

Alpha SureFire® Ultra™ HV Multiplex

Terbium SureFire® Ultra™ HV Assay Kit For multiplexing with AlphaLISA® SureFire® Ultra™ (ALSU) HV assay kits

Assay Kits Manual

Assay Points	Catalog #
100 (96 well format)	ALSU-XXXX-X-HV

This Manual is a generic manual for all the kits. For target-specific information, relating to Kit Specificity, Control Lysates and Representative Data, please refer to the Technical Data Sheet of the kit, also available from www.perkinelmer.com

For Research Use Only. Not for use in Diagnostic Procedures.

Note: See important kit disposal information on page 5 of this manual

For a full, electronic, version of this manual, please go to:
<http://www.perkinelmer.com/category/alpha-surefire-kits>

Alpha SureFire® Ultra™ HV Multiplex

Assay Principle

The Alpha SureFire Ultra Multiplex kits allow the rapid, sensitive, and quantitative detection of two phosphoprotein targets in each well of an assay plate. This Multiplex measurement is achieved by the use of two types of Alpha Acceptor beads that emit at distinct wavelengths (Terbium, 545nm and Europium, 615nm).

The two distinct Alpha Acceptor beads report their binding to distinct antigens through their association with specific assay antibodies, as indicated below.

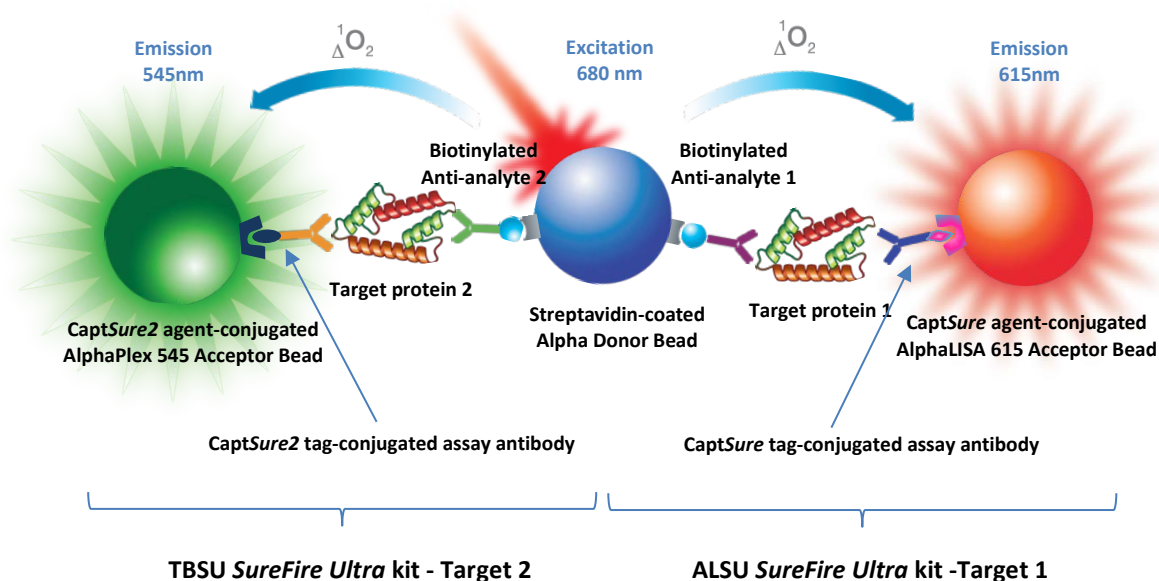
Terbium SureFire Ultra HV Assay kit for multiplexing with another AlphaLISA SureFire Ultra HV assay kit:

The TBSU assay kit contains reagents to be combined with those of an AlphaLISA SureFire Ultra HV assay kit (Cat No. ALSU-XXXX-X-HV) to allow multiplexed measurement of two phosphoprotein targets. Antibodies in each TBSU or ALSU HV kit measure a single target. Combining two such kits allows the highly flexible selection of two phosphoprotein targets to be measured.

The terbium “Alpha 545 Acceptor Beads” in the TBSU kit are conjugated with a “CaptSure2” agent, which binds a specific CaptSure2 tag on one of the assay antibodies. In combination with the biotinylated antibody provided, the 545 nm signal generated will report levels of the phosphoprotein of interest (Target 2). The TBSU kit is to be used in conjunction with the reagents in an ALSU kit, such that the ALSU kit will report that target of interest by Eu emission at 615nm (Target 1). This is shown diagrammatically below.

