

my name is
Hello
my name is

Same Great
Imaging Agent

IVIS® - Meet Your Reagents
IVISense™ and IVISbrite™

WE'VE RENAMED

Our Fluorescent & Bioluminescent Reagents

PerkinElmer's *in vivo* optical imaging reagents have been renamed to **IVISense** or **IVISbrite** – aligning better with our industry leading IVIS optical imaging instruments.

In addition, our new reagent names are more descriptive, helping you to easily select the right reagents for your imaging applications.

Q Have formulations of the products changed?

A **No.** The changes are in name only. Our *in vivo* imaging reagents maintain the same high-quality formulation and exquisite functionality as before, just as we maintain our commitment to stand behind our *in vivo* imaging reagents and support you in designing and executing your experiments.

Q Have part numbers changed?

A **No.** The changes are in name only. You can still order fluorescent and bioluminescent reagents using the same part numbers for easy reference.

Q The name on the vial label differs from the name on the package. How can I be sure a mistake was not made in preparing my order?

A You can verify that your order is correct by checking the part number and lot number on the vial label against the numbers on the package. If they match, then you can be confident that you have the correct reagents.

The **IVISense** brand represents our wide range of **fluorescent** probes, labels, and dyes.

The **IVISbrite** brand represents our **bioluminescent** reagents including tumor and bacterial cell lines, substrates, and lentiviral particles.

Fluorescent Agent Portfolio

Fluorescent Probes (Activatable)		
Cat #	Former Name	New Name
NEV11112	Cat B 680 FAST Fluorescent Imaging Agent	IVISense Cat B 680 FAST Fluorescent Probe
NEV11098	Cat B 750 FAST Fluorescent Imaging Agent	IVISense Cat B 750 FAST Fluorescent Probe
NEV11000	Cat K 680 FAST Fluorescent Imaging Agent	IVISense Cat K 680 FAST Fluorescent Probe
NEV10100	MMPsense 645 FAST Fluorescent Imaging Agent	IVISense MMP 645 FAST Fluorescent Probe
NEV10126	MMPsense 680 Fluorescent Imaging Agent	IVISense MMP 680 Fluorescent Probe
NEV10168	MMPsense 750 FAST Fluorescent Imaging Agent	IVISense MMP 750 FAST Fluorescent Probe
NEV11169	Neutrophil Elastase 680 FAST Fluorescent Imaging Agent	IVISense Neutrophil Elastase 680 FAST Fluorescent Probe
NEV10003	ProSense 680 Fluorescent Imaging Agent	IVISense Pan Cathepsin 680 Fluorescent Probe
NEV10001EX	ProSense 750 EX Fluorescent Imaging Agent	IVISense Pan Cathepsin 750 Fluorescent Probe
NEV11171	ProSense 750 FAST Fluorescent Imaging Agent	IVISense Pan Cathepsin 750 FAST Fluorescent Probe
NEV11079	ReninSense 680 FAST Fluorescent Imaging Agent	IVISense Renin 680 FAST Fluorescent Probe
Fluorescent Probes (Targeted)		
Cat #	Former Name	New Name
NEV11053	Annexin Vivo 750 Fluorescent Imaging Agent	IVISense Annexin-V 750 Fluorescent Probe
NEV10090	BombesinRSense 680 Fluorescent Imaging Agent	IVISense Bombesin Receptor 680 Fluorescent Probe
NEV10040	FolateRSense 680 Fluorescent Imaging Agent	IVISense Folate Receptor 680 Fluorescent Probe
NEV11070	HypoxiSense 680 Fluorescent Imaging Agent	IVISense Hypoxia CA IX 680 Fluorescent Probe
NEV10640	IntegriSense 645 Fluorescent Imaging Agent	IVISense Integrin Receptor 645 Fluorescent Probe
NEV10645	IntegriSense 680 Fluorescent Imaging Agent	IVISense Integrin Receptor 680 Fluorescent Probe
NEV10873	IntegriSense 750 Fluorescent Imaging Agent	IVISense Integrin Receptor 750 Fluorescent Probe
NEV10020EX	OsteoSense 680EX - 24 nmol Fluorescent Imaging Agent	IVISense Osteo 680 Fluorescent Probe
NEV10053EX	OsteoSense 750EX - 48 nmol Fluorescent Imaging Agent	IVISense Osteo 750 Fluorescent Probe
NEV11105	OsteoSense 800 Fluorescent Imaging Agent	IVISense Osteo 800 Fluorescent Probe
NEV10060	TLectinSense 680 Fluorescent Imaging Agent	IVISense Tomato Lectin 680 Fluorescent Probe
NEV10091	Transferrin-Vivo 750 Fluorescent Imaging Agent	IVISense Transferrin Receptor 750 Fluorescent Probe
760561	XenoLight RediJect 2-DG 750 Probe Explorer kit	IVISense 2-DG 750 Fluorescent Probe in RediJect™ Solution (1 vial)
760562	XenoLight RediJect 2-DG 750 Probe Standard kit	IVISense 2-DG 750 Fluorescent Probe in RediJect Solution (4 vials)
760567	XenoLight RediJect 2-DG 750 Control Dye	IVISense 2-DG 750 Fluorescent Control Dye in RediJect Solution (1 vial)
133397	XenoLight Rediject Bacterial Detection Probe Explorer kit	IVISense Bacterial 750 Fluorescent Probe in RediJect Solution (1 vial)
133398	XenoLight Rediject Bacterial Detection Probe Standard Kit	IVISense Bacterial 750 Fluorescent Probe in RediJect Solution (4 vials)
133399	XenoLight Rediject Bacterial Detection Probe Control Dye	IVISense Bacterial 750 Fluorescent Control Dye in RediJect Solution (1 vial)
Fluorescent Probes (Vascular)		
Cat #	Former Name	New Name
NEV10054EX	AngioSense 680 EX Fluorescent Imaging Agent	IVISense Vascular 680 Fluorescent Probe
NEV10011EX	AngioSense 750 EX Fluorescent Imaging Agent	IVISense Vascular 750 Fluorescent Probe
NEV10149	AngioSPARK 680	IVISense Vascular NP 680 Fluorescent Nanoparticles
NEV10150	AngioSPARK 750	IVISense Vascular NP 750 Fluorescent Nanoparticles
NEV11121	GastroSense 750 Fluorescent Imaging Agent	IVISense Gastrointestinal 750 Fluorescent Probe
NEV10117	Genhance 680 1mg Fluorescent Imaging Agent	IVISense Acute Vascular 680 Fluorescent Probe (1 mg)
NEV10130	Genhance 680 5mg Fluorescent Imaging Agent	IVISense Acute Vascular 680 Fluorescent Probe (5 mg)

