

HTRF® Europium 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.**



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

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 green 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.

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

Delay time	60µs
Integration time	200µs
Measurement time	1000ms
Dynamic range	Automatic

- **This reader only allows high performance HTRF measurement when assays are run**



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 μ s
Integration time	200 μ s
Measurement time	1000ms
Dynamic range	Automatic

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

