



HTRF

SpectraMax Detection Cartridge

HTRF® SpectraMax® Detection Cartridges for SpectraMax® i3 and Paradigm® Systems utilize a high-energy Xenon flash lamp for sensitive reading of HTRF reagents. This dual emission design on the SpectraMax Paradigm System gives the most accurate results while accelerating throughput 2X.

HTRF SpectraMax Detection Cartridge

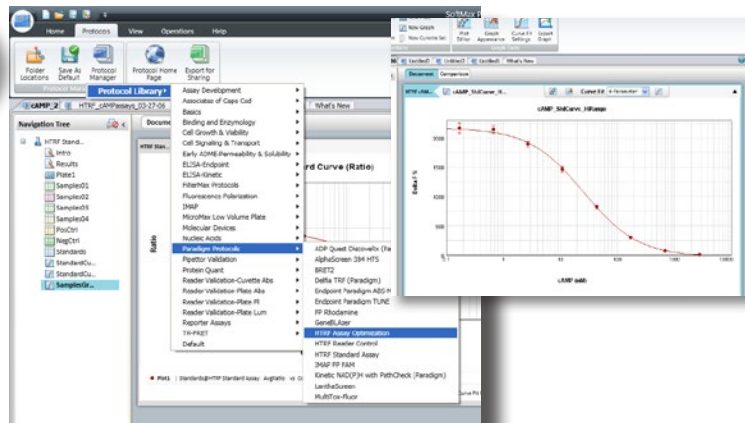
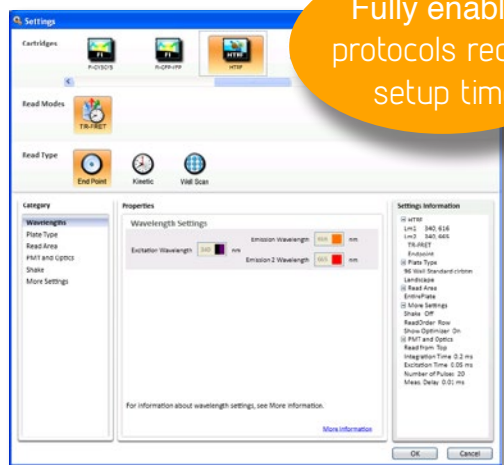
Description

Compatible systems	SpectraMax i3 and Paradigm Readers
Part number	0200-7011
Mounting	Top read
Cartridge size	2-Position
Light source	Xenon flash lamp
Excitation range	EX: HTRF
Emission range	EM1 616/10 nm EM2 665/10 nm

Certified for use with Cisbio's HTRF Europium Cryptate reagents.



Fully enabled protocols reduce setup time



SoftMax® Pro Software protocols not only offer preconfigured microplate reader settings but also contain predefined data reductions and calculations to instantly provide the results you seek.

Typical applications

The HTRF cartridge provides optimal performance for key applications for drug discovery, bioprocess development, and others.

- GPCR Assays
- Cytokine Assays
- cAMP Quantitation
- Human Ab Screening
- Kinase Assays

For more information and typical applications refer to <http://www.htrf.com>

HTRF is a registered trademark of CisBio. The fluorescence ratio associated with the HTRF readout is a correction method developed by CisBio and covered by the US patent 5,527,684 and its foreign equivalents, for which CisBio has granted a license. Its application is strictly limited to the use of HTRF reagents and technology, excluding any other TR-FRET technologies.

Contact Us

Phone: +1-800-635-5577
Web: www.moleculardevices.com
Email: info@moldev.com
Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1-800-635-5577
Brazil +55-11-3616-6607
China (Beijing) +86-10-6410-8669
China (Shanghai) +86-21-3372-1088
Germany 00800-665-32860

Japan (Osaka) +81-6-7174-8831
Japan (Tokyo) +81-3-6362-5260
South Korea +82-2-3471-9531
United Kingdom +44-118-944-8000

