

CISBIO STARTER PACKS



FAST, EASY, AND ACCURATE RESEARCH

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

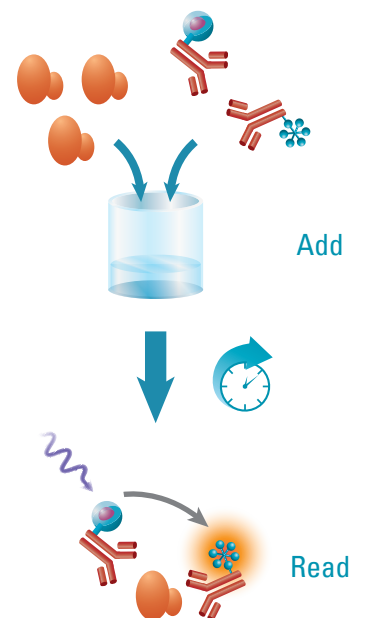
Welcome to HTRF!

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 μ L Detection
reagent 4 μ 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/

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 to find a list of our regional distributors

ORDERING INFORMATION

PRODUCT	PART#	PRICE
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	425 €
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#	PRICE
Human IFN gamma - 1 x 96 tests	62HIFNGPET	
Human IL1 beta - 1 x 96 tests	62HIL1BPET	
Human IL6 - 1 x 96 tests	62HIL06PET	383 €
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#	PRICE
HTRF 96 well low volume plate	66PL96001	12€
HTRF 96 well low volume plate - 5 plates	66PL96005	55 €
HTRF 96 well low volume plate - 25 plates	66PL96025	225 €
HTRF 96 well low volume plate - 100 plates	66PL96100	700 €