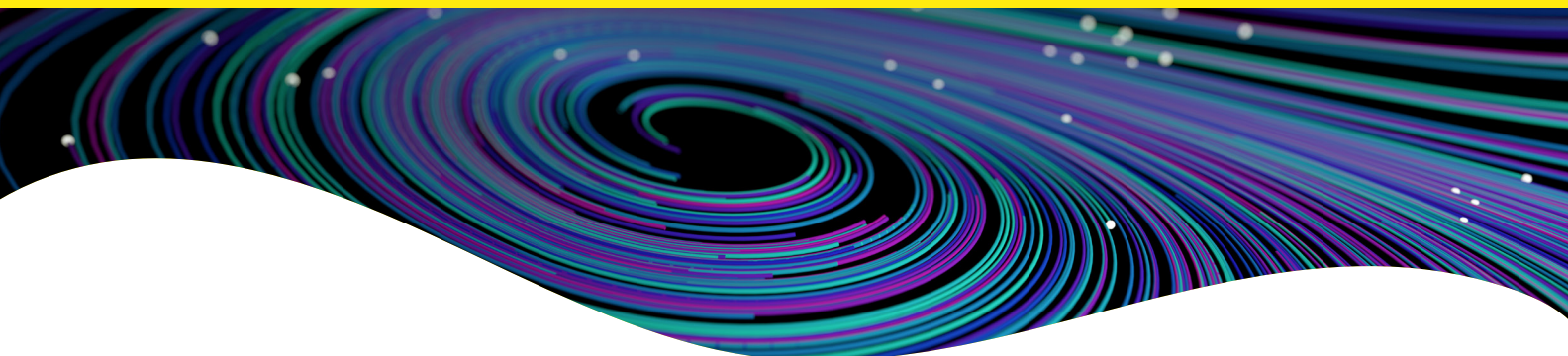


# HTRF setup recommendations for Spark 20M.



## HTRF Europium cryptate donor / red acceptor readout setup recommendations for Spark 20M

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 20M must be equipped with the HTRF™ module. Spark 20M 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.

Configurations : (Excitation/ Emission)	Configuration of head optical for excitation and emission selection			
	Filter/Filter (F/F)	Filter/Monochromator (F/M)	Monochromator/Filter (M/F)	Monochromator/Monochromator (M/M)
Excitation wavelength	320 nm/25 nm	320 nm/25 nm	320 nm/20 nm	320 nm/20 nm
Acceptor emission wavelength	665 nm/8 nm	665 nm/10 nm	665 nm/8 nm	665 nm/10 nm
Donor emission wavelength	620 nm/10 nm	620 nm/10 nm	620 nm/10 nm	620 nm/10 nm
Flashes	50	100	100	100
Lag time	100 µs	100 µs	100 µs	100 µs
Integration time	300 µs	300 µs	300 µs	300 µs
Mirror	510 dichroic	510 dichroic	510 dichroic	510 dichroic
Gain	Optimal gain	Optimal gain	Optimal gain	Optimal gain
Z optimization	Calculated from well giving the highest signal			
Plate color	Black and white	Black and white	Black and white	White plate only

## HTRF Terbium cryptate donor / green acceptor readout setup recommendations for Spark 20M

Two sequential measurements should be carried out at 620nm for the cryptate emission, and at 520nm 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.

The spark 20M must be equipped with the HTRF module. Spark 20M 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.

Configurations : (Excitation/ Emission)	Configuration of head optical for excitation and emission selection			
	Filter/Filter (F/F)	Filter/Monochromator (F/M)	Monochromator/Filter (M/F)	Monochromator/Monochromator (M/M)
Excitation wavelength	340 nm/35 nm	340 nm/35 nm	340 nm/25 nm	340 nm/25 nm
Acceptor emission wavelength	520 nm/10 nm	520 nm/10 nm	520 nm/10 nm	520 nm/10 nm
Donor emission wavelength	620 nm/10 nm	620 nm/10 nm	620 nm/10 nm	620 nm/10 nm
Flashes	50	100	100	100
Lag time	100 µs	100 µs	100 µs	100 µs
Integration time	300 µs	300 µs	300 µs	300 µs
Mirror	510 dichroic	510 dichroic	510 dichroic	510 dichroic
Gain	Optimal gain	Optimal gain	Optimal gain	Optimal gain
Z optimization	Calculated from well giving the highest signal			
Plate color	Black and white	Black and white	Black and white	<b>White plate only</b>

## HTRF Terbium cryptate donor / red acceptor readout setup recommendations for Spark 20M

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 20M must be equipped with the HTRF module. Spark 20M 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.

Configuration of head optical for excitation and emission selection				
Configurations : (Excitation/ Emission)	Filter/Filter (F/F)	Filter/Monochromator (F/M)	Monochromator/Filter (M/F)	Monochromator/Monochromator (M/M)
Excitation wavelength	340 nm/35 nm	340 nm/35 nm	340 nm/25 nm	340 nm/25 nm
Acceptor emission wavelength	665 nm/8 nm	665 nm/10 nm	665 nm/8 nm	665 nm/10 nm
Donor emission wavelength	620 nm/10 nm	620 nm/10 nm	620 nm/10 nm	620 nm/10 nm
Flashes	50	100	100	100
Lag time	100 µs	100 µs	100 µs	100 µs
Integration time	300 µs	300 µs	300 µs	300 µs
Mirror	510 dichroic	510 dichroic	510 dichroic	510 dichroic
Gain	Optimal gain	Optimal gain	Optimal gain	Optimal gain
Z optimization	Calculated from well giving the highest signal			
Plate color	Black and white	Black and white	Black and white	White plate only

