



pHSense™ setup recommendations for Revvity microplate readers.

pHSense™ readout for EnVision Nexus

EnVision™ Nexus™ is equipped with a specific optical device which enables the measurement of 620nm Europium emission.

pHSense readout can be achieved by EnVision Nexus after the installation of the HTRF/Delfia dedicated optical module which includes the optimizes excitation and emission filters and the dichroic mirror. The measurement conditions should then be set up in the instrument software according to these indications:

Setup	
Measurement mode	Standard
Measure from	Top
Measurement height	Must be adjusted
Measurement time	800 ms
Delay	1000 µs
Window time	400µs
Number of sequential windows	1
EnVision Nexus lamp	
Filter module	HTRF™/LANCE (330/665/620) Ref.: HH36794002 - Dual DELFLIA (330/615) Ref.: HH36794001 - Single
Energy source	Lamp
EnVision Nexus laser	
Filter module	HTRF™/LANCE (343/665/620) Ref.: HH36794003 - Dual DELFLIA (615) Ref.: HH36794006 - Single
Energy source	Laser

The pHSense assay operates with a single 620 nm detection mode. If a dual mirror is used, please ignore the 665 nm data.

pHSense readout for EnVision

EnVision microplate reader must be equipped with specific optical mirror and filters which enable the measurement of 620nm Europium emission.

Envision lamp Setup	
Cycle	2000 µs
Delay	1000 µs
Measurement height	Must be adjusted
Number of flash	100
Total time window	400 µs
Top mirror dual	
Top mirror	TRF Lance D407/D630 advanced dual mirror (#666) Ref.: 2105-4430 TRF LANCE/DELFIATM D400/D630 dual mirror (#662) Ref.: 2100-4160
Excitation filter	UV2 (TRF) 320 (#111) Ref.: 2100-5060 or UV (TRF) 340 (#101) Ref.: 2100-5010
Emission filter	APC 665 (#205) Ref.: 2100-5110
2 nd emission filter	CY5 620 (#118) Ref.: 2100-5240 or 615 (#203) Ref.: 2100-5090
Top mirror single	
Top mirror	TRF D407 advanced single mirror (#615) Ref.: 2105-4390 TRF LANCE/DELFIATM D400 single mirror (#412) Ref.: 2100-4170
Excitation filter	UV2 (TRF) 320 (#111) Ref.: 2100-5060 or UV (TRF) 340 (#101) Ref.: 2100-5010
Emission filter	CY5 620 (#118) Ref.: 2100-5240 or 615 (#203) Ref.: 2100-5090

EnVision microplate plate reader must be appropriately configured for pHSense readout by setting up the measurement conditions in the software according to these indications:

Envision laser Setup	
Cycle	16600 µs
Delay	1000 µs
Measurement height	Must be adjusted
Number of flash	50
Total time window	400 µs
Top mirror dual	
Top mirror	TRF LASER D407/D630 advanced dual mirror (#667) Ref.: 2105-4440 TRF LASER D400/D630 dual mirror (#446) Ref.: 2103-4290
Emission filter	APC 665 (#205) Ref.: 2100-5110
2 nd emission filter	CY5 620 (#118) Ref.: 2100-5240 or 615 (#203) Ref.: 2100-5090
Top mirror single	
Top mirror	TRF LASER D407 advanced single mirror (#616) Ref.: 2105-4400 TRF LANCE/DELFIATM D400 single mirror (#445) Ref.: 2103-4280
Emission filter	CY5 620 (#118) Ref.: 2100-5240 or 615 (#203) Ref.: 2100-5090

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

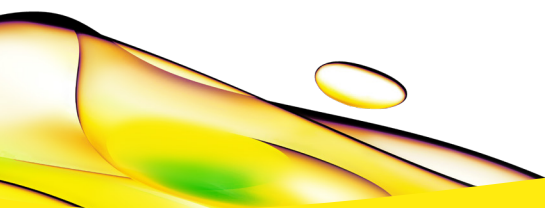
The pHSense assay operates with a single 620 nm detection mode. If a dual mirror is used, please ignore the 665 nm data.

pHSense readout for VICTOR Nivo

The VICTOR™ Nivo™ must be equipped with specific optical mirror and filters which enable the measurement of 620nm Europium emission.

The VICTOR Nivo reader must be appropriately configured for pHSense readout by setting up the measurement conditions in the software according to these indications:

Operation	TRF-EndPoint
Measurement type	Single label
Excitation filter	320/75 nm Ref.: HH35000947
Emission filter	620/10 nm Ref.: HH35000936 or 615/8 nm Ref.: HH35000935
Dichroic mirror	D400 Ref.: HH35000971
Delay time (μs)	1000
Emission time (μs)	400
Measurement direction	Top
Measurement time (ms)	500
Z-focus (mm)	Must be adjusted
Flash energy (mJ)	Low (10)



revvity