

NEW ASSAYS AND REAGENTS



NEW PRODUCTS AND APPLICATIONS TO MEET YOUR NEEDS

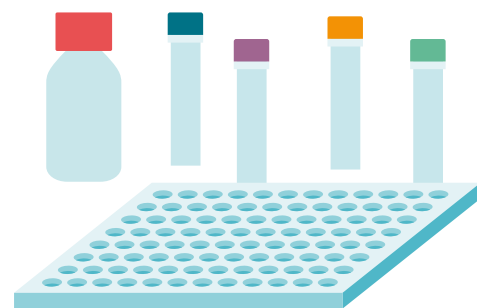
Here is our up-to-date listing of new products from the combined PerkinElmer-Cisbio reagent portfolio.

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

- DELFIA®,
- LANCE® *Ultra*™
- HTRF®
- AlphaLISA®
- CETSA®



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

NEW PRODUCTS



CELLULAR PHOSPHO-PROTEIN ASSAYS

HTRF® PRODUCTS

PART- 500 tests

FGFR1 phospho	64FGFR1Y6PEG
FGFR1 total	64FGFR1TPEG
FGFR2 phospho	64FGFR2Y6PEG
FGFR2 total	64FGFR2TPEG
FGFR3 phospho	64FGFR3Y6PEG
FGFR3 total	64FGFR3TPEG
FGFR4 phospho	64FGFR4Y6PEG
FGFR4 total	64FGFR4TPEG
Rb phospho S780	64RBS780PEG
Rb phospho S807/811	64RBS807PEG
Rb total	64RBTPEG
SMAD2 phospho S465/467	64SMAD2S5PEG
SMAD2 total	64SMAD2TPEG
Ubiquitin phospho S65	64UBIS65PEG
VEGFR2 total	64VEGFR2TPEG

PROTEIN PROTEIN INTERACTION ASSAYS

HTRF® PRODUCTS

PART- 500 tests

IL12/IL12RB1	64BDIL12PEG
IL23/IL12RB1	64BDPIL12PEG
IL23/IL23R	64BDPIL23PEG
IL6/IL6R	64BDIL6PEG
LAG3/MHCII	64LAG3PEG

CELLULAR THERMAL SHIFT ASSAYS

Alpha SureFire® CETSA® PRODUCTS

PART- 500 tests

STING	CETSA-STING-A500
-------	------------------

VIROLOGY ASSAYS

HTRF® PRODUCTS

PART- 500 tests

S1/ACE2	64LAG3PEG
SARS-Cov2 capsid kit	63ADK113PEG
SARS-Cov2 Spike S1	63ADK114PEG

ALPHALISA ASSAYS

AlphaLISA® PRODUCTS

PART- 500 tests

Cereblon Total	AL3139C
Phospho-Tau (T181)	AL3136C

HTRF products also available under 10 000 tests & bulk size

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

Alpha SureFire® CETSA® also available under 5,000 tests

CELLULAR PHOSPHO-PROTEIN ASSAYS

AlphaLISA® SureFire® Ultra™ PRODUCTS

PART- 500 tests

Phospho HDAC4 (S246)	ALSU-PHDAC-A500
Phospho Histone H2A.X (S139)	ALSU-PH2X-A500
Phospho MLKL (S358)	ALSU-PMLKL-B500
Phospho RET (panY)	ALSU-PRET-A500
Phospho RIPK1 (S166)	ALSU-PRIPK1-A500
Phospho RPS6 (S235/236)	ALSU-PS6R-B500
Phospho SLP76 (Y113)	ALSU-PSLP-B500
Phospho SLP76 (Y145)	ALSU-PSLP-C500
Phospho-PI3 Kinase p85 (Tyr458) (Mouse)	ALSU-PP13K-A500
Phospho-PLC G1 (S1248)	ALSU-PPLCG1-B500
Phospho-PLC G1 (Y1253)	ALSU-PPLCG1-C500
Phospho-PLC G1 (Y783)	ALSU-PPLCG1-A500
Phospho-PLC G2 (Y1197)	ALSU-PPLCG2-C500
Phospho-PLC G2 (Y1217)	ALSU-PPLCG2-D500
Phospho-PLC G2 (Y753)	ALSU-PPLCG2-A500
Phospho-PLC G2 (Y759)	ALSU-PPLCG2-B500
Total DAP12	ALSU-TDAP12-A500
Total H2A.X	ALSU-TH2X-A500
Total HDAC4	ALSU-THDAC-A500
Total IKKα	ALSU-TIKKA-A500
Total IRAK4	ALSU-TIRAK-A500
Total MLKL	ALSU-TMLKL-A500
Total PI3 Kinase p85	ALSU-TP13K-A500
Total PLC G1	ALSU-TPLCG1-A500
Total PLC G2	ALSU-TPLCG2-A500
Total RET	ALSU-TRET-A500
Total RIPK1	ALSU-TRIPK1-A500
Total RPS6	ALSU-TS6R-A500

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

Phospho (S235/236)/Total S6 RP Multiplex	MPSU-PTRPS6-B500
Phospho (S240/S244)/Total S6 RP Multiplex	MPSU-PTRPS6-B500
Phospho (S780)/Total Rb Multiplex	MPSU-PTRB-N500
Phospho (S807/811)/Total Rb Multiplex	MPSU-PTRB-M500
Phospho (T821/826)/Total Rb Multiplex	MPSU-PTRB-N500
p-STAT1 (Ser727) (Eu) + Total STAT1 (Tb)	MPSU-PTST1-M500
p-STAT1 (Tyr701) (Eu) + Total STAT1 (Tb)	MPSU-PTST1-K500

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