

HTRF setup recommendations for Infinite F200 PRO.



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	Measurement 2
Excitation filter	320 (25) nm Ref: 8C138	320 (25) nm Ref: 8C138
Emission filter	620 (10) nm Ref: 6F041	665 (8.5) nm Ref: 9E336
Mirror	Dichroic 510	Dichroic 510
Lag time	150 µs	150 µs
Integration time	500 µs	500 µs
Number of reads	10	10
Gain	Optimal	Optimal
Z	Can be calculated on the well giving the highest signal	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	Measurement 2
Excitation filter	340 (35) nm Ref: 9E095	340 (35) nm Ref: 9E095
Emission filter	620 (10) nm Ref: 6F041	520 (10) nm Ref: 7F90
Mirror	Dichroic 510	Dichroic 510
Lag time	150 µs	150 µs
Integration time	500 µs	500 µs
Number of reads	10	10
Gain	Optimal	Optimal
Z	Can be calculated on the well giving the highest signal	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.

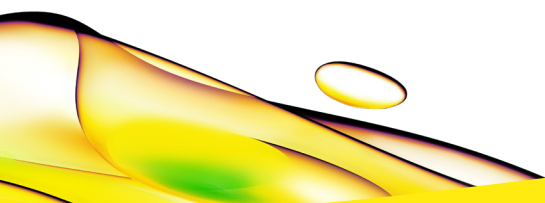
HTRF Terbium cryptate donor / red acceptor readout setup recommendations for Infinite F200 PRO

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	Measurement 2
Excitation filter	340 (35) nm Ref: 9E095	340 (35) nm Ref: 9E095
Emission filter	620 (10) nm Ref: 6F041	665 (8.5) nm Ref: 9E336
Mirror	Dichroic 510	Dichroic 510
Lag time	150 µs	150 µs
Integration time	500 µs	500 µs
Number of reads	10	10
Gain	Optimal	Optimal
Z	Can be calculated on the well giving the highest signal	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.



revvity