

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 (mirror and filters), which enables the 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/DELFI A D400/D630 dual mirror (#662) TRF advanced single mirror (#615) or TRF LANCE/DELFI A 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 via Optimization
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.



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 (mirror and filters), which enables the measurement of both 620 or 495 nm cryptate and 520nm acceptor emissions. The ratio* of the two fluorescence intensities 520/620 or 520/495 (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/DELFLIA D400 single mirror (#412)
Excitation filter	UV2 (TRF) 340 (#101)
Emission filter	Trf 520 (#275) or (#226)
2 nd emission filter	495 (#276) for simultaneous readouts (dual mirror) Note: CY5 620 (#118) or 615 (#203) can be used with dual mirror but 2 successive readouts should be done (520nm then 620nm or 615nm)
Measurement height	12 for 384sv – 7 for 96LV otherwise must be determined via Optimization
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.



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

Envision reader must be equipped with a specific optical device (mirror and filters), which enables the 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) 340 (Barcode #101)
Emission filter	APC 665 (Barcode #205)
2 nd emission filter	CY5 620 (Barcode #118 or #203)
Measurement height	12 for 384sv – 7 for 96LV otherwise must be determined via Optimization
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.



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

Envision reader must be equipped with a specific optical device (mirror and filters), which enables the 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 via Optimization
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



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

Envision reader must be equipped with a specific optical device (mirror and filters), which enables the measurement of both 620 nm or 495nm cryptate and 520 nm acceptor emissions. The ratio* of the two fluorescence intensities 520/620 or 520 /495 (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/DELFLIA D400 single mirror (#445)
Emission filter	Trf 520 (#275) or (#226)
2 nd emission filter	495 (#276) for simultaneous readouts (dual mirror) Note: CY5 620 (#118) or 615 (#203) can be used with dual mirror but 2 successive readouts should be done (520nm then 620nm or 615nm)
Measurement height	12 for 384sv – 7 for 96LV otherwise must be determined via Optimization
Cycle	16600µs
Delay	50µs
Number of flashes	20
Number of flashes for 2 nd detector	20
Total time of windows	400µs



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

Envision reader must be equipped with a specific optical device (mirror and filters), which enables the 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 via Optimization
Cycle	16600µs
Delay	50µs
Number of flashes	20
Number of flashes for 2 nd detector	20
Total time of windows	200µs

