

HTRF® Europium cryptate donor / Red 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 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.

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/D630 advanced dual mirror (#667) or TRF LASER D400/D630 dual mirror (#446) TRF LASER advanced single mirror (#616) or TRF LANCE/DELFI A D400 single mirror (#445)
Emission filter	APC 665 (#205)
2 nd 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	50 for white plate / 100 for black plate
Number of flashes for 2 nd detector	50 for white plate / 100 for black plate
Total time of windows	400µs

