

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

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 must be equipped with the HTRF® module. Spark 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.

Caution : Only the configuration Filter (Ex) / Filter (Em) is HTRF® compatible

Measurement 1

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

