

HTRF setup recommendations for RUBYstar.



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

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 μs
Integration time	400 μs
Number of flashes	20

Only BLACK plates must be used with Terbium cryptate

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