

# HTRF setup recommendations for CLARIOstar & CLARIOstar Plus.



## HTRF Europium cryptate donor / red acceptor readout setup recommendations for CLARIOstar & CLARIOstar Plus

CLARIOstar readers can be equipped with monochromators or filters and only filter version is HTRF certified.

CLARIOstar 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.

CLARIOstar 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	EX-TR
Emission filters	620 (10) nm 665 (10) nm
Dichroic mirror	LP-TR
Integration delay (lag time)	60 μs
Integration time	400 μs
Number of flashes	200
Optimal z-pos <sup>5</sup>	Volume and plate format dependent
Gain	2600 for 665 and 620 with black plate 2400 for 665 and 620 with white plate

## HTRF Terbium cryptate donor / green acceptor readout setup recommendations for CLARIOstar & CLARIOstar Plus

**CLARIOstar readers can be equipped with monochromators or filters and only filter version is HTRF certified.**

CLARIOstar 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.

CLARIOstar 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	EX-TR
Emission filters	620 (10) nm 520 (10) nm
Dichroic mirror	LP-TR
Integration delay (lag time)	60 µs
Integration time	400 µs
Number of flashes	200
Optimal z-pos <sup>§</sup>	Volume and plate format dependent
Gain	2600 for 665 and 620 with black plate 2400 for 665 and 620 with white plate

## HTRF Terbium cryptate donor / red acceptor readout setup recommendations for CLARIOstar & CLARIOstar Plus

**CLARIOstar readers can be equipped with monochromators or filters and only filter version is HTRF certified.**

CLARIOstar 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.d

CLARIOstar 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	EX-TR
Emission filters	620 (10) nm 665 (10) nm
Dichroic mirror	LP-TR
Integration delay (lag time)	60 $\mu$ s
Integration time	400 $\mu$ s
Number of flashes	200
Optimal z-pos <sup>§</sup>	Volume and plate format dependent
Gain	2600 for 665 and 620 with black plate 2400 for 665 and 620 with white plate

