与受体微珠，可检测蛋白一蛋白相互作用，或蛋白一核酸相互作用等多种测定系统。除以下相互作用方式外，其他的参考方法可在TechSupport ([www.perkinelmer.com](http://www.perkinelmer.com)) 中查询相关的文献列表。

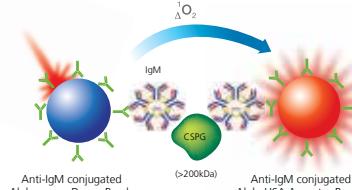
## p53/ HDM2 结合的低亲和性 (Protein/Protein 相互作用)

- AlphaScreen Streptavidin Donor Beads (#6760002)
- AlphaLISA Anti-GST Acceptor Beads (#AL110)



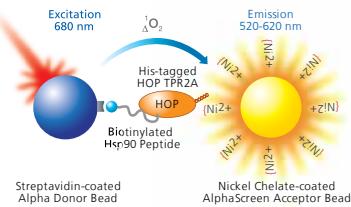
## 通过检测IgE的五聚体来测定 硫酸软骨素蛋白

- AlphaScreen Unconjugated Donor Beads (#6762013)
- AlphaLISA Anti-IgM Acceptor Beads (#AL130)



## 蛋白-蛋白相互作用的抑制剂测定

- AlphaScreen Nickel chelate detection kit (#6760619)

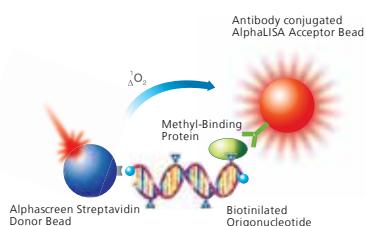


## Reference

Yi, F., Zhu, P., Southall, N., Inglese, J., Austin, C.P., Zheng, W. & Regan, L. An AlphaScreenbased high-throughput screen to identify inhibitors of Hsp90-cochaperone interaction. *J Biomol Screen* 14, 273-281 (2009). Dept. Molecular Biophysics & Biochemistry, Yale University NIH Chemical Genomics Center

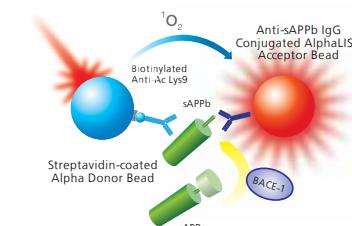
## DNA甲基转移酶测定 (DNA/Protein 相互作用)

- AlphaScreen Streptavidin Donor Beads (#6760002)
- AlphaLISA Unconjugated Acceptor Beads (#6772001)



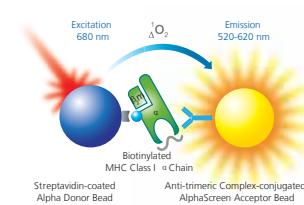
## 蛋白酶 (BACE-1) 测定

- AlphaScreen Streptavidin Donor Beads (#6760002)
- AlphaLISA Unconjugated Acceptor Beads (#6772001)



## MHC I 类, 以及 II 类多肽结合

- AlphaScreen Streptavidin Donor Beads (#6760002)
- AlphaLISA Unconjugated Acceptor Beads (#6772001)

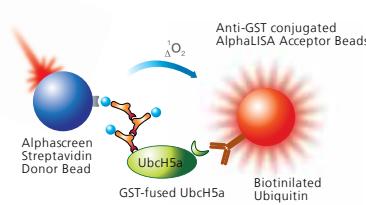


## Reference

Justesen, S., Harndahl, M., Lambeth, K., Nielsen, L.L., Roder, G. & Buus, S. Functional recombinant MHC class II molecules and hightthroughputpeptide-binding assays. *Immunome Res* 5: 2(2009). Harndahl, M., Justesen, S., Lambeth, K., Roder, G., Nielsen, M. & Buus, S. Peptide binding to HLA class I molecules: homogenous, hightthroughput screening and affinity assays. *J Biomol Screen* 14(2), 173-180(2009).

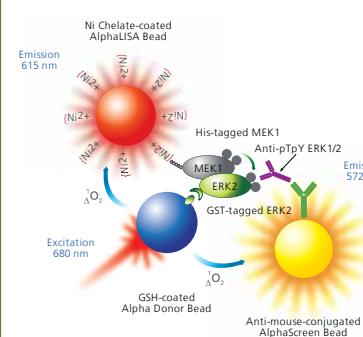
## 泛素测定 (E2的泛素化)

- AlphaScreen Streptavidin Donor Beads (#6760002)
- AlphaLISA Anti-GST Acceptor Beads (#AL110)



## 蛋白磷酸化及Total磷酸化同时检测

- AlphaScreen Mouse IgG detection kit (#6760606)
- Nickel chelate AlphaLISA Acceptor beads (#AL108)



