

HTRF[®] readout - Set up recommendations for PARADIGM[™] readers

To read HTRF[®], the PARADIGM[™] must be equipped with the PARADIGM[™] Cisbio HTRF[®] cartridge, which enables the simultaneous measurement of both the 620nm donor emission and the 665nm acceptor. A ratio of the two fluorescence intensities* (acceptor/donor) then allows the calculation of Delta F (%), i.e. the relative energy transfer rate for each data point.

HTRF[®] readout can be achieved by PARADIGM[™] readers after the installation of the HTRF[®] dedicated cartridge, which includes the optimised excitation and emission filters, the light source and the dichroic mirrors. The measurement conditions should then be set up in the DTX Multimode Detection software according to the following indications:

Setup

Cartridge	: PARADIGM [™] Cisbio HTRF [®]
Number of flashes	: 30
Delay (Lag time)	: 30 μ s
Integration time	: 500 μ s
Optimal z-position	: volume and plate format dependent. 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 CIS bio international with an application limited to the use of HTRF[®] reagents and technology, and for which CIS bio international has granted a licence to Beckman Coulter. The method is covered by the US patent 5,527,684 and its foreign equivalents.*