

HTRF Solutions for Cardiovascular Disease



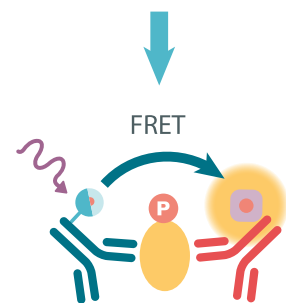
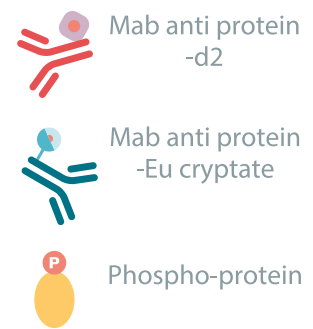
CARDIOVASCULAR

PHYSIOLOGICALLY RELEVANT RESULTS FOR FAST FLOWING RESEARCH

Built on the proven convenience and reliability of HTRF, Cisbio's scientists have developed an ample portfolio of kits and reagents catering to researchers studying hypertension, nitric oxide donors, atherosclerosis, cardiac hypertrophy and arrhythmia. cGMP, cAMP, IP-1, PCSK9, Phospho-Phospholamban, Apolipoproteins, cytokines are only the start of the Cisbio's all homogenous solutions for cardiovascular research.

It's your turn to benefit from HTRF

- Easy wash free mix-and-read protocol
- Minimized false negatives and repeat runs
- Consistent, adaptable, reliable data



HTRF sandwich assay principle

WHAT'S YOUR FOCUS? WHATEVER YOUR TARGET, HTRF IS YOUR SOLUTION

Our kits are available in larger formats (5,000 - 10,000 - 20,000 - 50,000 or 100,000). Check our website for details and availability.

<https://www.cisbio.com/drug-discovery/assay-kits-reagents>

CARDIOPROTECTION & CARDIOMIOCYTE GROWTH

Product	Tests	Cat. No.#
AKT Total kit	1x96	64NKTPET
AKT Total kit	500	64NKTPEG
AKT Phospho - S473 kit	1x96	62AKSPET
AKT Phospho - S473 kit	500	64AKSPEG
AKT Phospho - S308 kit	500	64AKTPEG
AKT1 Phospho - S473 kit	500	63ADK078PEG
AKT1 Total kit	500	63ADK079PEG
AKT2 Phospho S473	500	63ADK080PEG
AKT2 Total kit	500	63ADK081PEG
AKT3 Phospho S473 kit	500	63ADK082PEG
AKT3 Total kit	500	63ADK083PEG
mTOR Phospho-S2448 kit	500	64TORPEG
STAT1 Phospho-Y701 kit	500	63ADK026PEG
STAT1 Total kit	500	63ADK096PEG
STAT3 Phospho-Y705 kit	1x96	62AT3PET
STAT3 total kit	500	62NT3PEG
4EBP1 Phospho-T37/46 kit	500	64BP1PEG

CONTRACTION/RELAXATION & BLOOD PRESSURE

Product	Tests	Cat. No.#
cAMP - Gs HiRange kit	1,000	62AMPEB
cAMP - Gi Dynamic kit	1,000	62AM9PEB
cAMP - Gs Dynamic kit	1,000	62AM4PEB
IBMX - Phosphodiesterase Inhibitor	40 µL	62AMXADA
IBMX - Phosphodiesterase Inhibitor	500 µL	62AMXADC
Forskolin - Adenylyl Cyclase Activator	40 µL	62AMYADA
Forskolin - Adenylyl Cyclase Activator	800 µL	62AMYADC
IP-One - Gq kit	1,000	62IPAPEB

HEART FIBROSIS & CARDIOMIOCYTE APOPTOSIS

Product	Tests	Cat. No.#
SMAD1 Phospho-SS463/465	1x96	63ADK062PET
SMAD1 Phospho-S463/465	500	63ADK062PEG
SMAD1 Total kit	500	63ADK063PEG
SMAD3 Phospho-S423/425	500	63ADK025PEG
SMAD3 total kit	500	64ND3PEG
TGF Beta kit	500	62HTGFBPEG

TRANSPORT OF CHOLESTEROL & ATHEROSCLEROSIS

Product	Tests	Cat. No.#
Human Apolipoprotein A1 kit	500	64APAPEG
Human Apolipoprotein B kit	500	64APBPEG
Human Apolipoprotein C3 kit	500	63ADK001PEG
Human Apolipoprotein E kit	500	63ADK004PEG
Human PCSK-9 kit	500	63ADK050PEG

NO PATHWAY VASORELAXATION

Product	Tests	Cat. No.#
cGMP kit	500	62GM2PEG
VASP Phospho-S239 kit	500	63ADK065PEG
VASP Phospho-S157 kit	500	63ADK066PEG
VASP Total kit	500	63ADK067PEG

EXPERT SERVICES TO SATISFY YOUR EXACT NEEDS



In addition to providing ready-to-use solutions, Cisbio's team of experienced experts also works hand-in-hand with clients to create custom solutions.

From compound characterization, to small- and large-scale reagent preparation and full assay development, Cisbio is the ideal trusted partner to support all your needs.

Keep your therapeutic research moving forward with faster, more robust, physiologically relevant results. Call us or visit

<https://www.cisbio.com/drug-discovery/assay-kits-reagents>

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

