

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 [§]	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 [§]	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.

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 [§]	Volume and plate format dependent
Gain	2600 for 665 and 620 with black plate 2400 for 665 and 620 with white plate

