

HTRF[®] Europium cryptate donor / Red acceptor readout Setup recommendations for CLARIOstar

CLARIOstar is equipped with a specific optical device which enables the measurement of both 620 nm cryptate and 665 nm acceptor emissions. The ratio* of the two fluorescence intensities 665/620 (acceptor/donor) allows the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

CLARIOstar readers must be appropriately configured for HTRF[®] readout by setting up the measurement conditions in the software according to the following indications:

Setup	
Excitation filter	330 (10) nm
Emission filters	620 (10) nm 665 (10) nm
Integration delay (lag time)	60 μs
Integration time	400 μs
Number of flashes	200
Optimal z-pos [§]	Volume and plate format dependent
Gain	Select gain adjustment (plate) Required value 90%



**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.*