

# NEW HTRF® ASSAYS AND REAGENTS



## NEW HTRF® PRODUCTS AND APPLICATIONS TO MEET YOUR NEEDS

Here is our up-to-date listing of new products.

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 if the assay you are looking for is not in this list.

### **Get faster, more reliable results using HTRF assays**

- More consistent and reliable data from every well with physiological screening environment
- Greater specificity and high Z' scores for robust, reproducible results
- Fewer false negatives and repeat runs
- Adapt easily to all throughput requirements
- Simple wash-free mix-and-read protocol



HTRF kits are ready-to-use and include detection reagents, all necessary buffers and controls.

NEW PRODUCTS



## CELLULAR PHOSPHO-PROTEIN ASSAYS

PRODUCT	PART#
ACC total kit - 500 tests	64NACPEG
MKK4 phospho-S257 kit - 500 tests	64MK4CPEG
MKK4 total kit - 500 tests	64NK4CPEG
RB phospho-S780 kit - 500 tests	63ADK104PEG
RB phospho-S807/811 kit - 500 tests	63ADK105PEG
RB total kit - 500 tests	63ADK106PEG

10 000 tests & bulk size available

## CYTOKINE ASSAYS

PRODUCT	PART#
Human IL13 kit - 500 tests	62HIL13PEG
Human IL22 kit - 500 tests	62HIL22PEG

10 000 tests & bulk size available

## BIOMARKER ASSAY

PRODUCT	PART#
HMGB1 kit - 500 tests	62HMGPEG

10 000 tests & bulk size available

## BIOCHEMICAL PROTEIN-PROTEIN INTERACTION ASSAYS

PRODUCT	PART#
Human H232 STING binding kit	64BDSTGHPEG
Human AQ STING binding kit	64BDSTGQPEG
Human STING WT binding kit t	64BDSTGPEG

10 000 tests & bulk size available

## RECEPTOR BINDING ASSAYS - TAG-LITE®

PRODUCT	PART#
Adenosine A1 receptor labeled cells - 200 tests	C1TT1A1
Adenosine A2B receptor labeled cells - 200 tests	C1TT1A2B
Adenosine A3 receptor labeled cells - 200 tests	C1TT1A3
Adenosine A1 receptor red antagonist - 5 000 tests	L0067RED
Adenosine A2B receptor red antagonist - 5 000 tests	L0068RED
Adenosine A3 receptor green antagonist - 5 000 tests	L0069GRE
Plasmid Adenosine A1 receptor	PSNAPA1
Plasmid Adenosine A2B receptor	PSNAPA2B
Plasmid Adenosine A3 receptor	PSNAPA3
ETB stable cell line	C1SU1ETB

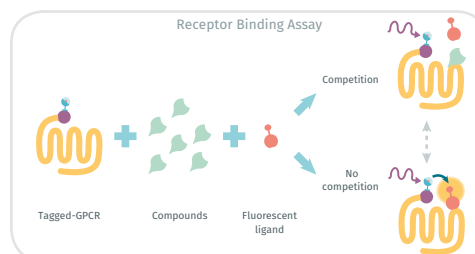
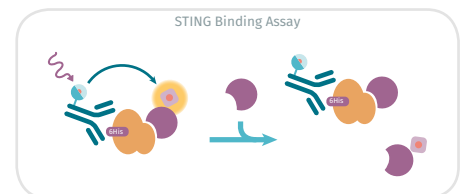
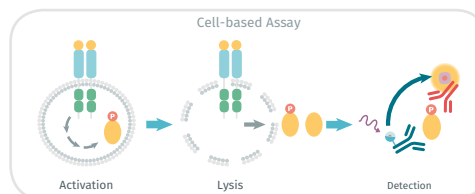
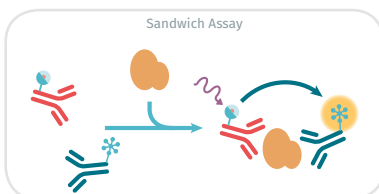
## PLATES AND TOP SEAL

PRODUCT	PART#
CulturPlate-96, White Opaque 96-well Microplate, Sterile and Tissue Culture Treated (50 plates)	6005680
OptiPlate-384, White Opaque 384-well Microplate (50 plates)	6007290
CulturPlate-384, White Opaque 384-well Microplate, Sterile and Tissue Culture Treated (50 plates)	6007680
ProxiPlate-384 Plus, White 384-shallow well Microplate (50 plates)	6008280
TopSeal-A PLUS (1 box)	6050185

Other packages available



## ASSAY FORMATS



### 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) to find a list of our regional distributors

