

HTRF[®] Tb readout - Set up recommendations for Synergy 4[™] readers

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. A ratio of the two fluorescence intensities* (acceptor/donor) then allows the calculation of Delta F (%), i.e. the relative energy transfer rate for each data point.

Synergy 4 readers must be appropriately configured for HTRF[®] Tb readout by setting up the measurement conditions in the Gen5[™] Reader Control and Data Analysis Software. In particular, these parameters should be entered as defined in the table below. The Synergy 4 must be equipped with the TRF module.

HTRF assays must be read using the filter-based detection mode only. The monochromator mode is not HTRF compatible.

USE WHITE PLATES ONLY

Measurement 1

| | |
|--------------------------|---|
| Excitation filter | : 340 (30) nm |
| Emission filter | : 620 (10) nm |
| Optics Position | : Top 400 nm |
| Number of Flashes | : 10 |
| Lag time | : 100µs |
| Integration time | : 300µs |
| Sensitivity | : Value to optimise on the well having the highest signal in order to reach 50000 counts. |
| Z | : Take the default value given in the software. |

Measurement 2

| | |
|--------------------------|---|
| Excitation filter | : 340 (30) nm |
| Emission filter | : 665 (8) nm |
| Optics Position | : Top 400 nm |
| Number of Flashes | : 10 |
| Lag time | : 100µs |
| Integration time | : 300µs |
| Sensitivity | : Value to optimise on the well having the highest signal in order to reach 50000 counts. |
| Z | : Take the default value given in the software. |

* The fluorescence ratio is a correction method developed by CIS bio international with an application limited to the use of HTRF reagents and technology, and for which CIS bio international has granted a licence to BioTek. The method is covered by the US patent 5,527,684 and its foreign equivalents.