

## 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

Excitation filter	340 (20) nm	Ref.: HF113
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	340 (20) nm	Ref.: HF113
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.