The protocol details how to combine these kits to allow multiplexed measurement of the desired targets.

Alpha SureFire® Ultra™ Multiplex Technology



General Information on the Terbium *SureFire® Ultra™ HV* Assay kit

The Terbium *SureFire Ultra HV* Assay (TBSU) kit is used to measure the phosphorylation of a target of interest plus another target of choice from the AlphaLISA *SureFire Ultra HV* (ALSU) list of kits. The TBSU kit must be used in conjunction with an ALSU HV kit, and can not be used alone to measure a single target.

Multiplexing is achieved through the use of two Alpha Acceptor beads. The first (545nm Tb) in this TBSU HV kit will report on the levels of the designated target of interest. The second (615nm Eu) Acceptor bead from an ALSU HV kit reports on the protein designed to be measured using that kit.

The Donor beads in this kit are identical to those in a standard AlphaLISA *SureFire Ultra HV* kit.

We recommend testing separately the signal generated in the two channels (545 and 615 nm) using the control lysates from each kit.

This kit has been formulated to provide optimal signal:background (i.e. S:B) assay windows, and to perform without interference in the presence of extraneous antibodies.

Kit-Specificity Information / Control Lysate Information / Representative Data

See Technical Data Sheet in assay kit box.

Unless otherwise indicated, the antibodies in each TBSU kit are identical to the antibodies to the same target in the corresponding ALSU kit, except for the CaptSure or CaptSure2 tag.

The assay utilizes the bead-based Alpha Technology, and requires an Alpha Technology-compatible plate reader capable of reading dual emission wavelengths. See www.perkinelmer.com/AlphaPlex for more information about the AlphaPlex technology and download the “AlphaPlex Quick Start Guide” and the “AlphaPlex Assay Development Guide” to find guidance about filters and mirrors selection, instrument protocol and channels crosstalk correction.

Note: The buffers in the Alpha *SureFire Ultra* Multiplex kits are the same formulation as the buffers in the AlphaLISA *SureFire Ultra* kits, but they have a different formulation compared to the buffers from the AlphaScreen *SureFire* kits, and buffers from the latter type of kit should not be interchanged.

Kit Contents

Kit Size	100 points
Reaction Buffer 3 - TBSU (Biotinylated anti-target antibody)	1 x 0.09 mL
Reaction Buffer 4 - TBSU (CaptSure2-tagged anti-target antibody)	1 x 0.09 mL
Alpha 545 CaptSure2 Acceptor Beads (2mg/mL in PBS plus 0.05% Proclin-300)	1 x 0.04 mL
Alpha Streptavidin Donor Beads (2mg/mL in PBS plus 0.05% Proclin-300)	1 x 0.04 mL
Positive Control Lysate	1 lyophilized tube to be re-dissolved with 250 µL H ₂ O

The above volumes supplied are in excess to the actual volume required to perform assay.

IMPORTANT: The components of this kit will be used in conjunction with the components of another standard AlphaLISA SureFire Ultra (ALSU) HV kit to provide multiplexing. On page 6 we detail how to set up the assay solutions from both kits.

Storage Conditions Upon Receipt

The kit should be placed at 4°C upon receipt. DO NOT freeze the kit buffers or beads. The Reaction buffers contain antibodies and freeze/thaw cycles can lead to a loss of activity. Alpha Donor Beads need to be stored at 4°C in the dark, and should be returned to the kit box after each use.

The Activation Buffer (use the one from the AlphaLISA SureFire Ultra kit) precipitates at 4°C. To re-dissolve, warm to 37°C and mix before each use. Alternatively, Activation buffer can be stored at room temperature with no loss in activity. All other components to be returned to 4°C after each use.

The Positive control lysate tube should be placed at -20°C or -80°C for long term storage.

See kit box label for expiry date.

Precautions

Only the Alpha Donor beads are light-sensitive. All the other assay reagents can be used under normal light conditions. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (filter #90 from LEE Filters (preferred) or Roscolux filters #389 from Rosco, or the equivalent) can be applied to light fixtures.

