

HTRF® Europium cryptate donor / Red acceptor readout Setup recommendations for Synergy™ NEO

Synergy NEO must be 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.

Synergy NEO readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Gen5™ Reader Control and Data Analysis Software. In particular, these parameters should be entered as defined in the table below

Setup

Top filter cube	EX 330 / LUM EM 620 / 665 / LUM
Light source	Xenon flash
Lamp energy	Low (faster)
Delay	150 µs
Data time collection	500 µs
Measurement data point	10
Read height	plate format dependant 8.5mm for 384 wells low volume
Read speed	Normal
Gain	Automatic gain adjustment Autoscale

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 Synergy[™] NEO

Synergy NEO must be 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.

Synergy NEO readers must be appropriately configured for HTRF[®] readout by setting up the measurement conditions in the Gen5[™] Reader Control and Data Analysis Software. In particular, these parameters should be entered as defined in the table below

Setup

Top filter cube	EX 340 / LUM EM 620 / 520 / LUM
Light source	Xenon flash
Lamp energy	Low (faster)
Delay	150 µs
Data time collection	500 µs
Measurement data point	10
Read height	plate format dependant 8.5mm for 384 wells low volume
Read speed	Normal
Gain	Automatic gain adjustment Autoscale

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 Synergy™ NEO

Synergy NEO must be 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.

Synergy NEO readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Gen5™ Reader Control and Data Analysis Software. In particular, these parameters should be entered as defined in the table below

Setup

Top filter cube	EX 340 / LUM EM 620 / 665 / LUM
Light source	Xenon flash
Lamp energy	Low (faster)
Delay	150 µs
Data time collection	500 µs
Measurement data point	10
Read height	plate format dependant 8.5mm for 384 wells low volume
Read speed	Normal
Gain	Automatic gain adjustment Autoscale

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

