

## HTRF® Terbium cryptate donor / Green acceptor readout Setup recommendations for Envision equipped in Laser

Envision reader must be 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 two fluorescence intensities 665/520 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

Envision reader must be appropriately configured for HTRF® readout by setting up the measurement conditions in the software according to the following indications:

### Setup

---

Top mirror	TRF LASER D407/D505 advanced dual mirror (#665) or TRF LASER D400/D505 dual mirror (#447)  TRF LASER advanced single mirror (#616) or TRF LANCE/DELFI A D400 single mirror (#445)
Emission filter	TRF 520 (#275)
2 <sup>nd</sup> emission filter	CY5 620 (#118) or 615 (#203)
Measurement height	12 for 384sv – 7 for 96LV otherwise must be determined
Cycle	16600µs
Delay	50µs
Number of flashes	20
Number of flashes for 2 <sup>nd</sup> detector	20
Total time of windows	400µs

