

HTRF® Terbium cryptate donor / Red acceptor readout Setup recommendations for Varioskan Lux reader

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.

Varioskan Lux readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the software according to the following indications:

Setup	
Excitation filter	334nm (40)nm
Emission filter Donor	620 (10) nm
Emission filter Acceptor	665 (10) nm
Delay time	60 µs
Integration time	200µs
Measurement time	1000ms
Dynamic range	Automatic

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

908-018-125



**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 BMG LABTECH. The method is covered by the US patent 5,527,684 and its foreign equivalents.*