

# NEW ASSAYS AND REAGENTS



## NEW PRODUCTS AND APPLICATIONS TO MEET YOUR NEEDS

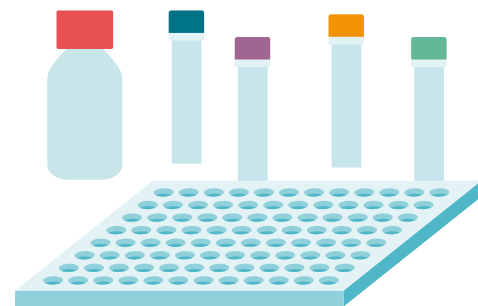
Here is our up-to-date listing of new products from the combined PerkinElmer-Cisbio reagent portfolio.

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®



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

NEW PRODUCTS



## CELLULAR PHOSPHO-PROTEIN ASSAYS

### HTRF® PRODUCTS PART- 500 tests

|                        |              |
|------------------------|--------------|
| SYK phospho-Tyr525/526 | 64SYKY525PEG |
| SYK total              | 64SYKTPEG    |

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

|                     |                 |
|---------------------|-----------------|
| ACC total           | ALSU-TACC-A500  |
| CDK2 total          | ALSU-TCDK2-A500 |
| JAK2 total          | ALSU-TJAK2-A500 |
| Raf1 phospho-Ser338 | ALSU-PRAF-A500  |
| TBK1 phospho-Ser172 | ALSU-PTBK1-A500 |
| TBK1 total          | ALSU-TTBK1-A500 |

## BIOCHEMICAL PROTEIN-PROTEIN INTERACTION ASSAYS

### HTRF® PRODUCTS PART- 500 tests

|                               |          |
|-------------------------------|----------|
| Human PD1 / PD-L1 binding kit | 64PD1PEG |
|-------------------------------|----------|

### AlphaLISA® PRODUCTS PART- 500 tests

|                     |         |
|---------------------|---------|
| Human CD28/CD86 kit | AL3131C |
| Cyno PD1/PDL1 kit   | AL583C  |

## BIOMARKERS ASSAYS

### HTRF® PRODUCTS PART- 500 tests

|                  |           |
|------------------|-----------|
| FcRn binding kit | 64FCRNPET |
|------------------|-----------|

### AlphaLISA® PRODUCTS PART- 500 tests

|                 |        |
|-----------------|--------|
| Mouse TREM2 kit | AL591C |
|-----------------|--------|

## CYTOKINE ASSAYS

### AlphaLISA® PRODUCTS PART- 500 tests

|                        |         |
|------------------------|---------|
| Human IL3 kit          | AL3134C |
| Mouse IL3 kit          | AL587C  |
| Mouse IFN $\gamma$ kit | AL586C  |

## TAG-LITE REAGENTS

### HTRF® PRODUCTS PART- 200 tests

|   |          |
|---|----------|
| Delta OpioidR expressing cells labeled with Terbium | C1TT1DOP |
| Kappa OpioidR expressing cells labeled with Terbium | C1TT1KOP |
| Mu OpioidR expressing cells labeled with Terbium    | C1TT1MOP |
| Opioid Receptor red antagonist - fluorescent probe  | L0005RED |

## BIOLOGICS BINDING ASSAYS

### HTRF® PRODUCTS PART- 200 tests

|   |            |
|---|------------|
| Mouse Hybridoma binding (GST antigen) kit | 64HYGSTPEG |
| Mouse Hybridoma binding (HIS antigen) kit | 64HYHISPEG |

## 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](http://www.cisbio.com) and [www.perkinelmer.com](http://www.perkinelmer.com)



cisbio

*HTRF products also available under 10 000 tests & bulk size  
AlphaLISA® SureFire® Ultra™ products also available under 10 000 - 50 000  
& 100 000 tests  
AlphaLISA® products also available under 100 - 5 000 tests & bulk size*