

HTRF setup recommendations for EnVision



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/DELTA 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) or TRF LASER advanced single mirror (#616) or TRF LANCE/DELFLIA 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/DELFI A 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/DELFLIA 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 |

