

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,



**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.*