## Reference

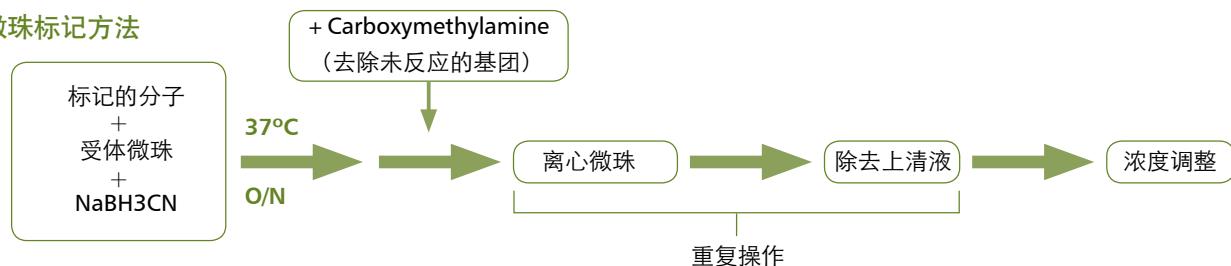
Arcand, M., Roby, P., Bossé, R., Lipari, F., Padrós, J., Beaudet, L., Marcil, A. & Dahan, S. Single-well monitoring of protein-protein interaction and phosphorylation dephosphorylation events. *Biochemistry* 49 (15), 3213-3215 (2010).

# Alpha Tool Box

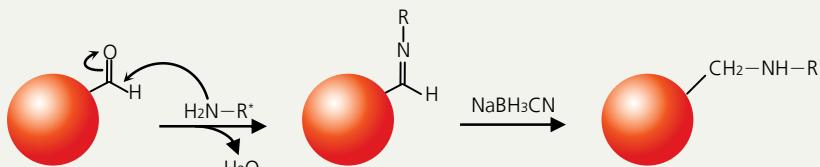
## 未标记的受体与供体微珠 (AlphaScreen/AlphaLISA/AlphaPlex 545)

未标记的微珠表面具有醛基 (-CHO) 活性基团，可以与标记分子上的氨基 (-NH2) 通过还原胺化反应相互结合。可用简单的步骤来进行蛋白微珠标记实验，再将未标记的蛋白分子通过洗涤和离心方法进行分离。

### 微珠标记方法



### 微珠表面结构式



\*R: 标记的分子, 如抗体

### 未标记的供体裸珠

产品名称	货号	检测点
AlphaScreen	1 mg	6762013
Unconjugated Donor beads	5 mg	6762011
	50 mg	6762012

数量: 1 mg 的量在终浓度为 20  $\mu\text{g}/\text{mL}$ 、25  $\mu\text{L}/\text{well}$  的反应体系下可测 2000 个实验点。

保存: + 4°C。

### AlphaScreen标记试剂盒

产品名称	货号	检测点
AlphaScreen Conjugation kit	2 mg each	6760000K

数量: 终浓度为 20  $\mu\text{g}/\text{mL}$ 、25  $\mu\text{L}/\text{well}$  的反应体系下可测 4,000 个实验点。

组分: 链霉亲和素包被的供体微珠, 无标签的 AlphaScreen 受体裸珠。

保存: + 4°C。

### 未标记的受体裸珠

产品名称	货号	检测点
AlphaScreen	1 mg	6762003
Unconjugated Donor beads	5 mg	6762001
	50 mg	6762002
AlphaLISA	1 mg	6772001
Unconjugated Acceptor beads	5 mg	6772002
	50 mg	6772003
AlphaPlex 545	1 mg	6782001
Unconjugated Acceptor beads	5 mg	6782002
	50 mg	6782003
AlphaPlex 645	1 mg	6792001
Unconjugated Acceptor beads	5 mg	6792002
	50 mg	6792003

数量: AlphaScreen 受体微珠 1 mg 的量, 在终浓度为 20  $\mu\text{g}/\text{mL}$ 、25  $\mu\text{L}/\text{well}$  的反应体系下可测 2,000 个实验点。

AlphaScreen 受体微珠 1 mg 的量, 在终浓度为 10  $\mu\text{g}/\text{mL}$ 、50  $\mu\text{L}/\text{well}$  的反应体系下可测 2,000 个实验点。

保存: + 4°C。

## 标记的供体微珠

### 标记的供体微珠

产品名称	货号	检测点
Streptavidin-coated Donor beads	1 mg	6760002S
	5 mg	6760002
	50 mg	6760002B
Glutathione-coated Donor beads	1 mg	6765300
	5 mg	6765301
	25 mg	6765302
Nickel Chelate Donor beads	1 mg	AS101D
	5 mg	AS101M
	25 mg	AS101R
Protein A Donor Beads	1 mg	AS102D
	5 mg	AS102M
	25 mg	AS102R
Anti-FLAG Donor Beads	1 mg	AS103D
	5 mg	AS103M
	25 mg	AS103R

