

HTRF® Europium cryptate donor / Red acceptor readout Setup recommendations for Spark using filters

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) to calculate the 665/620 signal ratio.

Spark must be optically equipped for HTRF® readout. Spark readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Spark Control™ software.

Note : Setup recommendations below are for Filter (Ex) / Filter (Em) configuration only.

Measurement 1

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

Measurement 2

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

For others configurations involving monochromator (FM,MF,MM), please contact Tecan to determine the version of your Spark reader (standard or enhanced). Then refer to other setups presented in the same web section.



HTRF® Terbium cryptate donor / Green acceptor readout Setup recommendations for Spark using filters

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 to calculate the 520/620 signal ratio.

Spark must be optically equipped for HTRF® readout. Spark readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Spark Control™ software.

Note : Setup recommendations below are for Filter (Ex) / Filter (Em) configuration only.

Measurement 1

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

Measurement 2

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

For others configurations involving monochromator (FM,MF,MM), please contact Tecan to determine the version of your Spark reader (standard or enhanced). Then refer to other setups presented in the same web section.



HTRF® Terbium cryptate donor / Red acceptor readout Setup recommendations for Spark using filters

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) to calculate the 665/620 signal ratio.

Spark must be optically equipped for HTRF® readout. Spark readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Spark Control™ software.

Note : Setup recommendations below are for Filter (Ex) / Filter (Em) configuration only.

Measurement 1

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

Measurement 2

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

For others configurations involving monochromator (FM,MF,MM), please contact Tecan to determine the version of your Spark reader (standard or enhanced). Then refer to other setups presented in the same web section.



**HTRF® Europium cryptate donor / Red acceptor readout
Setup recommendations for Spark (enhanced version) using
the monochromator**

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) to calculate the 665/620 signal ratio.

The spark must be optically equipped for HTRF® readout. Spark readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Spark Control™ software. In particular, these parameters should be entered as defined in the table below.

Note : For configuration involving monochromator, please contact Tecan to determine the version of your Spark reader : Standard or Enhanced. **Only Enhanced version is HTRF® compatible**

Configuration of optical head for excitation and emission selection			
Configuration : (Excitation/Emission)	Filter/Monochromator (F/M)	Monochromator/Filter (M/F)	Monochromator/Monochromator (M/M)
Excitation wavelength	320nm/20nm	320nm/25nm	320nm/25nm
Acceptor emission wavelength	665nm/10nm	665nm/8nm	665nm/10nm
Donor emission wavelength	620nm/10nm	620nm/10nm	620nm/10nm
Flashes	50		
Lag time	100µs		
Integration time	400µs		
Mirror	510 dichroic		
Gain	Optimal gain		
Z optimization	Calculated from well giving the highest signal		
Plate color	White plate only		



**HTRF® Terbium cryptate donor / Green acceptor readout
Setup recommendations for Spark (enhanced version) using
the monochromator**

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 to calculate the 665/620 signal ratio.

The spark must be optically equipped for HTRF® readout. Spark readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Spark Control™ software. In particular, these parameters should be entered as defined in the table below.

Note : For configuration involving monochromator, please contact Tecan to determine the version of your Spark reader : Standard or Enhanced. **Only Enhanced version is HTRF® compatible.**

Configuration of optical head for excitation and emission selection			
Configuration : (Excitation/Emission)	Filter/Monochromator (F/M)	Monochromator/Filter (M/F)	Monochromator/Monochromator (M/M)
Excitation wavelength	340nm/35nm	340nm/30nm	340nm/30nm
Acceptor emission wavelength	520nm/10nm	520nm/10nm	520nm/10nm
Donor emission wavelength	620nm/10nm	620nm/10nm	620nm/10nm
Flashes	50		
Lag time	100µs		
Integration time	400µs		
Mirror	510 dichroic		
Gain	Optimal gain		
Z optimization	Calculated from well giving the highest signal		
Plate color	White plate only		



**HTRF® Terbium cryptate donor / Red acceptor readout
Setup recommendations for Spark (enhanced version) using
the monochromator**

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) to calculate the 665/620 signal ratio.

The spark must be optically equipped for HTRF® readout. Spark readers must be appropriately configured for HTRF® readout by setting up the measurement conditions in the Spark Control™ software. In particular, these parameters should be entered as defined in the table below.

Note : For configuration involving monochromator, please contact Tecan to determine the version of your Spark reader : Standard or Enhanced. **Only Enhanced version is HTRF® compatible.**

Configuration of optical head for excitation and emission selection			
Configuration : (Excitation/Emission)	Filter/Monochromator (F/M)	Monochromator/Filter (M/F)	Monochromator/Monochromator (M/M)
Excitation wavelength	340nm/35nm	340nm/35nm	340nm/35nm
Acceptor emission wavelength	665nm/10nm	665nm/8nm	665nm/10nm
Donor emission wavelength	620nm/10nm	620nm/10nm	620nm/10nm
Flashes	50		
Lag time	100µs		
Integration time	400µs		
Mirror	510 dichroic		
Gain	Optimal gain		
Z optimization	Calculated from well giving the highest signal		
Plate color	White plate only		