Materials Required But Not Provided

Item	Suggested source	Catalog #	Size
1/2AreaPlate™ - 96 assay plate	PerkinElmer Inc.	6005560	50/box
TopSeal-A 384, clear adhesive sealing film	PerkinElmer Inc.	6050185	100/box
Envision®, Enspire® or EnSight™ Alpha-reader	PerkinElmer Inc.	-	-

Table : AlphaPlex Optics for EnVision Multilabel Reader – for complete information about how to set an AlphaPlex reading, please refer to the AlphaPlex Guides available at www.perkinelmer.com/AlphaPlex

	Description	Catalog #	Barcode	Recommendations
Mirrors	AlphaScreen	2101-4010	444	For Tb and Eu single and sequential reading ; not for Sm
	AlphaPlex Single Tb-Eu-Sm	2102-5910	605	Preferred mirror for all sequential AlphaPlex applications
	AlphaPlex Dual Tb-Eu	2102-5900	653	For simultaneous duplexing of Tb with Eu
Filters	AlphaScreen	2100-5710	244	Suitable for AlphaPlex single plexing, not for multiplexing
	Resorufine/ Amplex Red	2100-5570	124	Suitable for Tb single plexing and Tb/Eu duplexing.
	Europium	2100-5090	203	Preferred filter for all Eu applications and multiplexing
	AlphaPlex Tb	2100-5930	701	Preferred filter for all Tb applications and multiplexing

Important disposal information

The Lysis Buffer (5X) – *Ultra* in this kit contains Triton X-100, otherwise known as p-tert-octylphenol ethoxylate.

p-tert-octylphenol ethoxylate must be disposed of as Controlled Waste in accordance with Local Regulations.

Buffer Preparation and Subsequent Storage Conditions

1. Prepare sufficient Acceptor Mix from the ALSU AlphaLISA *SureFire Ultra HV* kit as you would for a standard single target assay, according to that kit manual.
2. The Terbium *SureFire Ultra HV* Assay kit reagents will be added to the pre-prepared ALSU Acceptor mix from step 1 to complete the Multiplex Acceptor Mix.
3. We recommend that you combine the Donor beads from both ALSU and TBSU kits into a single tube prior to use. Donor Mix will then be made up as indicated below.

Components	Preparation
1X Lysis Buffer Note: Lysis Buffer from ALSU kit	Dilute 5X Lysis buffer in MilliQ water to a final concentration of 1X For example: for 10 mL of 1X Lysis Buffer, add: 2 mL of 5X Lysis Buffer – <i>Ultra</i> to 8 mL MilliQ water. Discard unused 1X buffer.
<u>Multiplex Acceptor Mix</u> Prepare Acceptor Mix as normal from ALSU <i>SureFire Ultra HV</i> kit. <u>To this:</u> Add TBSU Reaction Buffer 3 (1:20) Add TBSU Reaction Buffer 4 (1:20) Vortex Add CaptSure2 Alpha 545 Acceptor Beads (1:50) Vortex See flowchart for table	Important: combine the reagents in the order indicated below <u>For example: for 20 samples (require minimum 300 µL of Multiplex Acceptor Mix):</u> Make up 300 µL of ALSU kit Acceptor Mix, according to that kit's protocol. To this ALSU Acceptor Mix add 15 µL of Reaction Buffer 3 and 15 µL of Reaction Buffer 4. Mix combined reagents by vortex. Then add 6µL CaptSure2 Alpha 545 Acceptor Beads. Mix again by vortex. The Multiplex Acceptor mix should be made up and used within 30min for best results. Excess mix should be discarded.
<u>Donor Mix*</u> Alpha Donor beads (1:25) (4 parts or 4%) Note: Dilution buffer from ALSU kit See flowchart for table * Prepare and use under low-light conditions.	Dilute Donor beads 25-fold in Dilution buffer <u>For example: for 20 samples (require minimum 300 µL of Donor Mix):</u> Add 12 µL Donor Beads to 288 µL of Dilution Buffer. Note: Dilution Buffer from ALSU kit. The Donor mix should be made up and used within 30min for best results. Excess mix should be discarded.
Positive control lysate	Reconstitute with 250µL water. Store at -20°C in single use aliquots and use within 3 months. Dilute as required.