产品名称	货号	检测点
Anti-mouse IgG Donor Beads	1 mg	AS104D
	5 mg	AS104M
	25 mg	AS104R
Anti-Rabbit IgG Donor Beads	1 mg	AS105D
	5 mg	AS105M
	25 mg	AS105R
Strep-Tactin Donor Beads	1 mg	AS106D
	5 mg	AS106M
	25 mg	AS106R
Anti-Digoxigenin Donor Beads	1 mg	AS107D
	5 mg	AS107M
	25 mg	AS107R

数量: 标记的 1 mg 微珠在终浓度为 20  $\mu\text{g}/\text{mL}$ 、25  $\mu\text{L}/\text{well}$  的反应体系中 (AlphaScreen 实验), 可测 2,000 个实验点。在终浓度为 40  $\mu\text{g}/\text{mL}$ 、50  $\mu\text{L}/\text{well}$  的反应体系中 (AlphaLISA 实验) 可测 500 个实验点。

保存: + 4°C。

## 标记的受体微珠 (AlphaScreen®/AlphaLISA/AlphaPlex 545) \*Protein A 标记

在没有直接抗待测物抗体的情况下，免疫球蛋白受体微珠可用于间接检测待测分子（左下表格）。抗融合标签抗体标记的受体微珠可与含有标签的待测蛋白相互作用（右下例子），或者可以检测生物素标记的抗体与标签标记的蛋白相互作用。

### 免疫球蛋白检测受体微珠

产品名称	货号	检测点
AlphaScreen受体微珠		
Protein A	5 mg	6760137M
AlphaScreen Acceptor beads	25 mg	6760137R
AlphaLISA(Eu)受体微珠		
Protein A	250 µg	AL101C
AlphaLISA Acceptor Beads	5 mg	AL101M
	25 mg	AL101R
Protein G	250 µg	AL102C
AlphaLISA Acceptor Beads	5 mg	AL102M
	25 mg	AL102R
Protein L	250 µg	AL126C
AlphaLISA Acceptor Beads	5 mg	AL126M
	25 mg	AL126R
Anti-Human IgG	250 µg	AL103C
AlphaLISA Acceptor Beads	5 mg	AL103M
	25 mg	AL103R
Anti-Human IgG1 (PK/PD)	250 µg	AL141C
AlphaLISA Acceptor Beads	5 mg	AL141M
	25 mg	AL141R
Anti-Human IgG4 (PK/PD)	250 µg	AL142C
AlphaLISA Acceptor Beads	5 mg	AL142M
	25 mg	AL142R
Anti-Human IgG1 (isotyping)	250 µg	AL153C
AlphaLISA Acceptor Beads	5 mg	AL153M
	25 mg	AL153R
Anti-Human IgG2 (isotyping)	250 µg	AL154C
AlphaLISA Acceptor Beads	5 mg	AL154M
	25 mg	AL154R
Anti-Human IgG3 (isotyping)	250 µg	AL155C
AlphaLISA Acceptor Beads	5 mg	AL155M
	25 mg	AL155R
Anti-Human IgG4 (isotyping)	250 µg	AL156C
AlphaLISA Acceptor Beads	5 mg	AL156M
	25 mg	AL156R
Anti-Mouse IgG	250 µg	AL105C
AlphaLISA Acceptor Beads	5 mg	AL105M
	25 mg	AL105R
Anti-Mouse IgM	250 µg	AL130C
AlphaLISA Acceptor Beads	5 mg	AL130M
	25 mg	AL130R
Anti-Mouse IgG1 (isotyping)	250 µg	AL157C
AlphaLISA Acceptor Beads	5 mg	AL157M
	25 mg	AL157R
Anti-Mouse IgG2a (isotyping)	250 µg	AL158C
AlphaLISA Acceptor Beads	5 mg	AL158M
	25 mg	AL158R
Anti-Mouse IgG2a (isotyping)	250 µg	AL159C
AlphaLISA Acceptor Beads	5 mg	AL159M
	25 mg	AL159R
Anti-Mouse IgG3 (isotyping)	250 µg	AL160C
AlphaLISA Acceptor Beads	5 mg	AL160M
	25 mg	AL160R
Anti-Mouse IgE (isotyping)	250 µg	AL161C
AlphaLISA Acceptor Beads	5 mg	AL161M
	25 mg	AL161R
Anti-Mouse IgM (isotyping)	250 µg	AL162C
AlphaLISA Acceptor Beads	5 mg	AL162M
	25 mg	AL162R
Anti-Mouse IgA	250 µg	AL162C
AlphaLISA Acceptor Beads	5 mg	AL163M
	25 mg	AL163R
Anti-Mouse total IgG	250 µg	AL164C
AlphaLISA Acceptor Beads	5 mg	AL164M
	25 mg	AL164R

