

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

RUBYstar is an HTRF® dedicated reader. Its specific optical design enables the simultaneous measurement of both 620 nm cryptate and 665 nm acceptor emissions. A ratio of the two fluorescence intensities* (acceptor/donor) allows the calculation of Delta F (%), i.e. the relative energy transfer rate for each data point.

The basic measurement conditions should be set up in the instrument software according to the following indications:

Setup	
Integration delay (lag time)	50 µs
Integration time	400 µs
Number of flashes	20



**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 BMG LABTECH. The method is covered by the US patent 5,527,684 and its foreign equivalents.*

HTRF® Terbium cryptate donor / Red acceptor readout Setup recommendations for RUBYstar

RUBYstar is an HTRF® dedicated reader. Its specific optical design enables the simultaneous measurement of both 620 nm cryptate and 665 nm acceptor emissions. A ratio of the two fluorescence intensities* (acceptor/donor) allows the calculation of Delta F (%), i.e. the relative energy transfer rate for each data point.

The basic measurement conditions should be set up in the instrument software according to the following indications:

Setup	
Integration delay (lag time)	50 μ s
Integration time	400 μ s
Number of flashes	20

Only BLACK plates must be used with Terbium cryptate!



**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 BMG LABTECH. The method is covered by the US patent 5,527,684 and its foreign equivalents.*