Dual target measurement using combined Terbium SureFire® Ultra™ HV Assay kit + AlphaLISA® SureFire® Ultra™ HV target assay kit

A. 2-Plate Assay - assay protocol for adherent cells

Cell Seeding

1. Seed cells (200 µL of cells per well) in 96 well tissue culture plates. Incubate at 37°C overnight in serum-containing media.

Cell Treatment

2. Remove culture media, and stimulate the cells with 50 µL agonists prepared in serum-free media. *(If testing antagonists, prior to stimulation remove culture medium and replace with 50 µL serum-free media containing antagonists. Return cells to 37°C incubator for desired time. 1 hour is often sufficient for signal transduction inhibitors, and 5-20 minutes for receptor antagonists).*

Note: Peptidic agonists and antagonists can often stick to plastic surfaces. To minimize this effect, dilute in serum-free media containing a suitable carrier protein (e.g. 0.1% protease-free BSA).

Lysate Preparation

3. To lyse cells, remove medium from wells, and add freshly prepared 1X Lysis Buffer (50-100 µL per well). Agitate on a plate shaker (~350 rpm) for 10 minutes at room temperature.
4. Take 30 µL of the lysate and transfer to a 96-well 1/2AreaPlate™ for assay. *Add 30 µL of Control lysates to separate wells. We recommend testing a serial dilution of Control lysate (eg 100, 50, 25, 12.5, 6.25 and 0% diluted in 1X Lysis Buffer).*

SureFire Ultra Assay

5. Add 15 µL of Acceptor Mix to wells. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature.
6. Add 15 µL of Donor Mix to wells under subdued light. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature in the dark.

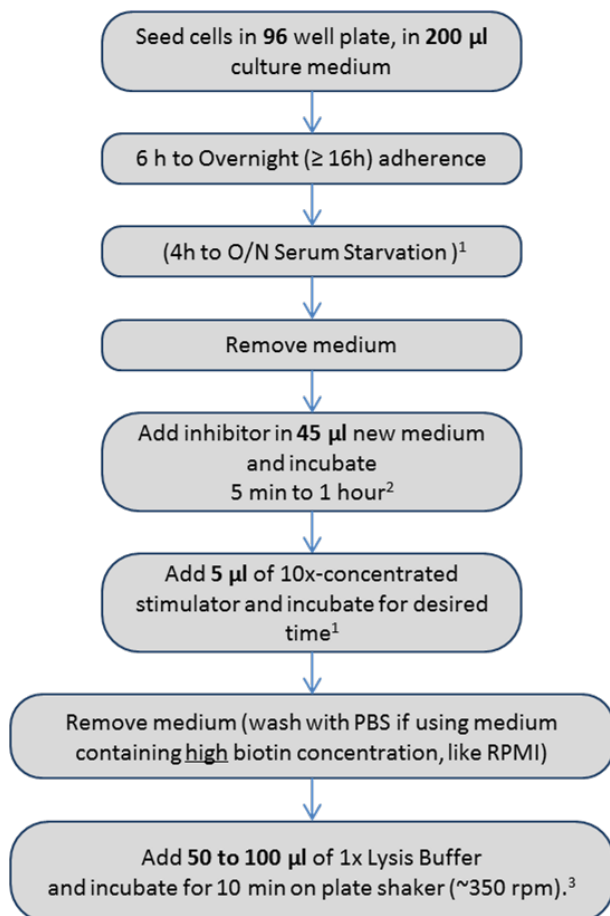
Note: Longer incubation may give greater sensitivity. Plates can be incubated overnight if required.

7. Read plate on an AlphaPlex Technology-compatible plate reader, using standard AlphaPlex settings (see above).

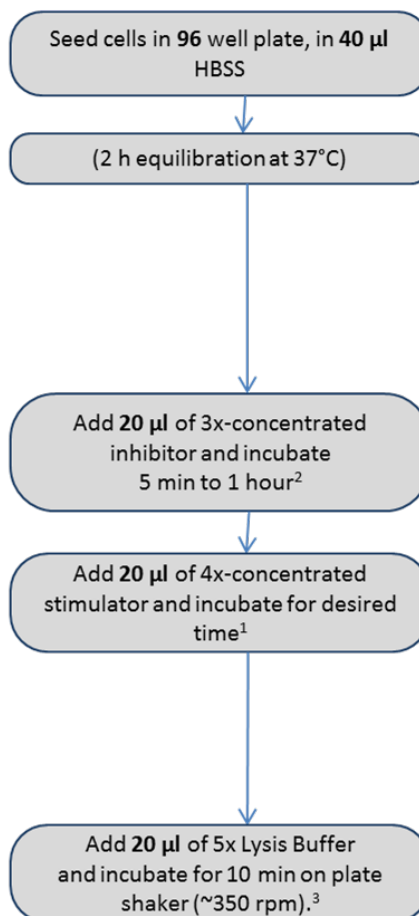
Alpha SureFire® Ultra™ HV Multiplex: 2-plates / 2-incubation assay flowchart