### 免疫球蛋白检测受体微珠 (续)

产品名称	货号	检测点
Anti-Mouse total IgG	250 µg	AL164C
AlphaLISA Acceptor Beads	5 mg	AL164M
	25 mg	AL164R
Anti-Rat IgG	250 µg	AL106C
AlphaLISA Acceptor Beads	5 mg	AL106M
	25 mg	AL106R
Anti-Rabbit IgG	250 µg	AL104C
AlphaLISA Acceptor Beads	5 mg	AL104M
	25 mg	AL104R
Anti-Goat IgG	250 µg	AL107C
AlphaLISA Acceptor Beads	5 mg	AL107M
	25 mg	AL107R
Anti-Chicken IgY	250 ug	AL313C
AlphaLISA Acceptor Beads	5 mg	AL313M
	25 mg	AL313R
Anti-Sheep IgG	250 µg	AL132C
AlphaLISA Acceptor Beads	5 mg	AL132M
	25 mg	AL132R
Anti-Bovine total IgG	250 µg	AL165C
AlphaLISA Acceptor Beads	5 mg	AL165M
	25 mg	AL165R
Anti-Bovine IgG 1	250 µg	AL166C
AlphaLISA Acceptor Beads	5 mg	AL166M
	25 mg	AL166R
Anti-Bovine IgG 2	250 µg	AL167C
AlphaLISA Acceptor Beads	5 mg	AL167M
	25 mg	AL167R
Anti-Bovine IgM	250 µg	AL168C
AlphaLISA Acceptor Beads	5 mg	AL168M
	25 mg	AL168R
Anti-Bovine IgA	250 µg	AL169C
AlphaLISA Acceptor Beads	5 mg	AL169M
	25 mg	AL169R
AlphaPlex 545(Tb)受体微珠		
Protein A	250 µg	AP101TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP101TB-M
	25 mg	AP101TB-R
Anti-Human IgG	250 µg	AP103TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP103TB-M
	25 mg	AP103TB-R
Anti-Rabbit IgG	250 µg	AP104TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP104TB-M
	25 mg	AP104TB-R
Anti-Mouse IgG	250 µg	AP105TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP105TB-M
	25 mg	AP105TB-R
AlphaPlex 645(Sm)受体微珠		
Protein A	250 µg	AP101SM-C
AlphaPlex 645 Acceptor Beads	5 mg	AP101SM-M
	25 mg	AP101SM-R
Anti-Human IgG	250 µg	AP103SM-C
AlphaPlex 645 Acceptor Beads	5 mg	AP103SM-M
	25 mg	AP103SM-R
Anti-Mouse IgG	250 µg	AP105SM-C
AlphaPlex 645 Acceptor Beads	5 mg	AP105SM-M
	25 mg	AP105SM-R

