

HTRF® Europium cryptate donor / Red acceptor readout Setup recommendations for Infinite® F200 PRO

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® F200 PRO must be equipped with the HTRF® module. Infinite® F200 PRO readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Tecan i-Control™ software. In particular, these parameters should be entered as defined in the table below.

Measurement 1

Excitation filter	320 (25) nm	Ref: 8C138
Emission filter	620 (10) nm	Ref: 6F041
Mirror	Dichroic 510	
Lag time	150 µs	
Integration time	500 µs	
Number of reads	10	
Gain	Optimal	
Z	Can be calculated on the well giving the highest signal	

Measurement 2

Excitation filter	320 (25) nm	Ref.: 8C138
Emission filter	665 (8.5) nm	Ref.: 9E336
Mirror	Dichroic 510	
Lag time	150 µs	
Integration time	500 µs	
Number of reads	10	
Gain	Optimal	
Z	Can be calculated on the well giving the highest signal	

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® F200 PRO

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.

The Infinite® F200 PRO must be equipped with the HTRF® module. Infinite® F200 PRO readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Tecan i-Control™ software. In particular, these parameters should be entered as defined in the table below:

Measurement 1

Excitation filter	340 (35) nm	Ref.: 9E095
Emission filter	620 (10) nm	Ref.: 6F041
Mirror	Dichroic 510	
Lag time	150 µs	
Integration time	500 µs	
Number of reads	10	
Gain	Optimal	
Z	Can be calculated on the well giving the highest signal	

Measurement 2

Excitation filter	340 (35) nm	Ref.: 9E095
Emission filter	520 (10) nm	Ref.: 7F90
Mirror	Dichroic 510	
Lag time	150 µs	
Integration time	500 µs	
Number of reads	10	
Gain	Optimal	
Z	Can be calculated on the well giving the highest signal	

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® F200 PRO

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® F200 PRO must be equipped with the HTRF® module. Infinite® F200 PRO readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the TECAN i-Control™ software. In particular, these parameters should be entered as defined in the table below:

Measurement 1

Excitation filter	340 (35) nm	Ref.: 9E095
Emission filter	620 (10) nm	Ref.: 6F041
Mirror	Dichroic 510	
Lag time	150 µs	
Integration time	500 µs	
Number of reads	10	
Gain	Optimal	
Z	Can be calculated on the well giving the highest signal	

Measurement 2

Excitation filter	340 (35) nm	Ref.: 9E095
Emission filter	665 (8.5) nm	Ref.: 9E336
Mirror	Dichroic 510	
Lag time	150 µs	
Integration time	500 µs	
Number of reads	10	
Gain	Optimal	
Z	Can be calculated on the well giving the highest signal	

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

