

HTRF setup recommendations for POLARstar OMEGA.



HTRF Europium cryptate donor / red acceptor readout setup recommendations for POLARstar OMEGA

POLARstar OMEGA is equipped with a specific optical device 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) allows the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

POLARstar OMEGA readers must be appropriately configured for HTRF™ readout by setting up the measurement conditions in the software according to the following indications:

Setup	
Excitation filter	337nm Ref.: 001-337TR (TR-Ex H)
Emission filters	620 (8.5) nm Ref.: 001-615TR 665 (10) nm Ref.: 001-665TR
Integration delay (lag time)	60 µs
Integration time	400 µs
Number of flashes	200
Optimal z-pos [§]	Volume and plate format dependant Adjustment to be done manually Select by default the following values: →10.5 for 384 well low-volume plate → 8.0 for 96 well half-area plate
Gain	2300 for 665 and 620

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 POLARstar OMEGA

POLARstar OMEGA is equipped with a specific optical device which enables the measurement of both 620 nm cryptate and 520 nm acceptor emissions. The ratio of the two fluorescence intensities 520/620 (acceptor/donor) allows the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

POLARstar OMEGA readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the software according to the following indications:

Setup	
Excitation filter	337 nm Ref.: 001-337TR (TR-Ex H)
Emission filters	620 (8.5) nm Ref.: 001-615TR 520 (10) nm Ref.: 001-520TR
Integration delay (lag time)	60 µs
Integration time	400 µs
Number of flashes	200
Optimal z-pos ⁵	Volume and plate format dependent Adjustment to be done manually Select by default the following values: → 10.5 for 384 well low-volume plate → 8.0 for 96 well half-area plate
Gain	2300 for 520 and 620

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

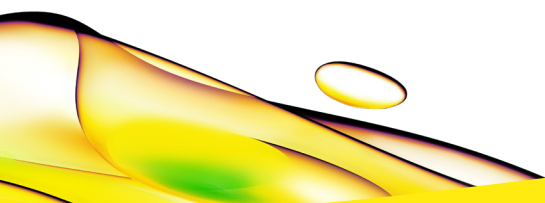
HTRF Terbium cryptate donor / red acceptor readout setup recommendations for POLARstar OMEGA

POLARstar OMEGA is equipped with a specific optical device, 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.

POLARstar OMEGA readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the software according to the following indications:

Setup	
Excitation filter	337 nm Ref.: 001-337TR (TR-Ex H)
Emission filters	620 (8.5) nm Ref.: 001-615TR 665 (10) nm Ref.: 001-665TR
Integration delay (lag time)	60 µs
Integration time	400 µs
Number of flashes	200
Optimal z-pos [§]	Volume and plate format dependent Adjustment to be done manually Select by default the following values: → 10.5 for 384 well low-volume plate → 8.0 for 96 well half-area plate
Gain	2300 for 665 and 620

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



revvity