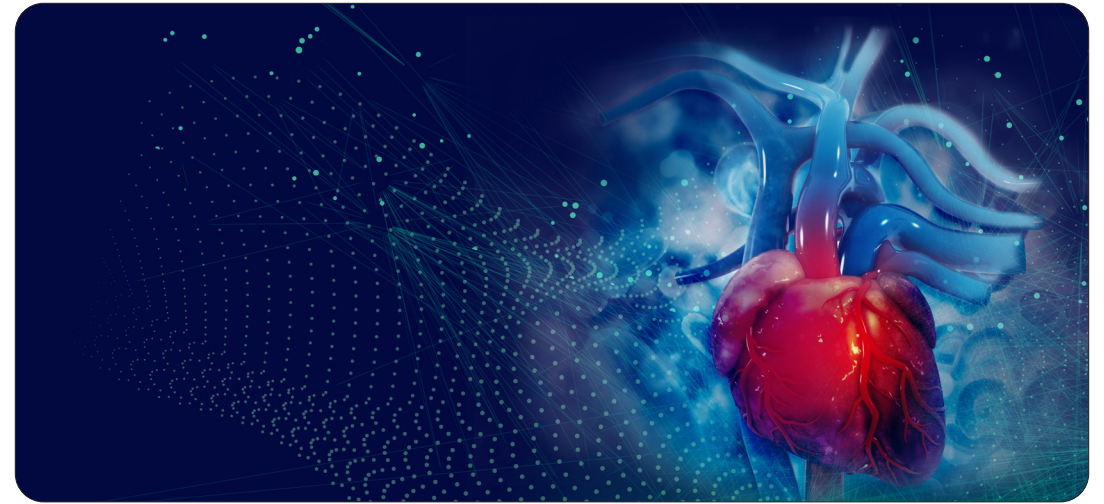
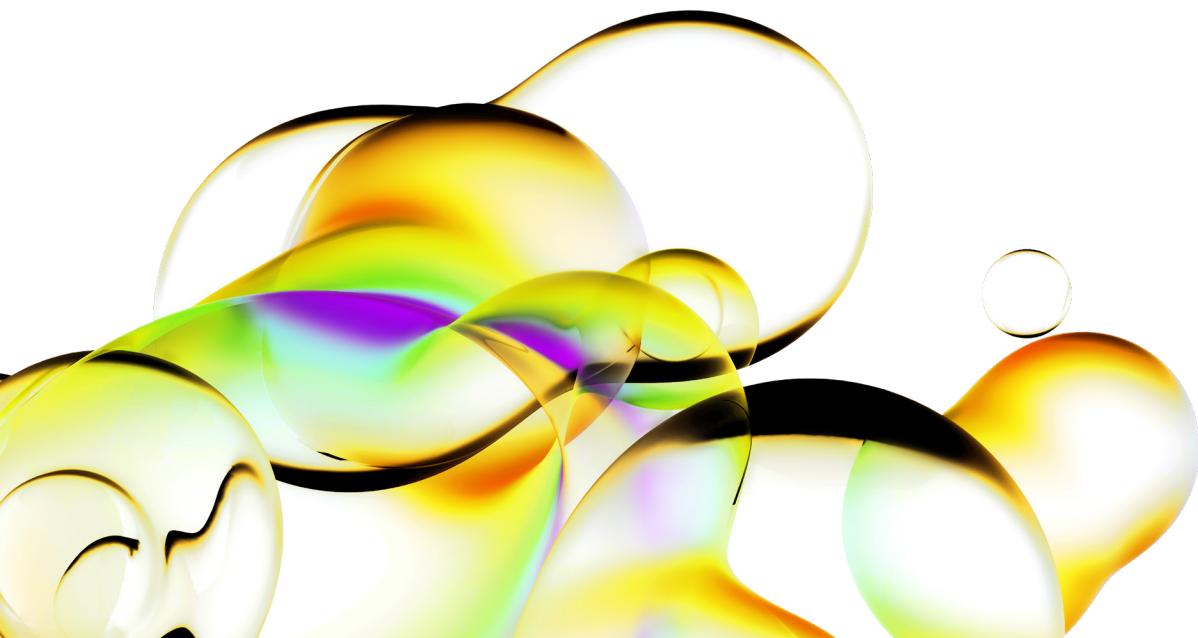




Solutions for cardiovascular disease.