Dual Targets

Adherent Cells

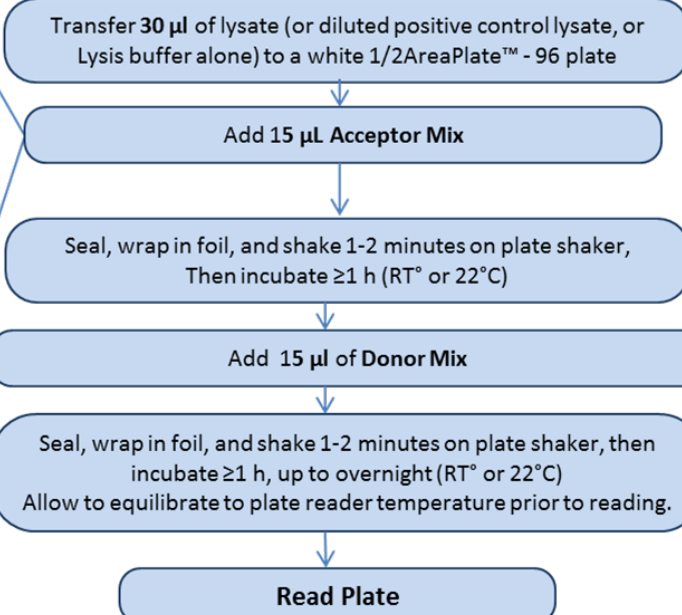


Suspension Cells



Multiplex Acceptor Mix:	Typical volumes (20 x 5 µL)	My volumes
ALSU Acceptor Mix	300	
TBSU Reaction Buffer 3	15 µL	
TBSU Reaction Buffer 4	15 µL	
CaptSure2 Alpha 545 Acceptor Beads	6 µL	

Donor Mix:	Typical volume (20 x 5 µL)	My volumes
Dilution Buffer	288 µL	
Donor Beads	12 µL	



¹ Depending on cell type and pathway analyzed.

² Depending on type of inhibitor used: 5 min is generally enough for receptor antagonists; more time is needed to block intracellular targets.

³ May stop and freeze lysates at -20°C if desired. If doing this, re-shake after thawing to ensure homogeneity of lysate before pipetting.

Dual target measurement using combined Terbium *SureFire® Ultra™ HV* Assay kit + AlphaLISA® *SureFire® Ultra™ HV* target assay kit

B. 1 Plate Assay - assay protocol for non-adherent cells, and for high-throughput applications.

Cell Seeding

1. Harvest cells by centrifugation, and re-suspend cells in HBSS at a suitable cell density. We recommend 10^7 cells/mL as a starting point. Seed 12 μ L of cells/well into a 96-well white opaque culture plate (eg 1/2AreaPlate™ - 96). Note: as engaging less cells per well can result in increased signal to background ratios, it is important to optimize this factor.
2. If using test agents/inhibitors, add 6 μ L/well of 4X inhibitors prepared in HBSS. Otherwise add 6 μ L/well of HBSS.

Note: Peptidic agonists and antagonists can often stick to plastic surfaces. To minimize this effect, dilute in HBSS containing a suitable carrier protein (e.g. 0.1% protease-free BSA).

3. Return cells to incubator at 37°C for 1-2 hours.

Cell Treatment

4. Treat cells with agonists/buffer by addition of 6 μ L/well of 4X agonist stock/buffer in HBSS containing 0.1% BSA. The final volume in the wells should be 24 μ L.

