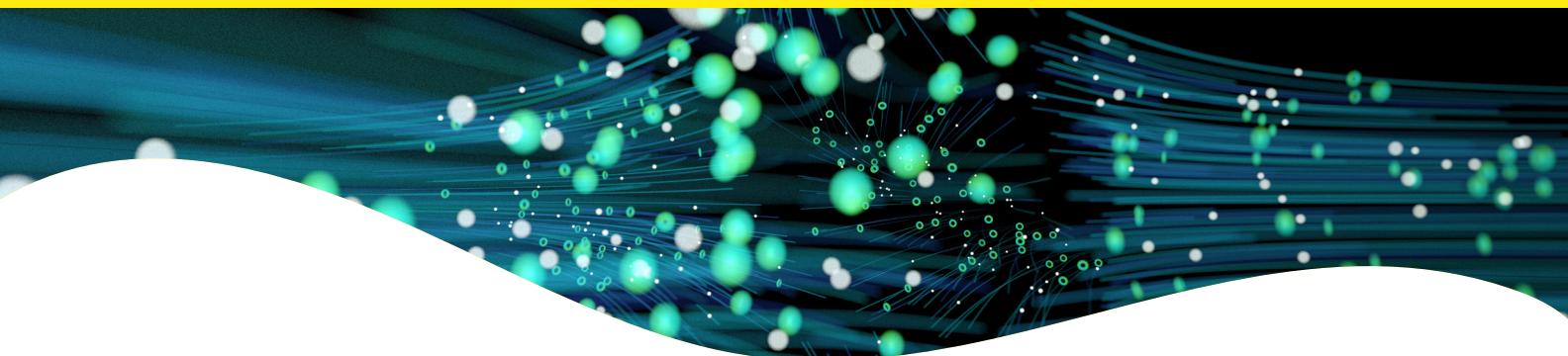


HTRF setup recommendations for SpectraMax Paradigm.



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 Revvity 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.

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 Revvity 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 Revvity 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 Volume and plate format dependant,

