

## HTRF<sup>®</sup> readout - Set up recommendations for RUBYstar

RUBYstar is an HTRF<sup>®</sup> dedicated reader. Its specific optical design enables the simultaneous measurement of both the 620 nm cryptate emission and the 665 nm XL665 emission. A ratio of the two fluorescence intensities\* (acceptor/donor) then 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 :

### Set up

---

Integration delay (lag time)	50 $\mu$ s
Integration time	400 $\mu$ s
Number of flashes	20

*\* The fluorescence ratio is a correction method developed by CIS bio international, whose application is limited to the use of HTRF<sup>®</sup> reagents and technology, and for which CIS bio international has granted a license to BMG LABTECH. The method is covered by the US patent 5,527,684 and its foreign equivalents.*