LanthaScreen[®] Terbium (Tb) Assay Setup Guide on the Berthold Technologies Tristar² S LB 942 Microplate Reader

The Berthold Technologies Tristar² S LB 942 Microplate Reader was tested for compatibility with LanthaScreen[®] Activity Assay, a TR-FRET assay from Thermo Fisher Scientific, using LanthaScreen[®] Fluorescein-labeled Poly GT (PV3610) and Tb-anti-GST Antibody (PV3550).

The following document is intended to demonstrate setup of this instrument for any Tb-based TR-FRET assay and provide representative data. For more detailed information and technical support of Thermo Fisher Scientific assays please call 1-800-955-6288 ext. 40266. For more detailed information and technical support of Berthold Technologies' instruments or software, please contact Berthold Technologies Bioanalytic at +49 7081-177-0 or www.berthold-bio.com.

A. Recommended Optics

	Wavelength (nm)	Berthold Technologies' Filters	Included in Filter Package	
Excitation	320	320/40 (ld. Nr 52733¹)	ld. Nr. 62772	
Emission 1	520	520/10uv (ld. Nr. 38836 ¹)	ld. Nr. 68492 or 62772	
Emission 2	495	495/10uv (ld. Nr. 39798 ¹)	ld. Nr. 68492	

¹ Although filters 52733, 38836 and 39798 are available separately, they need to be mounted in a special way to be suitable for LanthaScreen[®]. Use only the filters provided in filter packages 68492 or 62772. Do not disassemble, change or move the filters included in the LanthaScreen[®] filter slides, as that could render them unsuitable for LanthaScreen[®]. Do not use filters purchased separately.

Filter Package 62772 includes:

320/40, Id. Nr. 52733
340/26, Id. Nr. 54083*
620/10uv, Id. Nr. 47731*
665/7uv, Id. Nr. 52544*
520/10uv, ld. Nr. 38836

Filter Package 68492 includes

Excitation slide:	340/26, Id. Nr. 54083*
Emission slide:	520/10uv, ld. Nr. 38836
	495/10uv, ld. Nr. 39798

* Not used in this application.

For this assay the excitation slide from filter package 62772 and the emission slide from filter package 68492 are used.

- **Note**: Eu-based TR-FRET and Tb-based TR-FRET use different excitation filters for optimal performance.
- Note: Monochromator based detection is not recommend for TR-FRET assays.

B. Instrument Setup

The following instructions are provided for the MikroWin software. The ICE software is also compatible with LanthaScreen[®], and the same instrument settings can be easily programmed in ICE. Contact Berthold Technologies if you need support to program in ICE the instrument settings detailed here.

- 1. Make sure the plate reader is turned on and then open the MikroWin software on the computer.
- Click on Instrument >> Excitation Filter Slide. Check if the right filters are assigned to the right positions of the filter slides; if they are not, assign each filter to the corresponding position in the filter slide. Please follow the example below:

Excitation slide xD (from filter package 62772) Slot 1: 320/40 (HTRF Eu cryptate); usage: TRFluorescence Slot 2: 340/26 (HTRF Tb cryptate); usage: TRFluorescence

Excitation Filter Slide		\times
Unused Filters 355/40 (Umbelliferone) 405/10 (Absorbance) 492/10 (Absorbance) 620/10 (Absorbance)	All filters can be dragged from the Unused Filters list to an empty position in the slide, or from one position to an empty one or back to the list. Excitation Filter Slide Positions $320/4\Gamma \ 340/2F^2 \ 340/2F^2 \ 5$	
	Excitation filter slide xD 🗸	
Add Remove	Eject Slide OK Cancel	



3. Click on Instrument >> Emission Filter Slide. Check if the right filters are assigned to the right positions of the filter slides; if they are not, assign each filter to the corresponding position in the filter slide (add new filters and enter the settings below if needed). Please follow the example below:

Emission Slide mK (from filter package 68492) Slot 2: 495/10uv; usage: TRFluorescence Slot 3: 520/10uv; usage: TRFluorescence

Emission Filter Slide		×				
Unused Filters 460/25 (Umbelliferone) 480/20 Coelenterazin 530/25 eYFP 400/70 Deep Blue C 515/20 GFP2 565/70 (Alpha Screen) 510/60 ChromaGloGreen 610/lp ChromaGloRed	Filters can be dragged from the Unused Filters list an empty position in the wheel, or from one sition to an empty one or back to the list. ission Filter Slide Positions					
Add Remove	Eject Slide OK Cancel					

4. If you already have a pre-existing template for LanthaScreen[®], open it and use this document to review your settings; if you don't have yet any suitable template, click on Settings in the menu bar at the top portion of the window to start creating a new template.



5. A new window will open. Select the Plate type corresponding to the plate you are using and highlight the wells you most commonly will measure. If unsure about what plate type to select, contact Berthold Technologies for assistance.

