

HTRF setup recommendations for Mithras LB 940.



HTRF Europium cryptate donor / red acceptor readout setup recommendations for Mithras LB 940

The Mithras LB940 reader must be equipped with the TR-FRET reading module which includes the necessary optical components for HTRF™ readout. Two sequential readings at 620 nm and 665 nm emission wavelengths are performed. The ratio of the fluorescence intensities 665/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

The Mithras LB940 operating software comes with preset ready-to-use parameter files for HTRF measurements including the ratio calculation. The recommended settings are defined under the TR-Fluorescence protocol as described below:

	Measurement 1	Measurement 2
Excitation filter	D320 (40) Ref.: 52733	D320 (40) Ref.: 52733
Emission filter	D620 (TRF) (10) Ref.: 47731	D665 (TRF) (7.5) Ref.: 52544
Lamp energy	100	100
Cycle time	2000 µs	2000 µs
Delay time	50 µs	50 µs
Reading time	300 µs	300 µs
Counting time	1s Optimal	1s Optimal
Operation mode	by plate	by plate

HTRF Terbium cryptate donor / green acceptor readout setup recommendations for Mithras LB 940

The Mithras LB940 reader must be equipped with the TR-FRET reading module which includes the necessary optical components for HTRF readout. Two sequential readings at 620 nm and 520 nm emission wavelengths are performed. The ratio of the fluorescence intensities 520/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

The Mithras LB940 operating software comes with preset ready-to-use parameter files for HTRF measurements including the ratio calculation. The recommended settings are defined under the TR-Fluorescence protocol as described below:

	Measurement 1	Measurement 2
Excitation filter	D340 / 26 Ref.: 54083	D340 / 26 Ref.: 54083
Emission filter	D620 (TRF) (10) Ref.: 47731	D520 (TRF) (10) Ref.: 38836
Lamp energy	100	100
Cycle time	2000 μ s	2000 μ s
Delay time	50 μ s	50 μ s
Reading time	400 μ s	300 μ s
Counting time	1s Optimal	1s Optimal
Operation mode	By plate	By plate

HTRF Terbium cryptate donor / red acceptor readout setup recommendations for Mithras LB 940

The Mithras LB940 reader must be equipped with the TR-FRET reading module which includes the necessary optical components for HTRF readout. Two sequential readings at 620 nm and 665 nm emission wavelengths are performed. The ratio of the fluorescence intensities 665/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

The Mithras LB940 operating software comes with preset ready-to-use parameter files for HTRF measurements including the ratio calculation. The recommended settings are defined under the TR-Fluorescence protocol as described below:

	Measurement 1	Measurement 2
Excitation filter	D340 / 26 Ref.: 54083	D340 / 26 Ref.: 54083
Emission filter	D620 (TRF) (10) Ref.: 47731	D665 (TRF) (7.5) Ref.: 52544
Lamp energy	100	100
Cycle time	2000 μ s	2000 μ s
Delay time	50 μ s	50 μ s
Reading time	400 μ s	400 μ s
Counting time	1s Optimal	1s Optimal
Operation mode	By plate	By plate

