

## HTRF<sup>®</sup> Terbium cryptate donor / Green acceptor readout Setup recommendations for CLARIOstar

CLARIOstar is equipped with a specific optical device which enables the measurement of both 620 nm cryptate and 520 nm acceptor emissions. The ratio\* of the two fluorescence intensities 520/620 (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<sup>®</sup> readout by setting up the measurement conditions in the software according to the following indications:

Setup	
Excitation filter	330 (80) nm
Emission filters	620 (10) nm 520 (10) nm.
Integration delay (lag time)	60 µs
Integration time	400 µs
Number of flashes	200
Optimal z-pos <sup>§</sup>	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<sup>®</sup> 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.*