Fluorescent Probes (Vascular) - Continued		
Cat #	Former Name	New Name
NEV10118	Genhance 750 1mg Fluorescent Imaging Agent	IVISense Acute Vascular 750 Fluorescent Probe (1 mg)
NEV10177	Genhance 750 5mg Fluorescent Imaging Agent	IVISense Acute Vascular 750 Fluorescent Probe (5 mg)
NEV30000	GFR-Vivo 680 Fluorescent Imaging Agent	IVISense GFR 680 Fluorescent Probe
NEV10116	Superhance 680 Fluorescent Imaging Agent	IVISense Edema 680 Fluorescent Probe

Fluorescent Panels		
Cat #	Former Name	New Name
NEV20003	Fluorescent Imaging Agent Panel – Arthritis	IVISense Arthritis Fluorescent Imaging Panel
NEV20013	Fluorescent Imaging Agent Panel – Inflammation	IVISense Inflammation Fluorescent Imaging Panel
NEV20005	Fluorescent Imaging Agent Panel – Oncology	IVISense Oncology Fluorescent Imaging Panel
NEV20016	Fluorescent Imaging Agent Panel – Toxicology	IVISense Toxicology Fluorescent Imaging Panel
NEV20015	Fluorescent Imaging Agent Panel – Tumor Metabolism	IVISense Tumor Metabolism Fluorescent Imaging Panel
NEV20014	Fluorescent Imaging Agent Panel – Vascular	IVISense Vascular Fluorescent Imaging Panel
NEV20011	Fluorescent Imaging Agent Panel – Featured Probes Pack	IVISense Featured Probes Pack Fluorescent Imaging Panel
NEV20000	Fluorescent Imaging Agent Panel – Sample Pack	IVISense Sample Pack Fluorescent Imaging Panel

Fluorescent Dyes (645, 680, & 800)		
Cat #	Former Name	New Name
NEV11173	VivoTag 645 (1 mg)	IVISense 645 NHS Fluorescent Dye (1 mg)
NEV11174	VivoTag 645 (5 mg)	IVISense 645 NHS Fluorescent Dye (5 mg)
NEV11273	VivoTag 645-MAL (1 mg)	IVISense 645 MAL Fluorescent Dye (1 mg)
NEV11274	VivoTag 645-MAL (5 mg)	IVISense 645 MAL Fluorescent Dye (5 mg)
NEV11119	VivoTag 680 XL (1mg)	IVISense 680 NHS Fluorescent Dye (1 mg)
NEV11120	VivoTag 680 XL (5 mg)	IVISense 680 NHS Fluorescent Dye (5 mg)
NEV11219	VivoTag 680XL-MAL (1 mg)	IVISense 680 MAL Fluorescent Dye (1 mg)
NEV11220	VivoTag 680XL-MAL (5 mg)	IVISense 680 MAL Fluorescent Dye (5 mg)
NEV11107	VivoTag 800 (1 mg)	IVISense 800 NHS Fluorescent Dye (1 mg)
NEV11108	VivoTag 800 (5 mg)	IVISense 800 NHS Fluorescent Dye (5 mg)

