

LanthaScreen® Terbium Assay Setup Guide on the PerkinElmer VICTOR™ Multilabel Plate Reader

NOTE: The PerkinElmer VICTOR™ Multilabel Plate Reader was tested for compatibility with Invitrogen's LanthaScreen® Terbium-based TR-FRET Assay. The following document is intended to demonstrate setup of this instrument. For more detailed information and technical support of Invitrogen assays please call 1-800-955-6288, select option "3", then extension 40266. For more detailed information and technical support of PerkinElmer instruments or software, please call 1-800-762-4000 or by e-mail at productinfo@perkinelmer.com.

A. Recommended Optics

Invitrogen part number	wavelength (nm)	diameter (mm)
Excitation Filter ¹	340/60	15
Emission Filter 1 (PV00325)	495/10	25
Emission Filter 2 (PV00325)	520/25	25

¹ TRF excitation filter pre-installed

B. Instrument Setup

Protocol description

Protocol name LanthaScreen
Protocol number N/A
Name of the plate type NUNC 384
Number of wells in the plate 16 X 24
Height of the plate 14.4 mm
Offset of the wells 9.000 mm, 12.000 mm
Distance between wells 4.500 mm, 4.500 mm
Number of repeats 1
Delay between repeats 0 s
Measurement height 8.00 mm
Protocol notes

Name of the label tb
Label technology Time resolved fluorometry
Emission filter name 520
Emission filter slot B6
Second meas. emission filter name . 495
Second meas. emission filter slot . B7
Emission aperture Normal
Excitation filter D340
Delay 100 µs
Window time 200 µs
Cycle 1000 µs
Second window delay time 0 µs
Second window time 0 µs
Light integrator capacitors 1
Light integrator ref. level 144
Flash energy area High
Flash energy level 210
Flash absorbance measurement No
Beam Normal