

# HTRF Europium cryptate donor / Red acceptor readout Setup recommendations for PHERAstar FSX Lamp.



PERAstar FSX is equipped with a specific optical device, which enables the simultaneous measurement of both 620 nm cryptate and 665 nm acceptor emissions. 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.

HTRF™ readout can be achieved by PHERAstar FSX after the installation of the HTRF dedicated optical block which includes the optimized excitation and emission filters, the dichroic mirror and the beam splitter. The measurement conditions should then be set up in the instrument software according to the following indications:

Setup	
Optic module	HTRF 337/620/665 Ref: 906D1
Energy source	Lamp
Integration delay (lag time)	60 μs
Integration time	400 μs
Number of flashes	300
Optimal z-pos <sup>§</sup>	Volume and plate format dependent

<sup>§</sup>The focal height "z" is automatically calculated according to the plate format and the final working volume dispensed in the plate.