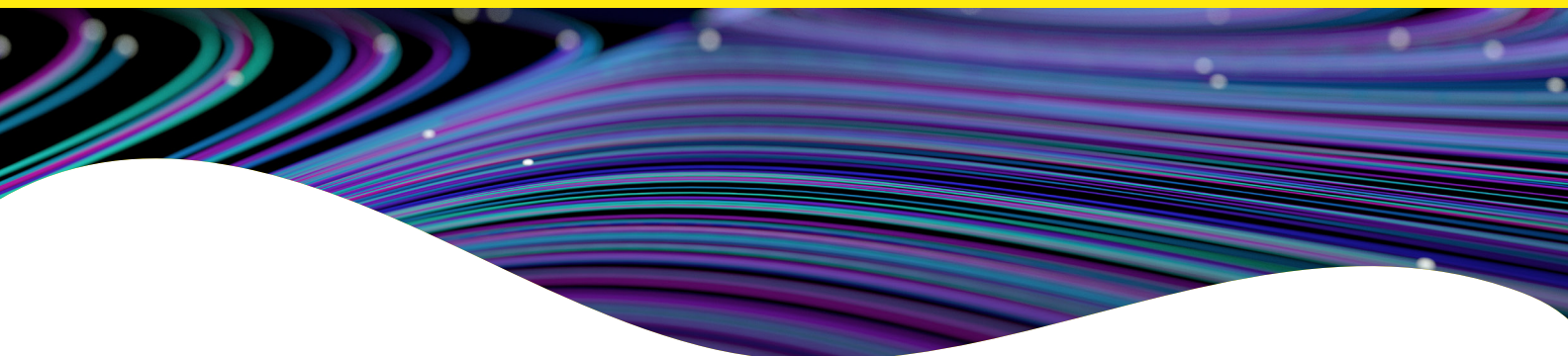


HTRF setup recommendations for Safire².



HTRF Europium cryptate donor / red acceptor readout setup recommendations for Safire²

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.

Safire²™ readers must be appropriately configured for HTRF™ readout by setting up the measurement conditions in the “multilabeling” function of Xfluor4 or Magellan software. In particular, these parameters should be entered as below. No special upgrade is required for HTRF readout, as it is a monochromator-based instrument:

	Measurement 1	Measurement 2
Excitation wavelength	317 nm	317 nm
Excitation bandwidth	20 nm	20 nm
Emission wavelength	620 nm	665 nm
Emission bandwidth	10 nm	10 nm
Number of reads	100	100
Lag time	60 μs	60 μs
Integration time	500 μs	500 μs
Gain	Optimal	Optimal
Z position	Optimal	Optimal

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 Safire²

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 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.

Safire² readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the "multilabeling" function of Xfluor4 or Magellan software. In particular, these parameters should be entered as below. No special upgrade is required for HTRF readout, as it is a monochromator-based instrument:

	Measurement 1	Measurement 2
Excitation wavelength	343 nm	343 nm
Excitation bandwidth	20 nm	20 nm
Emission wavelength	620 nm	520 nm
Emission bandwidth	10 nm	10 nm
Number of reads	100	100
Lag time	60 µs	60 µs
Integration time	500 µs	500 µs
Gain	Optimal	Optimal
Z position	Optimal	Optimal

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

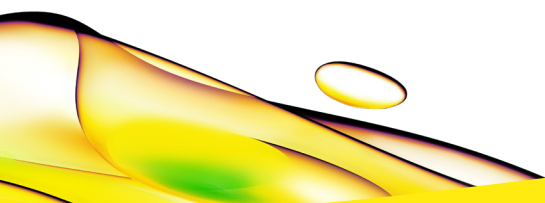
HTRF Terbium cryptate donor / red acceptor readout setup recommendations for Safire²

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.

Safire² readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the “multilabeling” function of Xfluor4 or Magellan software. In particular, these parameters should be entered as below. No special upgrade is required for HTRF readout, as it is a monochromator-based instrument:

	Measurement 1	Measurement 2
Excitation wavelength	340 nm	340 nm
Excitation bandwidth	20 nm	20 nm
Emission wavelength	620 nm	665 nm
Emission bandwidth	10 nm	10 nm
Number of reads	100	100
Lag time	60 µs	60 µs
Integration time	500 µs	500 µs
Gain	Optimal	Optimal
Z position	Optimal	Optimal

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



revvity