数量：AlphaScreen蛋白A受体微珠5 mg，在终浓度为20µg/mL、25 µL/well的反应体系中，可测10,000个点；

AlphaLISA 受体微珠 250 µg 、在终浓度为10µg/mL、50µL/well的反应体系中、可测500个点。

保存：+ 4°C。

# Alpha Tool Box

## 融合标签检测受体微珠

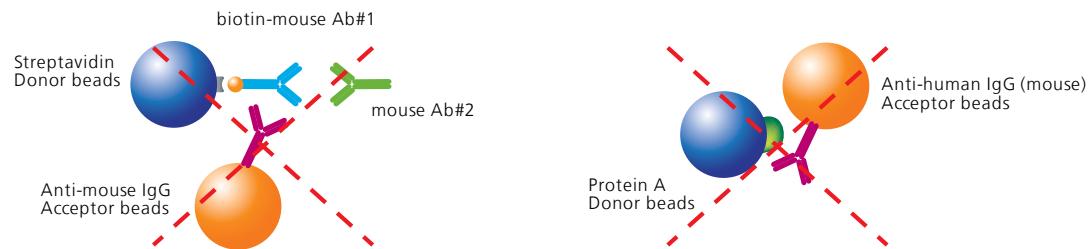
产品名称	货号	检测点
<b>AlphaLISA (Eu)受体微珠</b>		
Nickel Chelate	250 µg	AL108C
AlphaLISA Acceptor Beads	5 mg	AL108M
	25 mg	AL108R
Glutathione	250 µg	AL109C
AlphaLISA Acceptor Beads	5 mg	AL109M
	25 mg	AL109R
Anti-GST	250 µg	AL110C
AlphaLISA Acceptor Beads	5 mg	AL110M
	25 mg	AL110R
Anti-c-Myc	250 µg	AL111C
AlphaLISA Acceptor Beads	5 mg	AL111M
	25 mg	AL111R
Anti-FLAG	250 µg	AL112C
AlphaLISA Acceptor Beads	5 mg	AL112M
	25 mg	AL112C
Anti-DIG	250 µg	AL113C
AlphaLISA Acceptor Beads	5 mg	AL113M
	25 mg	AL113R
Streptavidin	250 µg	AL125C
AlphaLISA Acceptor Beads	5 mg	AL125M
	25 mg	AL125R
Anti-FITC	250 µg	AL127C
AlphaLISA Acceptor Beads	5 mg	AL127M
	25 mg	AL127R
Anti-His Tag	250 µg	AL128C
AlphaLISA Acceptor Beads	5 mg	AL128M
	25 mg	AL128R
Anti-V5	250 µg	AL129C
AlphaLISA Acceptor Beads	5 mg	AL129M
	25 mg	AL129R
Anti-MBP	250 µg	AL134C
AlphaLISA Acceptor Beads	5 mg	AL134M
	25 mg	AL134R
Strep-Tactin	250 µg	AL136C
AlphaLISA Acceptor Beads	5 mg	AL136M
	25 mg	AL136R
Anti-HA	250 µg	AL170C
AlphaLISA Acceptor Beads	5 mg	AL170M
新品	25 mg	AL170R
<b>AlphaPlex 545(Tb)受体微珠</b>		
Glutathione	250 µg	AP109TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP109TB-M
	25 mg	AP109TB-R
Streptavidin	250 µg	AP125TB-C
AlphaPlex 545 Acceptor beads	5 mg	AP125TB-M
	25 mg	AP125TB-R
Anti-FLAG	250 µg	AP112TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP112TB-M
	25 mg	AP112TB-R
Anti-6x His	250 µg	AP128TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP128TB-M
	25 mg	AP128TB-R
Anti-HA	250 µg	AP170TB-C
AlphaPlex 545 Acceptor Beads	5 mg	AP170TB-M
	25 mg	AP170TB-R
<b>AlphaPlex 645(Sm) 受体微珠</b>		
Streptavidin	250 µg	AP125SM-C
AlphaPlex 645 Acceptor Beads	5 mg	AP125SM-M
	25 mg	AP125SM-R
Anti-FLAG	250 µg	AP112SM-C
AlphaPlex 645 Acceptor Beads	5 mg	AP112SM-M
	25 mg	AP112SM-R
Anti-6x His	250 µg	AP128SM-C
AlphaPlex 645 Acceptor Beads	5 mg	AP128SM-M
	25 mg	AP128SM-R

## 糖链受体检测受体微珠

产品名称	货号	检测点
LCA (Lens culinaris agglutinin)	250 µg	AL104C
AlphaLISA Acceptor Beads <i>(Detection of fucosylated glycoproteins)</i>	5 mg	AL104M
	25 mg	AL104R

## Tips: 供体微珠与受体微珠的结合方式

受体微珠与供体微珠不一定以我们需要的方式结合。如在实验中若同时使用蛋白A偶联的供体微珠和hIgG偶联的受体微珠，会因为供体微珠与受体微珠直接相互作用而使得背景值非常高。



AlphaLISA Acceptor Bead	Alpha Donor Bead							
	Protein A	Anti-FLAG	Anti-Mouse	Anti-Rabbit	Strep-Tactin <sup>®</sup>	Ni chelate	GSH	Streptavidin
Anti-Chicken IgY	✗	✓	✓	✗	✓	✓	✓	✓
Anti-cMyc	✓	✓	✓	✓	✓	✓	✓	✓
Anti-Digoxigenin	✓	✓	✗	✓	✓	✓	✓	✓
Anti-FITC	✗	✓	✗	✓	✓	✓	✓	✓
Anti-FLAG	✓	✓	✗	✓	✓	✓	✓	✓
Anti-GFP	✓	✓	✗	✓	✓	✓	✓	✓
Anti-Goat IgG	✓	✓	✓	✗	✓	✓	✓	✓
Anti-GST	✓	✓	✓	✓	✓	✓	✓	✓
Anti-His	✓	✓	✗	✓	✓	✓	✓	✓
Anti-Human IgG	✗	✓	✗	✓	✓	✓	✓	✓
Anti-MBP	✓	✓	✓	✓	✓	✓	✓	✓
Anti-Mouse IgG	✓	✗	✓	✓	✓	✓	✓	✓
Anti-Mouse IgM	✓	✓	✗	✓	✓	✓	✓	✓
Anti-Rabbit IgG	✓	✓	✓	✓	✓	✓	✓	✓
Anti-Rat IgG	✓	✗	✓	✓	✓	✓	✓	✓
Anti-Sheep IgG	✗	✓	✗	✗	✓	✓	✓	✓
Anti-V5	✗	✓	✗	✓	✓	✓	✓	✓
Glutathione	✓	✓	✓	✓	✓	✓	✓	✓
Ni chelate	✓	✓	✓	✓	✓	✓	✓	✓
Protein A	✓	✓	✓	✓	✓	✓	✓	✓
Protein G	✓	✓	✗	✗	✓	✓	✓	✓
Protein L	✓	✓	✓	✓	✓	✗	✓	✓
Strep-Tactin <sup>®</sup>	✓	✓	✓	✓	✓	✓	✓	✓
Streptavidin	✓	✓	✓	✓	✓	✓	✓	✓