Physiologically relevant results for fast flowing research

Human cardiovascular diseases are a significant health problem worldwide due to the large number of people affected and the incomplete knowledge of the cardiovascular process's pathogenesis. Cardiovascular disorders can develop in several different diseases such as hypertension, hypercholesterol, atherosclerosis, heart fibrosis or strokes; all of which have ties with the regulation of blood pressure, fat transportation, artery diameter and the state of inflammation in arteries and vessels.

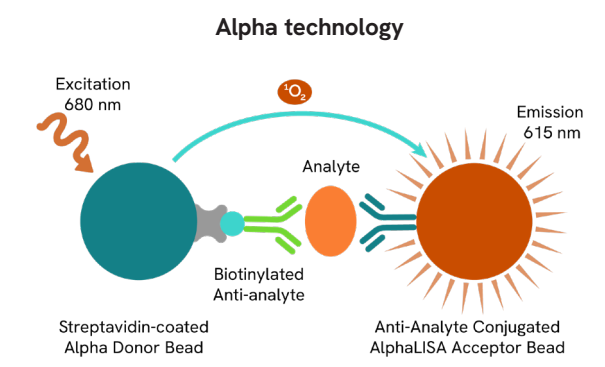
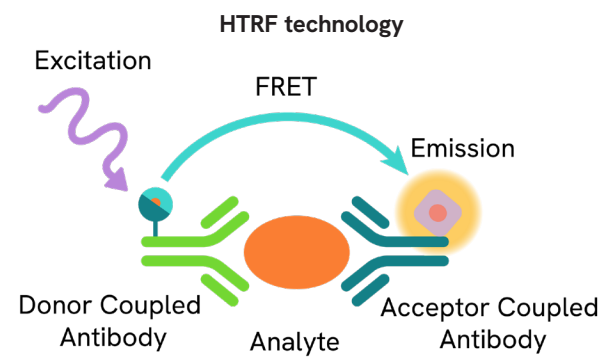
Alpha is a versatile, bead-based platform that enables you to assay the most complex samples in one well and with no wash steps. HTRF is a fast, sensitive, homogeneous, and ready to use assay platform with no wash steps.

Advance your research on Cardiovascular disorders and benefit from Revvity's broad offering of reagent technologies:

- Products covering all significant contributors to a cardiovascular disorder involved in cardiomyocyte health, vasorelaxation, artery inflammation and heart fibrosis.
- Easy-to-use Alpha, HTRF and LANCE Ultra assay technologies enable the cell-based detection of cytokines, biomarkers, and cell signaling proteins, providing fast and accurate results."

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.



Cardioprotection and cardiomyocyte growth

Product	Tests	HTRF Part#	Alpha Part#
AKT1, phospho-S473	500	63ADK078PEG	ALSU-PAKT1-C500
AKT1, Total	500	63ADK079PEG	ALSU-TAKT1-B500
AKT1/2/3, phospho-S473	500		ALSU-PAKT-B500
AKT1/2/3, phospho-T308	500		ALSU-PAKT-A500
AKT1/2/3, phospho-T450	500		ALSU-PAKT-C500
AKT1/2/3, Total	500		ALSU-TAKT-B500
mTOR, phospho-S2448	500	64TORPEG	ALSU-PMTOR-C500
mTOR, phospho-S2481	500		ALSU-PMTOR-B500
mTOR, Total	500		ALSU-TMTOR-B500
STAT1, phospho-S727	500		ALSU-PST1-B500
STAT1, phospho-Y701	500		ALSU-PST1-A500
STAT1, Total	500	63ADK096PEG	ALSU-TST1-A500
STAT3, phospho-Y705	500		ALSU-PST3-A500
STAT3, Total	500	64NT3PEG	ALSU-TST3-A500
STAT4, phospho-Y693	500		ALSU-PST4-A500
STAT4, Total (human)	500		ALSU-TST4-A500
STAT5, phospho-Y694/699	500		ALSU-PST5-B500
STAT5, Total	500		ALSU-TST5-A500
STAT6, phospho-Y641	500	64AT6PEG	ALSU-PST6-A500
STAT6, Total (human)	500		ALSU-TST6-A500
AKT, phospho-S473	500	64AKSPEG	
AKT, phospho-T308	500	64AKTPEG	MPSU-PAKT-E500

