

HTRF® Terbium cryptate donor / Green 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 520nm acceptor emissions. The ratio* of the two fluorescence intensities 520/620 (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 D407/D505 advanced dual mirror (#664) or TRF D400/D505 dual mirror (#657)
	TRF advanced single mirror (#615) or TRF LANCE/DELFI A D400 single mirror (#412)
Excitation filter	UV2 (TRF) 340 (#101)
Emission filter	TRF 520 (#275)
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.