Fluorescent Dyes (680 & 750 Self-Quenching)		
Cat #	Former Name	New Name
NEV10121	VivoTag-S 680 (1 mg)	IVISense 680 NHS Fluorescent Self-Quenching Dye (1 mg)
NEV10122	VivoTag-S 680 (5 mg)	IVISense 680 NHS Fluorescent Self-Quenching Dye (5 mg)
NEV10123	VivoTag-S 750 (1 mg)	IVISense 750 NHS Fluorescent Self-Quenching Dye (1 mg)
NEV10124	VivoTag-S 750 (5 mg)	IVISense 750 NHS Fluorescent Self-Quenching Dye (5 mg)
NEV11223	VivoTag-S 750 MAL (1 mg)	IVISense 750 MAL Fluorescent Self-Quenching Dye (1 mg)
NEV11224	VivoTag-S 750 MAL (5 mg)	IVISense 750 MAL Fluorescent Self-Quenching Dye (5 mg)

Fluorescent Labeling Kit		
Cat #	Former Name	New Name
NEV11118	VivoTag 680XL Protein Labeling Kit	IVISense 680 NHS Fluorescent Labeling Kit

Fluorescent Cell Labeling Dyes		
Cat #	Former Name	New Name
NEV12001	VivoTrack 680 (Sample Kit)	IVISense 680 Fluorescent Cell Labeling Dye (1 x 0.2mg)
NEV12000	VivoTrack 680 (Standard Kit)	IVISense 680 Fluorescent Cell Labeling Dye (5 x 0.2mg)
125964	XenoLight DiR	IVISense DiR 750 Fluorescent Cell Labeling Dye

Bioluminescent Reagent Portfolio

Bioluminescent Substrates		
Powders		
Cat #	Former Name	New Name
122799	XenoLight D-Luciferin Pottasium salt (1 g)	IVISbrite D-Luciferin Potassium Salt Bioluminescent Substrate (1g)
122799-5	XenoLight D-Luciferin Pottasium salt (5 g)	IVISbrite D-Luciferin Potassium Salt Bioluminescent Substrate (5 x 1g)
122799-10	XenoLight D-Luciferin Pottasium salt (10 g)	IVISbrite D-Luciferin Potassium Salt Bioluminescent Substrate (10 x 1g)
Injectables		
Cat #	Former Name	New Name
770504	XenoLight Rediject D-Luciferin (50 Injections)	IVISbrite D-Luciferin Bioluminescent Substrate in RediJect Solution
770505	XenoLight Rediject D-Luciferin Ultra (50 Injections)	IVISbrite D-Luciferin Ultra Bioluminescent Substrate in RediJect Solution
760506	XenoLight Rediject Coelenterazine h (50 Injections)	IVISbrite Coelenterazine h Bioluminescent Substrate in RediJect Solution
760535	XenoLight RediJect Inflammation probe Explorer kit	IVISbrite MPO 425 Chemiluminescent Probe in RediJect Solution (1 vial)
760536	XenoLight RediJect Inflammation probe Standard kit	IVISbrite MPO 425 Chemiluminescent Probe in RediJect Solution (4 vials)
Lentiviral Particles		
Cat #	Former Name	New Name
CLS960002	RediFect Red-Fluc-Puromycin	IVISbrite Red F-luc-Puromycin Lentiviral Particles
CLS960003	RediFect Red-Fluc-GFP	IVISbrite Red F-luc-GFP Lentiviral Particles
CLS960004	RediFect Green Renilla-Puromycin	IVISbrite Green Renilla-Puromycin Lentiviral Particles
Tumor Cell Lines		
Cat #	Former Name	New Name
BW128090	Bioware Brite 4T1-Red-FLuc-GFP	IVISbrite 4T1 Red F-luc-GFP Bioluminescent Tumor Cell line
BW124087	Bioware Brite 4T1-Red-FLuc	IVISbrite 4T1 Red F-luc Bioluminescent Tumor Cell line
BW119266	Bioware Brite A549-Red-Fluc	IVISbrite A549 Red F-luc Bioluminescent Tumor Cell line
BW124734	Bioware Brite B16F10-Red-FLuc	IVISbrite B16F10 Red F-luc Bioluminescent Tumor Cell line
BW125058	Bioware Brite BxPC3-Red-FLuc	IVISbrite BxPC3 Red F-luc Bioluminescent Tumor Cell line
BW124317	Bioware Brite Colo205-Red-FLuc	IVISbrite Colo205 Red F-luc Bioluminescent Tumor Cell line
BW134246	Bioware Brite GL261-Red-FLuc	IVISbrite GL261 Red F-luc Bioluminescent Tumor Cell line
BW124318	Bioware Brite HCT-116-Red-FLuc	IVISbrite HCT-116 Red F-luc Bioluminescent Tumor Cell line
BW134280	Bioware Brite HepG2-Red-FLuc	IVISbrite HepG2 Red F-luc Bioluminescent Tumor Cell line
BW128092	Bioware Brite HT1080-Red-FLuc	IVISbrite HT1080 Red F-luc Bioluminescent Tumor Cell line
BW124353	Bioware Brite HT-29-Red-FLuc	IVISbrite HT-29 Red F-luc Bioluminescent Tumor Cell line
BW124735	Bioware Brite K562-Red-FLuc	IVISbrite K562 Red F-luc Bioluminescent Tumor Cell line
BW119267	Bioware Brite LL/2-Red-FLuc	IVISbrite LL/2 Red F-luc Bioluminescent Tumor Cell line
BW125055	Bioware Brite LNCaP-Red-FLuc	IVISbrite LNCaP Red F-luc Bioluminescent Tumor Cell line
BW119262	Bioware Brite MCF7-Red-FLuc	IVISbrite MCF-7 Red F-luc Bioluminescent Tumor Cell line
BW124316	Bioware Brite NCI-H460-Red-FLuc	IVISbrite NCI-H460 Red F-luc Bioluminescent Tumor Cell line
BW128444	Bioware Brite PC3-Red-FLuc	IVISbrite PC3 Red F-luc Bioluminescent Tumor Cell line
BW133416	Bioware Brite PC3-Red-FLuc-GFP	IVISbrite PC3 Red F-luc-GFP Bioluminescent Tumor Cell line
BW119276	Bioware Brite SKOV3-Red-FLuc	IVISbrite SKOV3 Red F-luc Bioluminescent Tumor Cell line
BW124577	Bioware Brite U87MG-Red-FLuc	IVISbrite U87MG Red F-luc Bioluminescent Tumor Cell line

