

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

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

