

## HTRF® Terbium cryptate donor / Green acceptor readout Setup recommendations for Cytation 1

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.

Cytation 1 readers must be appropriately configured for HTRF® 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.

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

### Measurement 1

Excitation filter	340 (30) nm
Emission filter	620 (10) nm
Optics position	top 400
Number of flashes	10
Lag time	100 µs
Integration time	300 µs
Sensitivity	Read the plate, select the well with the highest signal, and set its value at 50,000 counts
Z	Select the default value given in the software

### Measurement 2

Excitation filter	340 (30) nm
Emission filter	520 (10) nm
Optics position	top 400
Number of flashes	10
Lag time	100 µs
Integration time	300 µs
Sensitivity	Read the plate, select the well with the highest signal, and set its value at 50,000 counts
Z	Select the default value given in the software

**This reader only allows high performance HTRF measurement when assays are run in WHITE plates.**

