

## HTRF<sup>®</sup> Terbium cryptate donor / Green acceptor readout Setup recommendations for SpectraMax<sup>®</sup> Paradigm<sup>®</sup>

To read HTRF<sup>®</sup>, the SpectraMax<sup>®</sup> Paradigm<sup>®</sup> must be first equipped with the SpectraMax<sup>®</sup> Paradigm<sup>®</sup> Cisbio HTRF<sup>®</sup> 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<sup>®</sup> readout can be achieved by SpectraMax<sup>®</sup> Paradigm<sup>®</sup> readers after the installation of the HTRF<sup>®</sup> 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<sup>®</sup> 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 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<sup>®</sup> 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.*