

HTRF® Terbium cryptate donor / Red acceptor readout Setup recommendations for Victor Nivo 5S/5F

Victor Nivo is declined under 2 versions (3S/3F for 3 technologies and 5S/5F for 5 technologies) and only 5S/5F model is equipped with TR-FRET module.

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.

Victor Nivo 5S/5F reader must be appropriately configured for HTRF® readout by setting up the measurement conditions in the software according to the following indications:

Setup	
Excitation filter	320/75
Emission filter 1	620/10
Emission filter 2	665/8
Dichroic mirror	D400
Delay time	70µs
Emission time	200µs
Flash energy	Low (10mJ)
Measurement time	500ms
Z -focus (mm)	12 for 384sv or 96LV

This reader only allows high performance HTRF measurement when assays are run in WHITE plates.