✓	You can use this pair
✗	This pair might give high background
✗	Not recommended

# Alpha Tool Box

## AlphaScreen IgG / 融合标签检测试剂盒

链霉亲和素包被的供体微珠和抗IgG抗体标记的受体微珠的AlphaScreen试剂盒。

### 融合标签检测试剂盒

产品名称	货号	检测点
GST detection kit	500	6760603C
	10,000	6760603M
	50,000	6760603R
DIG detection kit	500	6760604C
	10,000	6760604M
	50,000	6760604R
FITC detection kit	500	6760605C
	10,000	6760605M
	50,000	6760605R
c-myc detection kit	500	6760611C
	10,000	6760611M
	50,000	6760611R
HA (hemagglutinin) detection kit	500	6760612C
	10,000	6760612M
	50,000	6760612R
FLAG (M2) detection kit	500	6760613C
	10,000	6760613M
	50,000	6760613R
Ni chelate detection kit	500	6760619C
	10,000	6760619M
	50,000	6760619R
5μM Biotinylated-His	2.5 μmole	6760303M
5μM Biotinylated-GST	0.75 μmole	6760305M

### IgG检测试剂盒

产品名称	货号	检测点
Mouse IgG detection kit	500	6760606C
	10,000	6760606M
	50,000	6760606R
Rabbit IgG detection kit	500	6760607C
	10,000	6760607M
	50,000	6760607R
Protein A detection kit	500	6760617C
	10,000	6760617M
	50,000	6760617R

组份：10X buffer, Biotinylated-IgG or tag, Streptavidin Donor beads, Acceptor beads.

数量：在终浓度为20μg/mL、25μL/well体系的384孔中进行实验。

保存：+ 4°C。

## 其他微珠（干扰测定、加样精确度、仪器设备验证相关微珠）

### TruHits 试剂盒 ..... 含链霉亲和素包被供体微珠和生物素化受体

此试剂盒是为了评估受体微珠颜色淬灭，单线态氧在实验中的淬灭，某些化合物如生物素类似物分子对Alpha实验信号的干扰等因素。

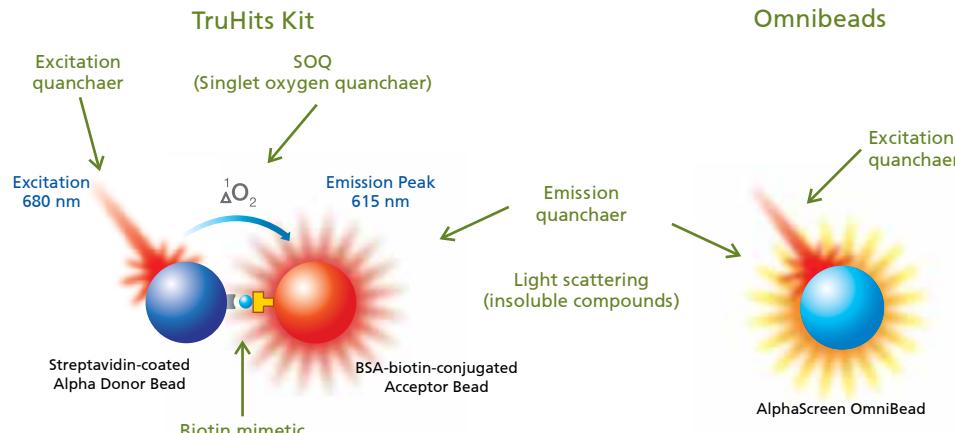
### Omnibeads ..... 该微珠上包被了产生Alpha信号所需的所有物质，激发光激发微珠后可直接产生信号。可验证仪器对每个孔的读取效率和精度，可通过Omnibeads校准仪器探头。

产品名称	货号	检测点
AlphaScreen TruHits Kit	1,000	6760627D
	10,000	6760627M
	1,000	AL900D
TruHits Kit	10,000	AL900M
	1,000	AP900D
	10,000	AP900M
Omnibeads	1,000	6760626D
	10,000	6760626M
	50,000	6760626R

数量：在试剂盒指定用量的情况下，使用384孔板的量。

保存：+ 4°C。

## 信号淬灭的种类



## 反应缓冲液

链霉亲和素包被的供体微珠和抗IgG抗体标记的受体微珠的AlphaScreen试剂盒。

Immunoassay Buffer .....	250mM HEPES, 1% Casein, 5%Triton X-100, 10mg/mL Dextran-500, 0.5% Proclin-300
Universal Buffer (5X) .....	5 × PBS, 0.5% BSA, 0.05% Proclin-300
HiBlock Buffer (10X) .....	250mM HEPES, 1% Casein, 5% Triton X-100, 10mg/mL Dextran-500, 0.5% Proclin-300, 5% gelatin, 5% BSA
Dissociation Buffer (3X) .....	AlphaLISA Residual Protein A kit (AL287)
NaCl Buffer (5X) .....	AlphaLISA CCL2/MCP1 kit (AL244) AlphaLISA Mouse Albumin kit (AL508)

