

## HTRF<sup>®</sup> Europium cryptate donor / Red acceptor readout Setup recommendations for SpectraMax iD5<sup>®</sup>

Two sequential measurements should be carried out: at 620 nm for the cryptate emission, and at 665 nm for the specific signal emitted by the acceptor (XL665 or d2). The ratio of the two fluorescence intensities 665/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

SpectraMax iD5 is a hybrid monochromator and filter-based and **only filter-based coupled with "Enhanced TRF" module is compatible**. The measurement conditions should be set up in the SoftMax<sup>®</sup> Pro software according to the following indications:

### Setup

|                                      |  |
|--------------------------------------|--|
| TRF module                           | Must be equipped with Enhanced TRF module (cartridge) in replacing of Standard TRF module  |
| Detection mode                       | Filter module  |
| Excitation filter (bandwidth)        | 340nm (70nm)   |
| Acceptor emission filter (bandwidth) | 665nm (10nm)   |
| Donor emission filter (bandwidth)    | 616nm (10nm)   |
| Number of flashes                    | 50   |
| Excitation time                      | 0.05ms (fixed value)   |
| Measurement delay                    | 0.1ms  |
| Integration time                     | 0.6ms  |
| Read height                          | Volume and plate format dependant.<br>Must be optimized before each new configured measurement using the labware optimization procedure of the software. |



## HTRF® Terbium cryptate donor / Red acceptor readout Setup recommendations for SpectraMax iD5®

Two sequential measurements should be carried out: at 620 nm for the cryptate emission, and at 665 nm for the specific signal emitted by the acceptor (XL665 or d2). The ratio of the two fluorescence intensities 665/620 (acceptor/donor) enables the calculation of Delta F (%) which represents the relative energy transfer rate for each sample.

SpectraMax iD5 is a hybrid monochromator and filter-based and **only filter-based coupled with "Enhanced TRF module is compatible"**. The measurement conditions should be set up in the SoftMax® Pro software according to the following indications:

### Setup

|                                      |  |
|--------------------------------------|--|
| TRF module                           | Must be equipped with Enhanced TRF module (cartridge) in replacing of Standard TRF module  |
| Detection mode                       |  |
|                                      | Filter module  |
| Excitation filter (bandwidth)        |  |
| Acceptor emission filter (bandwidth) | 340nm (70nm)   |
| Donor emission filter (bandwidth)    | 665nm (10nm)   |
|                                      | 616nm (10nm)   |
| Number of flashes                    |  |
| Excitation time                      | 30   |
|                                      | 0.05ms (fixed value)   |
| Measurement delay                    | 0.02ms   |
| Integration time                     | 0.2ms  |
| Read height                          | Volume and plate format dependant.<br>Must be optimized before each new configured measurement using the labware optimization procedure of the software. |

