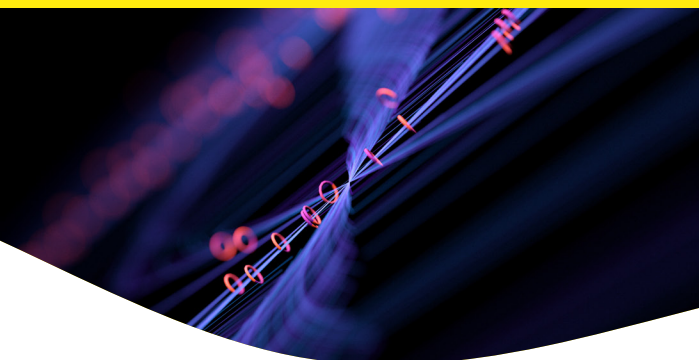


HTRF setup recommendations for Sense.



HTRF Europium cryptate donor / red acceptor readout setup recommendations for Sense

Two sequential measurements should be carried out: at 620 nm for the cryptate emission, and at 665 nm for the specific signal emitted by the acceptor (XL665 or d2). 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.

Sense 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	330 nm (80) nm
Emission filter Donor	620 (10) nm
Emission filter Acceptor	665 (7.5) nm
Lamp power	Super
Number of flashes	200
Delay	150 μs
Length	400 μs
Read mode	Top
Mirror	Automatic
Excitation aperture	Use default
Aperture Donor	8 for 384sv plate. For other plate type click on "Use default"
Aperture Acceptor	8 for 384sv plate. For other plate type click on "Use default"
Focus:	8 for 384sv plate. For other plate type click on "Use default"
Discriminator voltage	Use default
PMT voltage	Use default

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 Sense

Two sequential measurements should be carried out: at 620 nm for the cryptate emission, and at 520 nm for the specific signal emitted by the green acceptor. 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.

Sense 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	330 nm (80) nm
Emission filter Donor	620 (10) nm
Emission filter Acceptor	520 (10) nm
Lamp power	Super
Number of flashes	200
Delay	150 µs
Length	400 µs
Read mode	Top
Mirror	Automatic
Excitation aperture	Use default
Aperture Donor	8 for 384sv plate. For other plate type click on "Use default"
Aperture Acceptor	8 for 384sv plate. For other plate type click on "Use default"
Focus:	8 for 384sv plate. For other plate type click on "Use default"
Discriminator voltage	Use default
PMT voltage	Use default

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 Sense

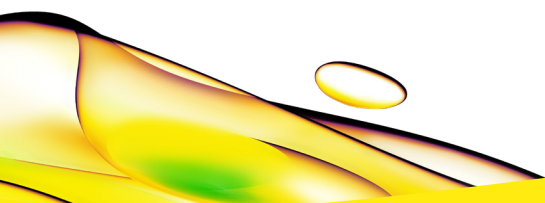
Two sequential measurements should be carried out: at 620 nm for the cryptate emission, and at 665 nm for the specific signal emitted by the acceptor (XL665 or d2).

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.

Sense 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	330 nm (80) nm
Emission filter Donor	620 (10) nm
Emission filter Acceptor	665 (7.5) nm
Lamp power	Super
Number of flashes	200
Delay	150 μ s
Length	400 μ s
Read mode	Top
Mirror	Automatic
Excitation aperture	Use default
Aperture Donor	8 for 384sv plate. For other plate type click on "Use default"
Aperture Acceptor	8 for 384sv plate. For other plate type click on "Use default"
Focus:	8 for 384sv plate. For other plate type click on "Use default"
Discriminator voltage	Use default
PMT voltage	Use default

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



revvity