产品名称	货号	检测点
AlphaLISA	10 mL	AL000C
Immunoassay Buffer (10X)	100 mL	AL000F
AlphaLISA	10 mL	AL001C
Universal Buffer (5X)	100 mL	AL001F
AlphaLISA	10 mL	AL003C
Lysis Buffer (5X)	100 mL	AL003F
AlphaLISA	10 mL	AL004C
HiBlock Buffer (10X)	100 mL	AL004F
AlphaLISA	10 mL	AL006C
Dissociation Buffer (3X)	100 mL	AL006F
AlphaLISA	10 mL	AL007C
NaCl Buffer (5X)	100 mL	AL007F

## 微孔板

AlphaPlate .....	消除白色微孔板孔与孔之间的串扰专门用于Alpha 实验检测的灰白色微孔板。
1/2 AreaPlate .....	孔的表面积设计为正常96孔板的一半。
CulturePlate .....	适用于基于细胞学水平的实验。具有很好的内部表面处理和γ射线灭菌。适用于贴壁细胞的培养。
TopSeal-A .....	防止样品的蒸发。建议密封在微孔板上。不会影响到信号值，可放置一段时间来进行检测。

产品名称	材质	规格	货号
AlphaPlate-384	聚苯乙烯	50 块	6005350
		200 块	6005359
AlphaPlate-384, shallow well	聚苯乙烯	50 块	6008350
AlphaPlate-1536	聚苯乙烯	200 块	6008359
1/2 AreaPlate-96 (白)	聚苯乙烯	50块	6004350
		200块	6005560
		50块	6005569
CulturePlate-96 (白)	聚苯乙烯	160块	6005680
		200块	6005689
		50块	6007680
CulturePlate-384 (白)	聚苯乙烯	160块	6007688
		200块	6007689
CulturePlate-1536 (白)	聚苯乙烯	50块	6004680
TopSeal-A 96 / 384	聚苯乙烯	100片	6050185

## 定制服务

### 受体配体微珠的标记服务

- 可将靶蛋白，配体或抗体标记到微珠上。
- 至少需标记1 mg。
- 可直接提供标记分子。

## 自定义检测服务

### 成功案例

- Anti-g8 Antibodies
- pACC
- Phospho-PERK
- CNTO 1959
- PAG (gestetion protein bovine)
- Phospho-PLC1 (Y783)
- Dimethyl H3 Peptide
- PCSK9
- Phospho-Tyr ERK
- DOT1 组蛋白甲基化
- Phospho-eEF2
- E1/E2 + CD81 相互作用
- FoxP3
- Phospho-elf2 α
- Serotonin
- Insulin Receptor 自磷酸化
- Phospho-LRP6
- SIGMA-1 Receptor/BIP Chaperone
- KIR
- Phospho-Merlin
- LRRK2
- Phospho-Myosine Phosphatase

# Alpha 检测设备

Alpha 检测需要使用配备了激光（发射 680 nm 激发光）的专用设备。本公司提供 3 种能够以高灵敏度和高精度检测 Alpha 信号的 Alpha 检测设备。

\* 有关各设备的详细信息请咨询本公司。

## EnVision

EnVision 是所有检测技术中灵敏度最高并具有极快的检测速度和高通量的一款独一无二的微孔板检测仪，其优异的灵活性和扩展性确保能适用于各种检测应用。用户可选择所需的检测功能进行配置。本系统还可用于 AlphaPlex 检测。

标准配置	选件
<ul style="list-style-type: none"><li>• 吸收光（滤光器）</li><li>• 荧光强度（滤光器、顶部/底部检测）</li><li>• 发光</li><li>• 荧光偏振</li><li>• 时间分辨荧光</li></ul>	<ul style="list-style-type: none"><li>• <b>Alpha 检测</b>（标准或 HTS）</li><li>• <b>AlphaPlex 检测</b>（标准）</li><li>• 高灵敏度发光</li><li>• 吸收光（单色器）</li><li>• 荧光强度（单色器）</li><li>• 377 nm 激光</li><li>• 温度控制</li><li>• 分液器</li><li>• 载板架（20 张/50 张）</li><li>• 条形码读取器</li><li>• 符合 Part11 的软件</li></ul>