Lysate Preparation

5. To lyse the cells, add 6 μ L/well of 5X Lysis Buffer. *Add 30 μ L of Control lysates to separate wells. We recommend testing a serial dilution of Control lysate (eg 100, 50, 25, 12.5, 6.25 and 0% diluted in 1X Lysis Buffer).*

SureFire Ultra Assay

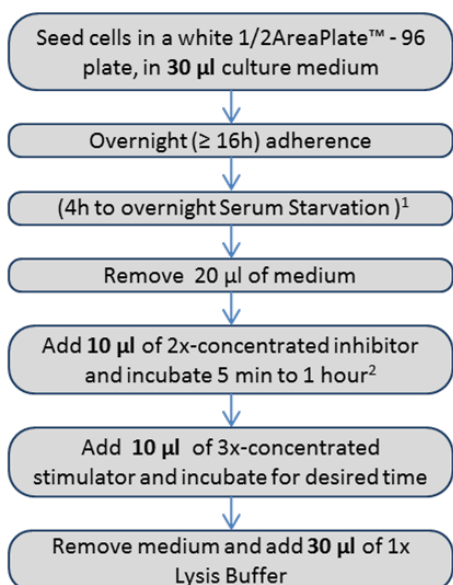
6. Add 15 μ L of Acceptor Mix to wells. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature.
7. Add 15 μ L of Donor Mix to wells under subdued light. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature in the dark.

Note: Longer incubation may give greater sensitivity. Plates can be incubated overnight if required.

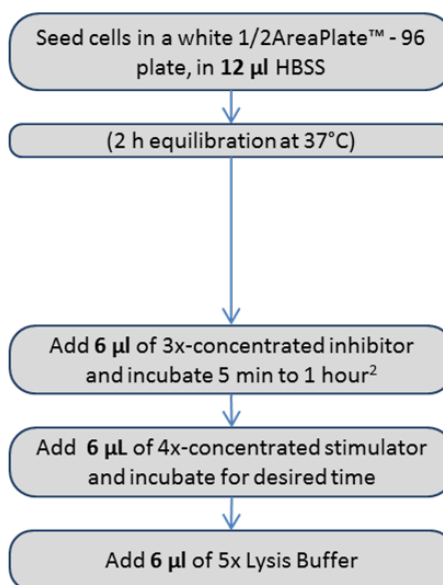
8. Read plate on an AlphaPlex Technology-compatible plate reader, using standard AlphaPlex settings (see above).

Alpha SureFire® Ultra™ HV Multiplex: 1-plate / 2-incubation assay flowchart Dual Targets

Adherent Cells



Suspension Cells



Seal and incubate for 10 min on plate shaker (~350 rpm).³

In control wells, add 30 µL positive control lysate dilution or lysis buffer alone.

Multiplex Acceptor Mix:	Typical volumes (20 x 5 µL)	My volumes
ALSU Acceptor Mix	300	
TBSU Reaction Buffer 3	15 µL	
TBSU Reaction Buffer 4	15 µL	
CaptSure2 Alpha 545 Acceptor Beads	6 µL	

Add 15 µL Acceptor Mix

Seal, wrap in foil, and shake 1-2 minutes on plate shaker, Then incubate ≥1 h (RT° or 22°C)

Donor Mix:	Typical volume (20 x 5 µL)	My volumes
Dilution Buffer	288 µL	
Donor Beads	12 µL	

Add 15 µl of Donor Mix

Seal, wrap in foil, and shake 1-2 minutes on plate shaker, then incubate ≥1 h, up to overnight (RT° or 22°C)
Allow to equilibrate to plate reader temperature prior to reading.

Read Plate

¹ Depending on cell type and pathway analyzed.

² Depending on type of inhibitor used: 5 min is generally enough for receptor antagonists; more time is needed to block intracellular targets.

³ May stop and freeze lysates at -20°C if desired. If doing this, re-shake after thawing to ensure homogeneity of lysate before pipetting.

