

# HTRF Setup recommendations for EnVision Nexus

## EnVision Nexus Lamp

### HTRF® Europium cryptate donor / Red acceptor readout

EnVision® Nexus™ is 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.



HTRF® readout can be achieved by EnVision Nexus after the installation of the HTRF® dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup	
Measurement mode	Standard
Measure from	Top
Emission Channel	Dual - Simultaneous
Filter Module	HTRF®/LANCE (330/665/620) Ref.: HH36794002
Energy source	Lamp
Measurement Height	13 for 384sv otherwise must be optimized
Measurement Time	800 ms
Delay	60 µs
Window Time	400µs
Number of sequential windows	1

## HTRF® Terbium cryptate donor / Red acceptor readout

EnVision Nexus is 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.

HTRF® readout can be achieved by EnVision Nexus after the installation of the HTRF® dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup	
Measurement mode	Standard
Measure from	Top
Emission Channel	Dual - Simultaneous
Filter Module	HTRF®/LANCE (330/665/620) Ref.: HH36794002
Energy source	Lamp
Measurement Height	13 for 384sv otherwise must be optimized
Measurement Time	800 ms
Delay	60 µs
Window Time	200 µs
Number of sequential windows	1

## HTRF® Terbium cryptate donor / Green acceptor readout

EnVision Nexus is equipped with a specific optical device, which enables the simultaneous measurement of both 620 nm cryptate and 520 nm 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.

HTRF® readout can be achieved by EnVision Nexus after the installation of the HTRF® dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup	
Measurement mode	Standard
Measure from	Top
Emission Channel	Dual - Simultaneous
Filter Module	HTRF®/LANCE (330/620/515) Ref.: HH36794004
Energy source	Lamp
Measurement Height	13 for 384sv otherwise must be optimized
Measurement Time	800 ms
Delay	60 µs
Window Time	200 µs
Number of sequential windows	1

## EnVision Nexus Laser

### HTRF® Europium cryptate donor / Red acceptor readout

EnVision Nexus is 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.

HTRF® readout can be achieved by EnVision Nexus after the installation of the HTRF® dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup	
Measurement mode	Standard
Measure from	Top
Emission Channel	Dual - Simultaneous
Filter Module	HTRF®/LANCE (343/665/620) Ref.: HH36794003
Energy source	Laser
Measurement Height	13 for 384sv otherwise must be optimized
Measurement Time	800 ms
Delay	60 µs
Window Time	400µs
Number of sequential windows	1

### HTRF® Terbium cryptate donor / Red acceptor readout

EnVision Nexus is 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.

HTRF® readout can be achieved by EnVision Nexus after the installation of the HTRF® dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup	
Measurement mode	Standard
Measure from	Top
Emission Channel	Dual - Simultaneous
Filter Module	HTRF®/LANCE (343/665/620) Ref.: HH36794003
Energy source	Laser
Measurement Height	13 for 384sv otherwise must be optimized
Measurement Time	800 ms
Delay	60 µs
Window Time	200µs
Number of sequential windows	1

## HTRF® Terbium cryptate donor / Green acceptor readout

EnVision Nexus is equipped with a specific optical device, which enables the simultaneous measurement of both 620 nm cryptate and 520 nm 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.

HTRF® readout can be achieved by EnVision Nexus after the installation of the HTRF® dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup	
Measurement mode	Standard
Measure from	Top
Emission Channel	Dual - Simultaneous
Filter Module	HTRF®/LANCE (343/620/515) Ref.: HH36794005
Energy source	Laser
Measurement Height	13 for 384sv otherwise must be optimized
Measurement Time	800 ms
Delay	60 µs
Window Time	200µs
Number of sequential windows	1