## EnSight

EnSight 是可配置吸收光和荧光单色器、高灵敏度发光以及 Alpha 检测等功能的灵活多模式微孔板检测仪。本系统还可用于微孔成像检测。

标准配置	选件
<ul style="list-style-type: none"><li>• 吸收光（滤光器）</li><li>• 条形码读取器</li><li>• 温度控制</li></ul>	<ul style="list-style-type: none"><li>• <b>Alpha 检测</b></li><li>• 吸收光（单色器）</li><li>• 荧光强度（单色器）</li><li>• 时间分辨荧光</li><li>• 高灵敏度发光</li><li>• 无标记</li><li>• 微孔成像</li></ul>



## EnSpire

EnSpire 是一款性能卓越、界面友好的多模式微孔板检测仪。我们提供 2 种型号，一种是 Alpha 检测专用型号，一种是除 Alpha 以外还可配备吸光、荧光、时间分辨荧光、发光、无标记技术的型号。

型号	标准配置	选件
EnSpire Alpha	<ul style="list-style-type: none"><li>• <b>Alpha 检测</b></li><li>• 条形码读取器</li></ul>	<ul style="list-style-type: none"><li>• 载板架（20 张/50 张）</li><li>• 符合 Part11 的软件</li></ul>
EnSpire	<ul style="list-style-type: none"><li>• 吸收光（滤光器）</li><li>• 条形码读取器</li></ul>	<ul style="list-style-type: none"><li>• <b>Alpha 检测</b></li><li>• 高灵敏度发光</li><li>• 吸收光（单色器）</li><li>• 荧光强度（单色器）</li><li>• 荧光强度底部检测</li><li>• 时间分辨荧光</li><li>• 无标记</li><li>• 载板架（20 张/50 张）</li><li>• 分液器</li><li>• 温度控制</li><li>• 符合 Part11 的软件</li></ul>



## 生命科学技术中心

第一步扫一扫二维码或点击以下链接

<http://www.perkinelmer.com/tools/resourcelibrary#/resources>



第二步选择并勾选所需

Home > Resources

### Resources

Use this section to access product documents, tools, and forms that are helpful in purchase decisions, ordering, and getting the most from your PerkinElmer products.

Search Product Documents Tools and Forms

#### Search

Search

- Application Brief
- Booklet
- Case Study
- Data Sheet
- Flyer
- Manual
- Poster
- Product Information Bulletin
- Specification Sheet
- White Paper

Application Note

Brochure

Catalog

FAQ

Guide

Material Safety Data Sheets

Poster (Scientific)

Product Note

Technical Note

Article

Calendar

Certificate of Analysis

Field Application Report

Handbook

Newsletter

Product Info

Resources

Training



点击!

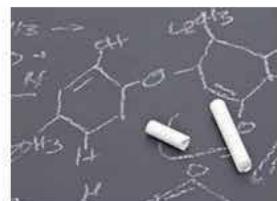
Submit

第三步：点击“Submit”获取相关信息

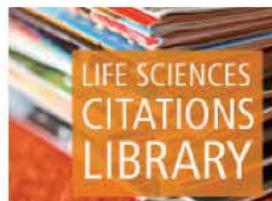
CoA (Certificates of analysis) 下载



技术资料下载



文献库检索



聆听 Webinar





## 珀金埃尔默企业管理（上海）有限公司

### 上海总公司

地址：上海浦东新区张江高科  
技园区张衡路1670号  
电话：+86 21-6064 5888  
传真：+86 21-6064 5999  
邮编：201203

### 北京分公司

地址：北京朝阳区酒仙桥路14号院  
兆维工业园甲2号楼1层东侧单元  
电话：+86 10 8434 8999  
传真：+86 10 8434 8988  
邮编：100015

### 成都分公司

地址：成都市高新区西芯大道5号  
汇都总部园6栋3楼  
电话：+86 28-8785 7220  
传真：+86 28-8785 7221  
邮编：610016

### 武汉分公司

地址：武汉市武昌区临江大道96号  
武汉万达中心写字楼1808-1809室  
电话：+86 27-8891 3055  
传真：+86 27-8891 3380  
邮编：430062

### 西安分公司

地址：陕西省西安市高新区锦业路69号  
创业研发园A座1009室  
电话：+86 029-8129 2671  
传真：+86 029-8129 2126  
邮编：710077

### 广州分公司

地址：广州市荔湾区芳村大道  
下市直街1号信义会馆12号  
电话：+86 20-3789 1888  
传真：+86 20-3789 1899  
邮编：510370

中文网址：[www.perkinelmer.com.cn](http://www.perkinelmer.com.cn)  
客户服务电话：[800 820 5046](tel:8008205046)

要获取全球办事处的完整列表，请访问 [www.perkinelmer.com>ContactUs](http://www.perkinelmer.com>ContactUs)  
版权所有 ©2016 PerkinElmer, Inc. 保留所有权利。PerkinElmer® 是 PerkinElmer, Inc. 的注册商标。所有其它商标均为其各自所有者的财产。  
所有数据仅供参考，PerkinElmer 拥有最终解释权。