Supplementary Buffers and Beads

If using the standard protocol, sufficient amounts of buffers and beads are provided in the kits. However in case the standard protocol would be modified, more buffers or beads may be needed. In this case, you can order additional buffers and beads using the following catalog numbers:

Item	Suggested source	Catalog #	Size
Lysis Buffer (5X) - <i>Ultra</i>	PerkinElmer Inc.	ALSU-LB-10mL	10mL
	PerkinElmer Inc.	ALSU-LB-100mL	100mL
Activation Buffer - <i>Ultra</i>	PerkinElmer Inc.	ALSU-AB-10mL	10mL
	PerkinElmer Inc.	ALSU-AB-100mL	100mL
Dilution Buffer - <i>Ultra</i>	PerkinElmer Inc.	ALSU-DB-10mL	10mL
	PerkinElmer Inc.	ALSU-DB-100mL	100mL
Alpha Streptavidin Donor Beads -2mg/mL	PerkinElmer Inc.	ALSU-ASDB-0.06mL	60µL
	PerkinElmer Inc.	ALSU-ASDB-1.2mL	1.2mL
	PerkinElmer Inc.	ALSU-ASDB-6mL	6mL

Useful Links

For FAQ and troubleshooting, please go to:

www.perkinelmer.com or www.perkinelmer.com/AlphaPlex

For a complete list of AlphaLISA *SureFire Ultra* and Alpha *SureFire Ultra* Multiplex kits, please go to:

www.perkinelmer.com/category/alpha-surefire-kits or www.tgrbio.com

For technical support please go to: www.perkinelmer.com/ASK

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

This product is not for resale or distribution except by authorized distributors.

LIMITED WARRANTY: PerkinElmer, Inc. warrants that, at the time of shipment, the products sold by it are free from defects in material and workmanship and conform to specifications which accompany the product. PerkinElmer Inc. makes no other warranty, express or implied with respect to the products, including any warranty of merchantability or fitness for any particular purpose. Notification of any breach of warranty must be made within 60 days of receipt unless provided in writing by PerkinElmer Inc. No claim shall be honored if the customer fails to notify PerkinElmer Inc. within the period specified. The sole and exclusive remedy of the customer for any liability of PerkinElmer Inc. of any kind including liability based upon warranty (express or implied whether contained herein or elsewhere), strict liability contract or otherwise is limited to the replacement of the goods or the refunds of the invoice price of goods. PerkinElmer Inc. shall not in any case be liable for special, incidental or consequential damages of any kind.

List of Available TBSU HV kits

Assay	Catalogue Numbers (100 points)
p-4E-BP1 (T37/46)	TBSU-P4EBP-A-HV
p-Akt1/2/3 (T308)	TBSU-PAKT-A-HV
p-Akt1/2/3 (S473)	TBSU-PAKT-B-HV
Total Cofilin	TBSU-TCOF-A-HV
p-CREB (S133)	TBSU-PCREB-A-HV
p-EGF Receptor (Y1068)	TBSU-PEGFR-A-HV
p-eIF2 α (S51)	TBSU-PEIF2-B-HV
p-eIF4E (S209)	TBSU-PEIF4-A-HV
p-ERK1/2 (T202/Y204)	TBSU-PERK-A-HV
p-GSK3 β (S9)	TBSU-PGS3B-A-HV
p-IGF-1 Receptor β (Y1135/1136)	TBSU-PIGFR-B-HV
p-IKK α (S176/180)	TBSU-PIKKA-A-HV
p-Insulin Receptor β (Y1150/1151)	TBSU-PINR-A-HV
p-JNK1/2/3 (T183/Y185)	TBSU-PJNK-A-HV
p-MEK1 (S218/222)	TBSU-PMEK1-A-HV
p-mTOR (S2448)	TBSU-PMTOR-A-HV
p-NF- κ B p65 (S536)	TBSU-PNFKB-A-HV
p-p38 MAPK (T180/Y182)	TBSU-PP38-B-HV
p-p70 S6K (T389)	TBSU-PP70-A-HV
p-Ribosomal Protein S6 (S240/244)	TBSU-PS6R-A-HV
p-SLP-76 (S376)	TBSU-PSLP-A-HV
p-SMAD1 (S463/465)	TBSU-PSM1-A-HV
p-SMAD3 (S423/425)	TBSU-PSM3-A-HV
p-STAT1 (S727)	TBSU-PST1-B-HV
p-STAT1 (Y701)	TBSU-PST1-A-HV
p-STAT3 (Y705)	TBSU-PST3-A-HV
p-STAT4 (Y693)	TBSU-PST4-A-HV
p-STAT5 (Y694/699)	TBSU-PST5-B-HV
p-STAT6 (Y641)	TBSU-PST6-A-HV
p-SYK (Y525/526)	TBSU-PSYK-A-HV
p-VEGF Receptor 2 (Y1175)	TBSU-PVGFR-A-HV