

## HTRF® Europium cryptate donor / Red acceptor readout Setup recommendations for Synergy™ NEO2

Synergy NEO2 must be equipped with a specific optical device, which enables the simultaneous measurement of both 620 nm cryptate and 665 nm acceptor emissions. 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.

Synergy NEO2 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

### Setup

Top filter cube	EX 330 / LUM EM 620 / 665 / LUM
Light source	Xenon flash
Lamp energy	Low (faster)
Delay	150 µs
Data time collection	500 µs
Measurement data point	10
Read height	plate format dependant 8.5mm for 384 wells low volume
Read speed	Normal
Gain	Automatic gain adjustment Autoscale

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



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