Bacterial Cell Lines		
Cat #	Former Name	New Name
119223	Xen14 <i>Escherichia coli</i>	IVISbrite <i>Escherichia coli</i> Xen14 Bioluminescent Bacterial Strain
119237	Xen19 <i>Listeria monocytogene</i>	IVISbrite <i>Listeria monocytogene</i> Xen19 Bioluminescent Bacterial Strain
119236	Xen44 <i>Proteus mirabilis</i>	IVISbrite <i>Proteus mirabilis</i> Xen44 Bioluminescent Bacterial Strain
119228	Xen05 <i>Pseudomonas aeruginosa</i>	IVISbrite <i>Pseudomonas aeruginosa</i> Xen05 Bioluminescent Bacterial Strain
119229	Xen41 <i>Pseudomonas aeruginosa</i>	IVISbrite <i>Pseudomonas aeruginosa</i> Xen41 Bioluminescent Bacterial Strain
119235	Xen33 <i>Salmonella typhimunium</i>	IVISbrite <i>Salmonella typhimunium</i> Xen33 Bioluminescent Bacterial Strain
119239	Xen08.1 <i>Staphylococcus aureus</i>	IVISbrite <i>Staphylococcus aureus</i> Xen8.1 Bioluminescent Bacterial Strain
119240	Xen29 <i>Staphylococcus aureus</i>	IVISbrite <i>Staphylococcus aureus</i> Xen29 Bioluminescent Bacterial Strain
119242	Xen31 <i>Staphylococcus aureus</i>	IVISbrite <i>Staphylococcus aureus</i> Xen31 Bioluminescent Bacterial Strain
119243	Xen36 <i>Staphylococcus aureus</i>	IVISbrite <i>Staphylococcus aureus</i> Xen36 Bioluminescent Bacterial Strain
119244	Xen40 <i>Staphylococcus aureus</i>	IVISbrite <i>Staphylococcus aureus</i> Xen40 Bioluminescent Bacterial Strain
119231	Xen27 <i>Shigella dysenteriae</i>	IVISbrite <i>Shigella dysenteriae</i> Xen27 Bioluminescent Bacterial Strain
119232	Xen24 <i>Yersinia enterocolitica</i>	IVISbrite <i>Yersinia enterocolitic</i> Xen24 Bioluminescent Bacterial Strain
119233	Xen25 <i>Yersinia enterocolitica</i>	IVISbrite <i>Yersinia enterocolitic</i> Xen25 Bioluminescent Bacterial Strain

HAVE MORE QUESTIONS?

Contact Global Tech Support

Global.TechSupport@perkinelmer.com

Visit Us

www.perkinelmer.com/category/in-vivo-imaging-reagents

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2021, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.