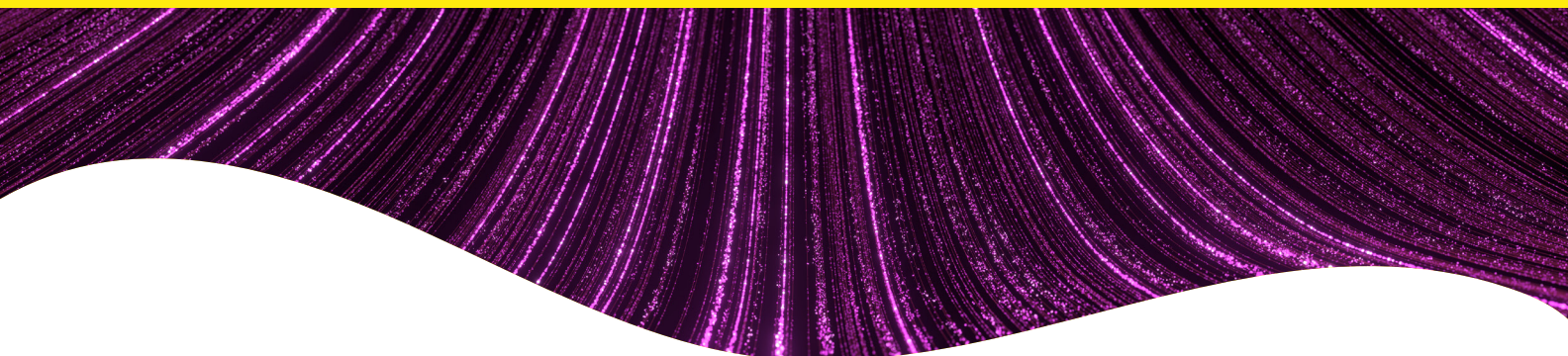


HTRF setup recommendations for Varioskan Lux reader.



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	334 nm (40) nm
Emission filter donor	620 (10) nm
Emission filter acceptor	665 (10) nm
Delay time	60 μs
Integration time	200 μs
Measurement time	1000 ms
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	60 µs
Integration time	200 µs
Measurement time	1000 ms
Dynamic range	Automatic

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

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	334 nm (40) nm
Emission filter donor	620 (10) nm
Emission filter acceptor	665 (10) nm
Delay time	60 µs
Integration time	200 µs
Measurement time	1000 ms
Dynamic range	Automatic

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

