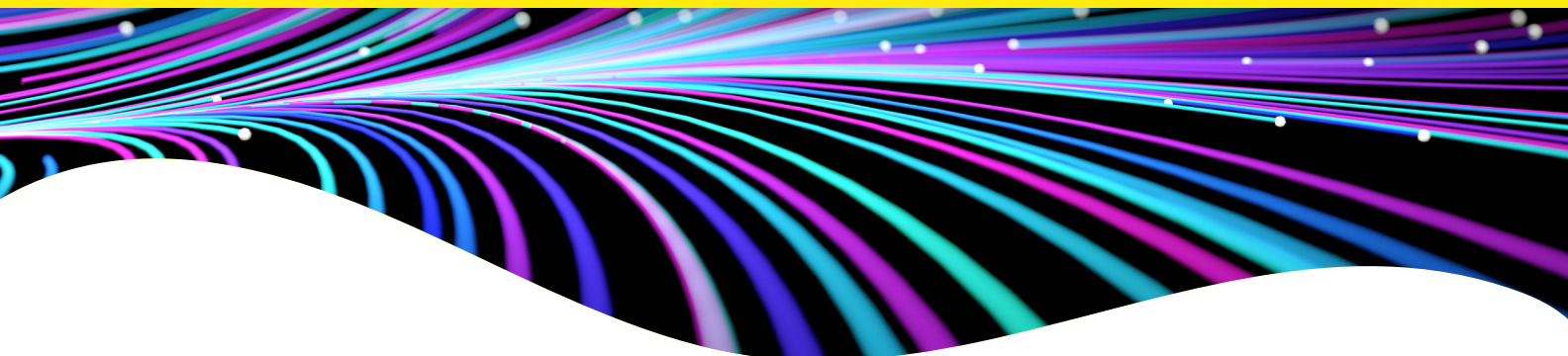


HTRF setup recommendations for Infinite M1000.



HTRF Europium cryptate donor / red acceptor readout setup recommendations for Infinite M1000

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.

Infinite® M1000 readers must be appropriately configured for HTRF™ readout by setting up the measurement conditions in the “multilabeling” function of Tecan i-Control™ 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:

	Label 1	Label 2
Measurement	Fluorescence intensity	Fluorescence intensity
Excitation wavelength	317 nm/20 nm	317 nm/20 nm
Emission wavelength	665 nm/10 nm	620 nm/10 nm
Mode	Top	Top
Flashes	Mode 2 [100Hz] : 100	Mode 2 [100Hz] : 100
Gain	Optimal	Optimal
Z position	Must be calculated from the well giving the highest signal	Must be calculated from the well giving the highest signal
Lag time	60 µs	60 µs
Integration time	500 µs	500 µs

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 Infinite M1000

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.

Infinite M1000 readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the “multilabeling” function of Tecan i-Control 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:

	Label 1	Label 2
Measurement	Fluorescence intensity	Fluorescence intensity
Excitation wavelength	340 nm/20 nm	340 nm/20 nm
Emission wavelength	520 nm/10 nm	620 nm/10 nm
Mode	Top	Top
Flashes	Mode 2 [100Hz] : 100	Mode 2 [100Hz] : 100
Gain	Optimal	Optimal
Z position	Must be calculated from the well giving the highest signal	Must be calculated from the well giving the highest signal
Lag time	60 µs	60 µs
Integration time	500 µs	500 µs

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 Infinite M1000

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.

Infinite M1000 readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the “multilabeling” function of Tecan i-Control 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:

	Label 1	Label 2
Measurement	Fluorescence intensity	Fluorescence intensity
Excitation wavelength	340 nm/20 nm	340 nm/20 nm
Emission wavelength	665 nm/10 nm	620 nm/10 nm
Mode	Top	Top
Flashes	Mode 2 [100Hz] : 100	Mode 2 [100Hz] : 100
Gain	Optimal	Optimal
Z position	Must be calculated from the well giving the highest signal	Must be calculated from the well giving the highest signal
Lag time	60 µs	60 µs
Integration time	500 µs	500 µs

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

