

NEW ASSAYS AND REAGENTS



NEW PRODUCTS AND APPLICATIONS TO MEET YOUR NEEDS

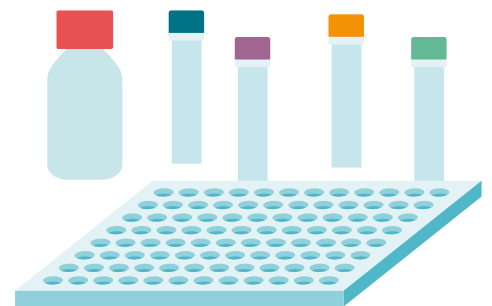
Here is our up-to-date listing of new products from the combined PerkinElmer-Cisbio reagent portfolio.

All assays are designed for small-scale studies up to large, full-library screens.

Please contact your local account manager to check availability, or to learn more.

Get faster, more reliable results using the many assay technologies we have available

- HTRF®
- AlphaLISA®
- AlphaLISA® *SureFire® Ultra™*



All kits are ready-to-use, plate-based solutions.

NEW PRODUCTS



NEW PRODUCTS LAUNCHES JANUARY 2021

BIOMARKERS ASSAYS

AlphaLISA® PRODUCTS **PART- 500 tests**

Glucagon Detection kit	AL3140C
SARS-CoV2 Spike S1 kit	AL3142C

HTRF® PRODUCTS **PART- 500 tests**

HIV p24 kit	64P24PEG
HO-1 kit	64HOX1PEG

CELLULAR PHOSPHO-PROTEIN ASSAYS

AlphaLISA® SureFire® Ultra™ PRODUCTS **PART- 500 tests**

Phospho RSK1 (Thr359/Ser363) Detection kit	ALSU-PRSK1-B500
Phospho RSK1 (Ser380) Detection kit	ALSU-PRSK1-C500
Phospho MLKL (Ser345) mouse Detection kit	ALSU-PMLKL-A500
Total MLKL (mouse) Detection kit	ALSU-TMLKL-B500
Phospho RSK1 (Thr573) Detection kit	ALSU-PRSK1-A500
Total RSK1 Detection kit	ALSU-TRSK1-A500
Phospho DAP12 (pan Tyr) Detection kit	ALSU-PDAP12-A500

HTRF® PRODUCTS **PART- 500 tests**

Cereblon total kit - Standard	64CRBNCDA
Insulin receptor B kit	63ADK016PEG
RIPK1 phospho S166 kit	64RIPK1S1PEG
RIPK1 total kit	64RIPK1TPEG

GPCRS BINDING ASSAYS

HTRF® PRODUCTS **PART- 500 tests**

AP2 total kit	64AP2TPEB
B-Arrestin 1 total kit	64BAR1TPEB
B-Arrestin 2 recruitment kit	62BDBAR2PEB
B-arrestin 2 total kit	64BAR2TPEB
GTP Gi binding kit	62GTPPET
Stabilization buffer 1 - 200 mL	62SB1FDF
Stimulation buffer 4	64SB4FDF
Wash buffer 1 - 200 mL	62WB1FDF

TAG-LITE ASSAYS

HTRF® PRODUCTS **PART- 200 tests**

pSNAP-AT2 Angiotensin receptor (Tag-lite plasmid) kit	PSNAPAT2
pSNAP-HTR4 Serotonin receptor (Tag-lite plasmid) kit	PSNAPHTR4

PROTEIN PROTEIN INTERACTION ASSAYS

HTRF® PRODUCTS **PART- 500 tests**

KRAS G12C GTP binding kit	64BDKRASGPEG
KRAS G12C/SOS1 binding kit	64KRASG12PEG
KRAS WT GTP binding kit	64BDKRASWPEG
KRAS WT/SOS1 binding kit	64KRASWTPEG

TOOLBOX

AlphaScreen PRODUCTS **PART- 1mg**

Anti-6xHis Alpha Donor Beads	AS116D
------------------------------	--------

HTRF products also available under 10 000 tests & bulk size

AlphaLISA® SureFire® Ultra™ products also available under 100 - 10 000 - 50 000 tests



NEW PRODUCTS LAUNCHES MARCH 2021

BIOMARKERS ASSAYS

AlphaLISA® PRODUCTS	PART
AlphaLISA alpha-SMA Kit	AL3143C
AlphaLISA IP-One kit	AL3145D

HTRF® PRODUCTS	PART
Human Amyloid b 1-42 Detection kit	62B42PEG
Protein Disaggregation Detection kit	64DAGGRPET

PROTEIN-PROTEIN INTERACTION ASSAYS

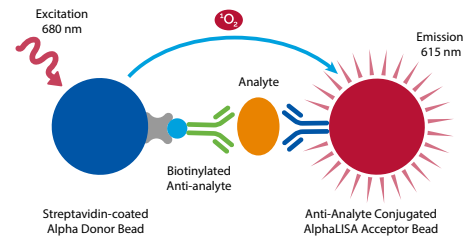
AlphaLISA® PRODUCTS	PART
AlphaLISA STING WT Binding Kit	AL3144C

CELLULAR PHOSPHO-PROTEIN ASSAYS

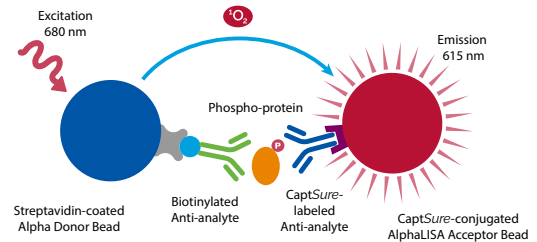
AlphaLISA® SureFire® Ultra™ PRODUCTS	PART
Beclin-1 Total Detection kit	ALSU-TBEC-A500
CHK2 pT68 Detection kit	ALSU-PCHK2-A500
CHK2 Total Detection kit	ALSU-TCHK2-A500
IRF3 pS386 Detection kit	ALSU-PIRF3-A500
IRF3 Total Detection kit	ALSU-TIRF3-A500

HTRF® PRODUCTS	PART
GAPDH Housekeeping Detection kit	64GAPDHPEG
STING pS365 (mouse) Detection kit	64MSTGPEG
STING total (mouse) Detection kit	64NMSTGPEG

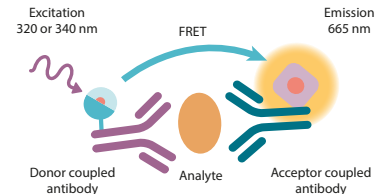
AlphaLISA® technology



AlphaLISA® SureFire® Ultra™ technology



HTRF® technology



HTRF products also available under 10 000 tests & bulk size

AlphaLISA® SureFire® Ultra™ products also available under 100 - 10 000 - 50 000 tests

AlphaLISA® products are also available as 100, 5000 tests & bulk sizes

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86 21 5018 9880 Japan +81-(0)43-306-8712

Visit www.cisbio.com and www.perkinelmer.com



cisbio