

HTRF setup recommendations for Infinite F Plex and F Nano+.



HTRF Europium cryptate donor / red acceptor readout setup recommendations for Infinite F Plex and F Nano+

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.

The Infinite® F Plex and F Nano+ must be equipped with the HTRF™ module. Infinite F version readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the Tecan i-Control™ software. These parameters should be entered as defined in the table below.

	Measurement 1	Measurement 2
Excitation filter	320(25) nm Ref: HF370	320(25) nm Ref: HF370
Emission filter	620(20) nm Ref: HJ471	665(8) nm Ref: HB045
Mode	Top	Top
Mirror	Dichroic 510	Dichroic 510
Number of flashes	25	25
Settle time	0	0
Gain	Optimal	Optimal
Lag time	150 µs	150 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 F Plex and F Nano+

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.

The Infinite F Plex and F Nano+ must be equipped with the HTRF module. Infinite F version readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the TECAN i-Control software. These parameters should be entered as defined in the table below:

	Measurement 1	Measurement 2
Excitation filter	340(20) nm Ref.: HF113	340(20) nm Ref.: HF113
Emission filter	620(20) nm Ref.: HJ471	665(8) nm Ref.: HB045
Mode	Top	Top
Mirror	Dichroic 510	Dichroic 510
Number of flashes	25	25
Settle time	0	0
Gain	Optimal	Optimal
Lag time	150 µs	150 µs
Integration time	500 µs	500 µs

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

