

HTRF® Europium cryptate donor / Red acceptor readout Setup recommendations for Infinite® M1000

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.

Infinite® M1000 readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the "multilabeling" function of Tecan 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:

	Label 1	Label 2
Measurement	Fluorescence intensity	Fluorescence intensity
Excitation wavelength	317nm / 20nm	317nm / 20nm
Emission wavelength	665nm / 10nm	620nm / 10nm
Mode	Top	Top
Flashes	Mode 2 [100Hz] : 100	Mode 2 [100Hz] : 100
Gain	Optimal	Optimal
Z position	Must be calculated from the well giving the highest signal	Must be calculated from the well giving the highest signal
Lag time	60µs	60µs
Integration time	500µs	500µs

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



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