

HTRF[®] Tb readout - Set up recommendations for Infinite[®] M1000 readers

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). A ratio of the two fluorescence intensities* (acceptor/donor) then allows the calculation of Delta F (%), i.e. the relative energy transfer rate for each data point.

Infinite[®] M1000 readers must be appropriately configured for HTRF[®] readout by setting up the measurement conditions in the "multilabeling" function of i-Control software. In particular, these parameters should be entered as below. No special upgrade is required for HTRF[®] readout, as it is a monochromator-based instrument.

Measurement 1

Excitation wavelength : 340 nm
Excitation bandwidth : 20 nm
Emission wavelength : 620 nm
Emission bandwidth : 10 nm
Number of reads : 100
Lag time : 60 µs
Integration time : 500 µs
Optimal Gain
Optimal Z position

Measurement 2

Excitation wavelength : 340 nm
Excitation bandwidth : 20 nm
Emission wavelength : 665 nm
Emission bandwidth : 10 nm
Number of reads : 100
Lag time : 60 µs
Integration time : 500 µs
Optimal Gain
Optimal Z position

WHITE plates only can be used on this reader!

** The fluorescence ratio is a correction method developed by CIS bio international with an application limited to the use of HTRF[®] reagents and technology, and for which CIS bio international has granted a licence to Tecan. The method is covered by the US patent 5,527,684 and its foreign equivalents.*