

HTRF® Terbium cryptate donor / Green acceptor readout Setup recommendations for PHERAstar FSX Laser

PHERAstar FSX is equipped with a specific optical device, which enables the simultaneous measurement of both 620 nm cryptate and 520 nm acceptor emissions. The ratio* of the fluorescence intensities 520/620 (acceptor/donor) allows the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

HTRF® readout can be achieved by PHERAstar FSX after the installation of the HTRF® dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup

| | |
|------------------------------|-----------------------------------|
| Optic module | HTRF® 337/620/520 Ref.: 1010B1 |
| Energy source | Laser |
| Integration delay (lag time) | 60 µs |
| Integration time | 400 µs |
| Number of flashes | 60 |
| Optimal z-pos § | Volume and plate format dependent |

§The focal height "z" is automatically calculated according to the plate format and the final working volume dispensed in the plate.



**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 BMG LABTECH.. The method is covered by the US patent 5,527,684 and its foreign equivalents.*