Product	Tests	HTRF Part#	Alpha Part#
AKT, Total	500	64NKTPEG	
AKT2, phospho-S473	500	63ADK080PEG	
AKT2, Total	500	63ADK081PEG	
AKT3, phospho-S473	500	63ADK082PEG	
AKT3, Total	500	63ADK083PEG	
STAT1, phospho-Y701	500	63ADK026PEG	
STAT3, phospho-Y705	500	62AT3PEG	
STAT5, phospho-Y694	500	64AT5PEG	
STAT6, Total (human; mouse)	500	64STAT6TPEG	
AKT1/2/3, phospho-S473 & AKT1, Total	500		MPSU-PTAKT-M500
AKT1/2/3, phospho-S473 & ERK 1/2, phospho-T202/Y205	500		MPSU-PAKER-K500

Contraction/relaxation and blood pressure

Product	Tests	HTRF Part#	Alpha Part#
cAMP - Gs HiRange	1,000	62AMPEB	
cAMP - Gi Dynamic	1,000	62AM9PEB	
cAMP - Gs Dynamic	1,000	62AM4PEB	
IBMX - Phosphodiesterase Inhibitor	40 µL	62AMXADA	
Forskolin - Adenylyl Cyclase Activator	40 µL	62AMYADA	
IP-One - Gq	1,000	62IPAPEB	AL3145D

Heart fibrosis and cardiomyocyte apoptosis

Product	Tests	HTRF Part#	Alpha Part#
TGF-b1	500	62HTGFBPEG	AL336C
TGF-b1, Biotin Free (human)	500		AL361C
SMAD1, phospho-S463/465	500	63ADK062PEG	ALSU-PSM1-A500
SMAD1, Total	500	63ADK063PEG	ALSU-TSM1-A500
SMAD2, phospho-S465/467	500	64SMAD2S5PEG	ALSU-PSM2-A500
SMAD2, Total	500	64SMAD2TPEG	ALSU-TSM2-A500
SMAD3, phospho-S423/425	500	63ADK025PEG	ALSU-PSM3-A500

Product	Tests	HTRF Part#	Alpha Part#
SMAD3, Total	500	64ND3PEG	ALSU-TSM3-A500
SMAD4, phospho-T277		64SMAD4T7PEG	
SMAD4, Total	500	64SMAD4TPEG	ALSU-TSM4-A500

Transport of cholesterol and atherosclerosis

Product	Tests	HTRF Part#	Alpha Part#
Apolipoprotein A1 (human)	500	64APAPEG	AL389C
Apolipoprotein B (human)	500	64APBPEG	AL392C
Apolipoprotein C3 (human)	500	63ADK001PEG	
Apolipoprotein E (human)	500	64APOEPEG	AL395C
PCSK9 (human)	500	63ADK050PEG	AL270C

Vasorelaxation

Product	Tests	HTRF Part#
cGMP	500	62GM2PEG
eNOS, phospho-S1177	500	64ENOSS7PEG
eNOS, Total	500	64ENOSTPEG
Phospholamban, phospho-S16 (human)	500	64PLN16PEG
Phospholamban, phospho-T17 (human)	500	63ADK074PEG
Phospholamban, Total (human)	500	63ADK075PEG
VASP, phospho-S157	500	63ADK066PEG
VASP, phospho-S239	500	63ADK065PEG
VASP, Total	500	63ADK067PEG

Expert services to satisfy your exact needs

- In addition to providing ready-to-use solutions, Revvity’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, Revvity 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 [Revvity.com](https://www.revvity.com).

