

## HTRF<sup>®</sup> Terbium cryptate donor / Red acceptor readout Setup recommendations for GENios Pro<sup>™</sup>

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.

GENios Pro<sup>™</sup> readers must be appropriately configured for HTRF<sup>®</sup> readout by setting up the measurement conditions in the "multilabeling" function of Xfluo4 software. In particular, these parameters should be entered as defined in table below following the installation of the Tecan HTRF<sup>®</sup> upgrade kit (Tecan #B122175) on GENios Pro<sup>™</sup>.

### Measurement 1

Excitation filter	340 (20) nm	Ref.: 30000405
Emission filter	620 (10) nm	Ref.: 30002292
Mirror	Dichroic 3 (e.g. FI 96)	
Number of flashes	10	
Lag time	150 µs	
Integration time	500 µs	
Gain	Optimal	
Z position	Optimal	

### Measurement 2

Excitation filter	340 (20) nm	Ref.: 30000405
Emission filter	665 (8.5) nm	Ref.: 30007518
Mirror	Dichroic 3 (e.g. FI 96)	
Number of flashes	10	
Lag time	150 µs	
Integration time	500 µs	
Gain	Optimal	
Z position	Optimal	



*\*The fluorescence ratio is a correction method developed by Cisbio Bioassays with an application limited to the use of HTRF<sup>®</sup> 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.*