

HTRF[®] Europium cryptate donor / Red acceptor readout Setup recommendations for SpectraMax i3[®]

To read HTRF[®], the SpectraMax i3[®] must be first equipped with the SpectraMax i3[®] 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 i3[®] 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
Number of flashes	30
Integration delay (lag time)	70 μ s
Integration time	400 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.



**The fluorescence ratio is a correction method developed by Cisbio Bioassays with an application limited to the use of HTRF[®] reagents and technology, and for which Cisbio Bioassays has granted a licence to Molecular Devices. The method is covered by the US patent 5,527,684 and its foreign equivalents.*

HTRF® Terbium cryptate donor / Green acceptor readout Setup recommendations for SpectraMax i3®

To read HTRF®, the SpectraMax i3® must be first equipped with the SpectraMax i3® 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 i3® 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
Number of flashes	30
Integration delay (lag time)	70 µs
Integration time	400 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



**The fluorescence ratio is a correction method developed by Cisbio Bioassays with an application limited to the use of HTRF® reagents and technology, and for which Cisbio Bioassays has granted a licence to Molecular Devices. The method is covered by the US patent 5,527,684 and its foreign equivalents.*

HTRF® Terbium cryptate donor / Red acceptor readout Setup recommendations for SpectraMax i3®

To read HTRF®, the SpectraMax i3® must be first equipped with the SpectraMax i3® 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 i3® 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
Number of flashes	30
Integration delay (lag time)	70 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,



**The fluorescence ratio is a correction method developed by Cisbio Bioassays with an application limited to the use of HTRF® reagents and technology, and for which Cisbio Bioassays has granted a licence to Molecular Devices. The method is covered by the US patent 5,527,684 and its foreign equivalents.*