

HTRF SOLUTIONS FOR AUTOIMMUNITY



AUTOIMMUNE DISEASE

Built on the proven reliability and efficiency of HTRF, Cisbio has developed an ample portfolio of products to assist with your autoimmune disease research. Whether your focus is lupus, multiple sclerosis, psoriasis, inflammatory bowel disease, rheumatoid arthritis, or autoimmune arthritis, Cisbio is able to provide you with a wide range of solutions for your drug discovery.

Receive faster, accurate results with HTRF

- Easy wash free mix-and-read protocol
- Minimized false negatives and repeat runs
- Consistent, adaptable, reliable data
- Truly homogeneous time-saving technology
- Broad dynamic range with high sensitivity

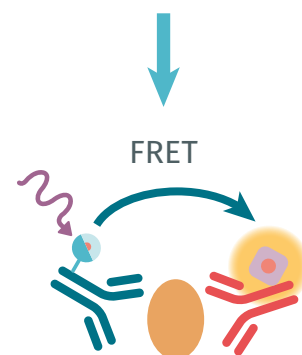
Products covering all major inflammation pathways

- STING -cGAS pathway
- NF-kB pathway
- JAK-STAT pathway
- BTK pathway
- ZAP-70 pathway

 Mab anti protein-d2

 Mab anti protein-Eu cryptate

 Target protein



**cisbio**
INTERACTION IS EVERYTHING

HTRF SOLUTIONS FOR LUPUS, MULTIPLE SCLEROSIS, INFLAMMATORY BOWEL DISEASE, RHEUMATOID ARTHRITIS, AUTOIMMUNE ARTHRITIS, AND PSORIASIS

CYTOKINES

Product	Tests	Cat. No.#
IFN beta kit	500	62HIFNBPEG
IFN alpha kit	500	62HIFNAPEG
Human IFN gamma kit	500	62HIFNGPEG
Mouse IFN gamma kit	500	62MIFNGPEG
Human TNF alpha kit	500	62HTNFAPEG
Cynomolgus TNF alpha kit	500	62CTNFAPEG
Mouse TNF alpha kit	500	62MTNFAPEG
Human IL-2 kit	500	62HIL02PEG
Human IL-4 kit	500	62HIL04PEG
Human IL-5 kit	500	62HIL05PEG
Human IL-6 kit	500	62HIL06PEG
Human IL-10 kit	500	62HIL10PEG
Human IL-12 kit	500	62HIL12PEG
Human IL-17 kit	500	62HIL17PEG
Mouse IL-2 kit	500	62MIL02PEG
Mouse IL-4 kit	500	62MIL04PEG
Mouse IL-5 kit	500	62MIL05PEG
Mouse IL-6 kit	500	62MIL06PEG
Mouse IL-10 kit	500	62MIL10PEG
Mouse IL-12 kit	500	62MIL12PEG
Mouse IL-17 kit	500	62MIL17PEG
Human CCL2 kit	500	62HCCL2PEG
Mouse CCL2 kit	500	62MCCL2PEG
Human CCL3 kit	500	62HCCL3PEG
Human CCL4 kit	500	62HCCL4PEG
Human CCL5 kit	500	62HCCL5PEG
Human CCL20 kit	500	62HCCL20PEG
Human CXCL1 kit	500	62HCXC1PEG
Human CXCL10 kit	500	62HCXC10PEG

KINASE (BIOCHEMICAL)

Product	Tests	Cat. No.#
KinEASE STK discovery kit	1,000	62ST0PEB
KinEASE-TK kit	1,000	62TK0PEB

PPI

Product	Tests	Cat. No.#
PD1/PDL1 binding assay kit	500	64ICP01PEG
CTLA4/B7-1 binding assay kit	500	64ICP04PEG
CTLA4/B7-2	500	64ICP05PEG
LAG3/MHCI binding assay kit	500	64ICP03PEG
CD47/SIRPa binding assay kit	500	62ICP02PEG
Human STING WT binding kit	500	64BDSTGPEG
Human H232 STING binding kit NEW	500	64BDSTGHPEG
Human AQ STING binding kit NEW	500	64BDSTGQPEG

PHOSPHO-PROTEINS

Product	Tests	Cat. No.#
Alpha tubulin kit	500	64ATUBPEG
AKT1 phospho-S473	500	63ADK078PEG
AKT1 total kit	500	63ADK079PEG
AKT2 phospho-S473	500	63ADK080PEG
AKT2 total kit	500	63ADK081PEG
AKT3 phospho-S473	500	63ADK082PEG
AKT3 total kit	500	63ADK083PEG
BTK phospho-Y223	500	63ADK017PEG
BTK total kit	500	63ADK064PEG
BAD phospho-S112	500	64BADPEG
CREB phospho-S133	500	64CREPEG
CREB total kit	500	63ADK052PEG
Advanced ERK phospho-T202/Y204	500	64AERPEG
ERK total kit	500	64NRKPEG
GSK3 alpha phospho-S21 kit	500	64GPAPEG
GSK3 beta phospho-S9 kit	500	64GPBPEG
GSK3 alpha total kit	500	64GTAPEG
GSK3 beta total kit	500	64GTBPEG
Human STING binding kit	500	64BDSTGPEG
IRF3 kit	500	6FRF3PEG
IKK β phospho-S177/181 kit	500	64KKBPEG
IKK β total kit	500	63ADK09PEG
JNK phospho-S2448 kit	500	64JNKPEG
mTOR phospho-S2448 kit	500	64TORPEG
NF-kB phospho-S172 kit	500	64NFBPEG
NF-kB total kit	500	64NFTPEG
p38 phospho-T180/Y182 kit	500	64P38PEG
SLP-76 phospho-S376 kit	500	63ADK076PEG
SLP-76 total kit	500	63ADK077PEG
STAT1 phospho-Y701 kit	500	63ADK026PEG
STAT1 total kit	500	63ADK096PEG
STAT3 phospho-Y705	500	62AT3PEG
STAT3 total kit	500	62AT3PEG
STAT5 phospho-Y694 kit	500	64AT5PEG
STING phospho-S366 kit	500	64STGPEG
STING total kit	500	64NTGEG
TBK1 total kit	500	62NTBPEG
TBK1 phospho-S172 kit	500	64TBKPEG
ZAP-70 phospho-Y319 kit	500	64ZAPPEG
ZAP-70 total kit	500	64ZATPEG

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

