

HTRF[®] Europium cryptate donor / Red acceptor readout Setup recommendations for SpectraMax[®] Paradigm[®]

To read HTRF[®], the SpectraMax[®] Paradigm[®] must be first equipped with the SpectraMax[®] Paradigm[®] Cisbio HTRF[®] cartridge, which enables the simultaneous measurement of both 620 nm donor and 665 nm acceptor emissions. The ratio of the two fluorescence intensities 665/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

HTRF[®] readout can be achieved by SpectraMax[®] Paradigm[®] readers after the installation of the HTRF[®] dedicated cartridge, which includes the optimized excitation and emission filters, the light source and the dichroic mirrors. The measurement conditions should then be set up in the SoftMax[®] Pro software according to the following indications:

Setup

Cartridge	HTRF Detection cartridge 0200-7011
Number of flashes	30
Integration delay (lag time)	30 μ s
Integration time	500 μ s
Optimal z-position	Volume and plate format dependant. Must be optimized before each new configured measurement using the labware optimization procedure of the software.



HTRF[®] Terbium cryptate donor / Green acceptor readout Setup recommendations for SpectraMax[®] Paradigm[®]

To read HTRF[®], the SpectraMax[®] Paradigm[®] must be first equipped with the SpectraMax[®] Paradigm[®] Cisbio HTRF[®] cartridge, which enables the simultaneous measurement of both 620 nm donor and 520 nm acceptor emissions. The ratio of the two fluorescence intensities 520/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

HTRF[®] readout can be achieved by SpectraMax[®] Paradigm[®] readers after the installation of the HTRF[®] dedicated cartridge, which includes the optimized excitation and emission filters, the light source and the dichroic mirrors. The measurement conditions should then be set up in the SoftMax[®] Pro software according to the following indications:

Setup

Cartridge	TR-FRET SpectraMax Paradigm Ex 340/80 – EM1 520/15- EM2 620/10
Number of flashes	30
Integration delay (lag time)	30 μ s
Integration time	500 μ s
Optimal z-position	Volume and plate format dependant, Must be optimized before each new configured measurement using the labware optimization procedure of the software Volume and plate format dependant



HTRF® Terbium cryptate donor / Red acceptor readout Setup recommendations for SpectraMax® Paradigm®

To read HTRF®, the SpectraMax® Paradigm® must be first equipped with the SpectraMax® Paradigm® Cisbio HTRF® cartridge, which enables the simultaneous measurement of both 620 nm donor and 665 nm acceptor emissions. The ratio of the two fluorescence intensities 665/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

HTRF® readout can be achieved by SpectraMax® Paradigm® readers after the installation of the HTRF® dedicated cartridge, which includes the optimized excitation and emission filters, the light source and the dichroic mirrors. The measurement conditions should then be set up in the SoftMax® Pro software according to the following indications:

Setup

Cartridge	HTRF Detection Cartridge 0200-7011
Number of flashes	30
Integration delay (lag time)	30 µs
Integration time	500 µs
Optimal z-position	Volume and plate format dependant. Must be optimized before each new configured measurement using the labware optimization procedure of the software Volume and plate format dependant,

