

# CISBIO STARTER PACKS

CISBIO STARTER PACK



## FAST, EASY, AND ACCURATE RESEARCH

Are you ready to discover a world with shorter time to result, easier experiments and quantitative data every time?

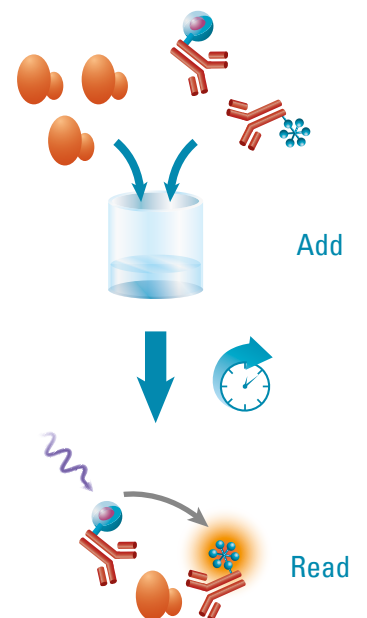
Welcome to HTRF!

Explore cell function inside out, discover our starter packs including everything you need to set-up your HTRF assay now.!

### Browse our starter packs

- 96 well kits including plate
- 500 well kits for up scale
- Plates available separately
- Step by step videos and data analysis

Sample  
16  $\mu$ L      Detection  
reagent 4  $\mu$ L



## PHOSPHO ASSAYS

### EXPLORE THE MAJOR PATHWAYS

- < 2 hours at bench
- Quantitative
- Specific
- Sample variety (cell culture, tissues homogenates ...)
- Normalization with Total assay

## CYTOKINE ASSAYS

### QUANTIFY EASILY CELL CYTOKINE RELEASE

- No wash protocol
- < 2 hours at bench
- Quantitative
- 4PL regression

## CISBIO LOW VOLUME 96 WELL PLATES

- Small sample size (total assay 20 µL)
- Compatible with all Cisbio catalogue
- HTRF validated (intra & inter batch)
- Compatible with all HTRF readers

## EXPERT SERVICES TO SATISFY YOUR EXACT NEEDS



In addition to providing ready-to-use kits with all the necessary reagents, Cisbio's team of experienced experts works hand-in-hand with its clients to create custom solutions. When you have specific requirements for cellular phospho-proteins, cytokines or other signaling proteins, Cisbio is the ideal trusted partner to develop the assays you need.

Keep your therapeutic research moving forward with faster, more robust, physiologically relevant results. Call us or visit [www.cisbio.com/](http://www.cisbio.com/)

### 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

### ORDERING INFORMATION

PRODUCT	PART#
ACC phospho-S79 kit - 1 x 96 tests	64ACCPET
AKT phospho-S473 kit - 1 x 96 tests	64AKSPET
AKT total kit - 1 x 96 tests	64NKTPET
AMPK phospho-T172 kit - 1 x 96 tests	64MPKPET
BAD phospho-S112 kit - 1 x 96 tests	64BADPET
CREB phospho-S133 kit - 1 x 96 tests	64CREPET
Advanced ERK phospho-T202/Y204 kit - 1 x 96 tests	64AERPET
ERK total kit - 1 x 96 tests	64NRKPET
NFkB phospho-S536 kit - 1 x 96 tests	64NFBPET
P38 phospho-T180/Y182 kit - 1 x 96 tests	64P38PET
P53 phospho-S15 kit - 1 x 96 tests	64P53PET
P53 total kit - 1 x 96 tests	64T53PET
SLP-76 phospho-S376 kit - 1 x 96 tests	63ADK076PET
SMAD1 phospho-S463/465 kit - 1 x 96 tests	63ADK062PET
STAT3 phospho-Y705 kit - 1 x 96 tests	62AT3PET

### ORDERING INFORMATION

PRODUCT	PART#
Human IFN gamma - 1 x 96 tests	62HIFNGPET
Human IL1 beta - 1 x 96 tests	62HIL1BPET
Human IL6 - 1 x 96 tests	62HIL06PET
Human IL8 - 1 x 96 tests	62HIL06PET
Human IL17 - 1 x 96 tests	62HIL178PET
Human TNF alpha - 1 x 96 tests	62HTNFAPET

### ORDERING INFORMATION

PRODUCT	PART#
HTRF 96 well low volume plate	66PL96001
HTRF 96 well low volume plate - 5 plates	66PL96005
HTRF 96 well low volume plate - 25 plates	66PL96025
HTRF 96 well low volume plate - 100 plates	66PL96100