

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

Excitation filter	320 (25) nm	Ref: HF370
Emission filter	620 (20) nm	Ref: HJ471
Mode	Top	
Mirror	Dichroic 510	
Number of flashes	25	
Settle time	0	
Gain	Optimal	
Lag time	150 µs	
Integration time	500 µs	

Measurement 2

Excitation filter	320 (25) nm	Ref.: HF370
Emission filter	665 (8) nm	Ref.: HB045
Mode	Top	
Mirror	Dichroic 510	
Number of flashes	25	
Settle time	0	
Gain	Optimal	
Lag time	150 µs	
Integration time	500 µs	

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

