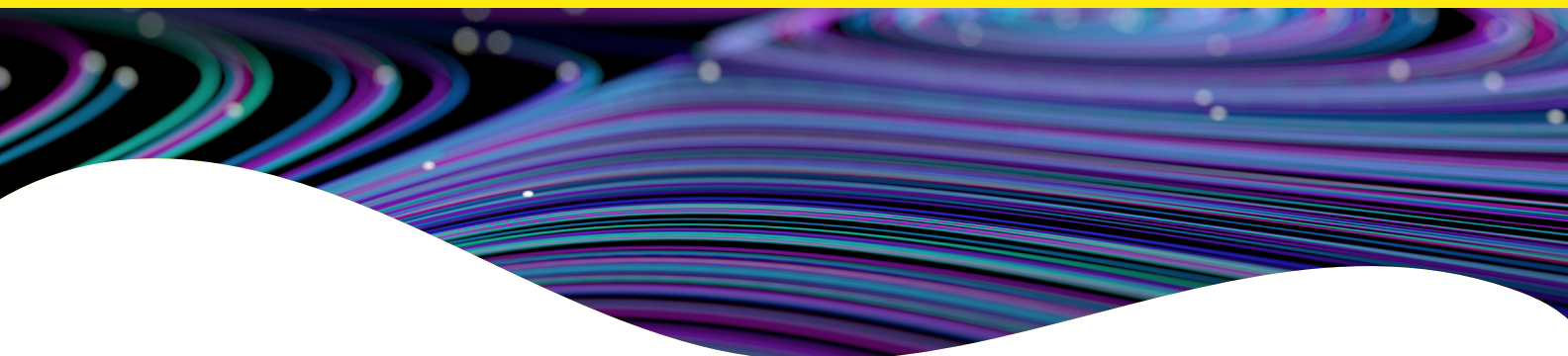


HTRF setup recommendations for Spark 10M filter / filter (FF) configuration.



HTRF Europium cryptate donor / red acceptor readout setup recommendations for Spark 10M Filter / Filter (FF) configuration

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.

The Spark 10M must be equipped with the HTRF™ module. Spark 10M readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the Tecan i-Control™ software. In particular, these parameters should be entered as defined in the table below.

	Measurement 1	Measurement 2
Excitation filter	320 (25) nm	320 (25) nm
Emission filter	620 (10) nm	665 (8) nm
Mirror	Dichroic 510	Dichroic 510
Lag time	100 µs	100 µs
Integration time	400 µs	400 µs
Flashes	75	75
Gain	Optimal gain	Optimal gain
Z	Can be calculated on the well giving the highest signal	Can be calculated on the well giving the highest signal

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

HTRF Terbium cryptate donor / green acceptor readout setup recommendations for Spark 10M Filter / Filter (FF) configuration

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 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.

The Spark 10M must be equipped with the HTRF module. Spark 10M readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the Tecan i-Control software. In particular, these parameters should be entered as defined in the table below.

	Measurement 1	Measurement 2
Excitation filter	340 (35) nm	340 (35) nm
Emission filter	620 (10) nm	520 (10) nm
Mirror	Dichroic 510	Dichroic 510
Lag time	100 µs	100 µs
Integration time	200 µs	200 µs
Flashes	75	75
Gain	Optimal gain	Optimal gain
Z	Can be calculated on the well giving the highest signal	Can be calculated on the well giving the highest signal

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

HTRF Terbium cryptate donor / red acceptor readout setup recommendations for Spark 10M Filter / Filter (FF) configuration

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.

The Spark 10M must be equipped with the HTRF module. Spark 10M readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the Tecan i-Control software. In particular, these parameters should be entered as defined in the table below.

	Measurement 1	Measurement 2
Excitation filter	340 (35) nm	340 (35) nm
Emission filter	620 (10) nm	665 (8) nm
Mirror	Dichroic 510	Dichroic 510
Lag time	100 µs	100 µs
Integration time	200 µs	200 µs
Flashes	75	75
Gain	Optimal gain	Optimal gain
Z	Can be calculated on the well giving the highest signal	Can be calculated on the well giving the highest signal

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

HTRF Europium cryptate donor / red acceptor readout setup recommendations for Spark 10M Monochromator/ Filter (MF) configuration

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.

The Spark 10M must be equipped with the HTRF module. Spark 10M readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the Tecan i-Control software. In particular, these parameters should be entered as defined in the table below.

	Measurement 1	Measurement 2
Excitation Monochromator	320 (20) nm	320 (25) nm
Emission filter	620 (10) nm	665 (8) nm
Mirror	Dichroic 510	Dichroic 510
Lag time	100 µs	100 µs
Integration time	400 µs	400 µs
Flashes	75	75
Gain	Optimal gain	Optimal gain
Z	Can be calculated on the well giving the highest signal	Can be calculated on the well giving the highest signal

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

HTRF Terbium cryptate donor / green acceptor readout setup recommendations for Spark 10M Monochromator / Filter (MF) configuration

Two sequential measurements should be carried out: at 620 nm for the cryptate emission, and at 520nm for the specific signal emitted by the 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.

The Spark 10M must be equipped with the HTRF module. Spark 10M readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the Tecan i-Control software. In particular, these parameters should be entered as defined in the table below.

	Measurement 1	Measurement 2
Excitation Monochromator	340 (35) nm	340 (35) nm
Emission filter	620 (10) nm	520 (10) nm
Mirror	Dichroic 510	Dichroic 510
Lag time	100 µs	100 µs
Integration time	200 µs	200 µs
Flashes	75	75
Gain	Optimal gain	Optimal gain
Z	Can be calculated on the well giving the highest signal	Can be calculated on the well giving the highest signal

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

HTRF Terbium cryptate donor / red acceptor readout setup recommendations for Spark 10M Monochromator/ Filter (MF) configuration

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.

The Spark 10M must be equipped with the HTRF module. Spark 10M readers must be appropriately configured for HTRF readout by setting up the measurement conditions in the Tecan i-Control software. In particular, these parameters should be entered as defined in the table below.

	Measurement 1	Measurement 2
Excitation Monochromator	340 (35) nm	340 (35) nm
Emission filter	620 (10) nm	665 (8) nm
Mirror	Dichroic 510	Dichroic 510
Lag time	100 μ s	100 μ s
Integration time	200 μ s	200 μ s
Flashes	75	75
Gain	Optimal gain	Optimal gain
Z	Can be calculated on the well giving the highest signal	Can be calculated on the well giving the highest signal

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

