

HTRF® Europium cryptate donor / Red acceptor readout Setup recommendations for Envision equipped in Flash lamp

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 Lance D407/D630 advanced dual mirror (#666) or TRF LANCE DELFIA D400/630 dual mirror (#662) TRF advanced single mirror (#615) or TRF LANCE/DELFIA D400 single mirror (#412) |
| Excitation filter | UV2 (TRF) 320 (#111) |
| 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 | 2000µs |
| Delay | 60µs |
| Number of flashes | 100 |
| Number of flashes for 2 nd detector | 100 |
| Total time of windows | 400µs |

This reader only allows high performance HTRF measurement when assays are run in WHITE plates.