Options	×
Samples Measurement	
Samples Measurement Plate parameter Plate type: Plate type: Berthold 96 - No:23300/23302 Plate Mode: Costar 24 3524 Plate Speed: NUNC Multidish 48 Measurement Order: BD Falcon 24-Multiwell Insert BD Falcon 24-Multidish 12 D Falcon 24-Multidish 12 NUNC Multidish 12 Terasaki 60 Terasaki 72 Terasaki 72	
A Petridish 35mm NUNC 384 Costar 96 clearbottom Costar 96 clear Greiner 96 Costar 96 HalfArea Optiplate 384 BD Falcon HTS FluoroBlok 24-Multiwell Berthold 96 clearbottom - No:24910/38840 Berthold 24 clearbottom - No:41081/41082	
F Petnaisn 6.3mm Berthold 96 sterile - No:51838/51839 G H OK Abbrechen	ehmen



6. Click on the Measurement tab and look for the TRF operation.

🜖 MikroWin - [Templa	te - Ma	trices - Calculatio	n]									_	
File Edit View Inst	rument	Installation F	Help	tings	Plate ID		Status Inf	ormation	Start	Print	Export		
Reader	iate M	Calculation Form	ula of Position A0	1: MEA		Add	Formula						
Instrument	#	1	2	3	4	5	6	7	8	9	10	11	12
Plate Layout Plate Layout Grouping	A	MEA	MEA	MEA	MEA	Options Sample	s Measurement				×	MEA	MEA
Overlay Matrices Controls	в	MEA	MEA	MEA	MEA	- Mea	surement Sequence:	Operations:	Ope	ration		MEA	MEA
Classifications Curvefitting Kinetics	с	MEA	MEA	MEA	MEA	5	 Dispense Lumin. Label K Lumin. Kinetic S Lumin. Scannir R Lumin. Repeat 	ng ed				MEA	MEA
Scanning Definitions Statistics	D	MEA	MEA	MEA	MEA		K Fluor. Label K Fluor. Kinetic S Fluor. Scanning R Fluor. Repeate S Fluor. Spectral	g d Scar				MEA	MEA
Compilation	E	MEA	MEA	MEA	MEA	Ope	ration Mode	> Bepeats: 1				MEA	MEA
Export Configuration	F	MEA	MEA	MEA	MEA	Tem	perature Control Temperature	25	°Celsius			MEA	MEA
	G	MEA	MEA	MEA	MEA		Keep Plate after Mea	s.	ОК	Abbrechen	Übernehmen	MEA	MEA
	н	MEA	MEA	MEA	MEA	MEA	MEA	MEA	MEA	MEA	MEA	MEA	MEA
	н	Sample_	ID 2 Reade	er_Values	Results								4 Þ H
		_	_		Measure	ment not perfo	ormed Vali	d Assay		Untitled.dat		default.par	

- 7. Double click on TRF to insert a TRF measurement operation. A new window will appear. If desired, enter a Name for the measurement operation. Configure the settings as shown in the screenshot below:
 - Enter Counting Time: 1.00
 - Select Aperture: 3 (Rd 2)
 - Select Excitation Filter: 320/40 (HTRF Eu cryptate)*
 - Select Excitation Optic: 3 Wide Filter 0.45mm
 - Select Emission Filter: 495/10uv*
 - Enter Timing settings: Cycle Time 5000, Delay Time 100, Reading Time 300
 - Check Second Measurement
 - Select Excitation Filter: 320/40 (HTRF Eu cryptate)*
 - Select Emission Filter: 520/10uv*

When finished, click OK.

* The name of the filters in the software sometimes does not match the LanthaScreen[®] naming conventions, and sometimes filters named as "Tb cryptate" are mentioned in a Eu assay, or the other way around. This is not an error; filter naming was designed for HTRF[®] assays, but for LanthaScreen[®] different filter combinations are sometimes chosen for the best performance.

FRF Label				×
Name:	LanthaScreen Tb		1	ОК
Use Delay before Re	ading (0,1 - 600 s)	0.0	[Cancel
Counting Time:	1.00	(0.05 - 600 s)		
Counter:	Photon	O Current		
Sensitivity:	O Low (500 V)	Medium (700 V)	O High (1000) V)
	O Manual (500 V	. 1000 V)	700	
Lamp Energy:	100		1 1 1 1	
Use	Filters	O Monochromator		
Aperture:	3 - Rd 2		~	•
Excitation Filter:	320/40 (HTRF Eu c	ryptate) - Slot xD1	~	•
Excitation Optic:	3 - Wide Filter 0.45r	nm	~	•
Emission Filter:	495/10uv	- Slot mK2	~	•
Cycle Time: Delay Time: Reading Time: Flashes per well::	5000 100 300 200	(2000 - 10000 μs) (0 - 4560 μs) (20 - 4760 μs)		
Second Measuremen	t			
Excitation Filter:	320/40 (HTRF Eu c	ryptate) - Slot xD1	~	•
Emission Filter:	520/10uv	- Slot mK3	~	•
Operation Mode:	O By plate	By well		



8. To save the template, click on File in the main menu, then Template and Save as. Browse to the desired folder, enter the desired filename and click OK.

🍨 Speichern unter									×
← → • ↑ <mark> </mark>	« Mik	kroWin 2010 → Pa	araTriStar2S → TR-	-Fluorescence		√ Č	"TR-Fluorescence	" durchsuch	٩
Organisieren 🔻	Neuer	Ordner							?
📌 Schnellzugriff 🛄 Desktop	А	Name	^	Es wurden	Änderungsdatum keine Suchergebnisse	Typ gefunden.	Größe		
➡ Downloads	A A A								
Lantha LanthaScreer	n								
Screenshots									
💣 Netzwerk									
Datei <u>n</u> ame: Datei <u>t</u> yp:	Lantha Mikro\	a <mark>Screen.par</mark> Win Para File (*.pai	r)						~
∧ Ordner ausblende	en						<u>S</u> peichern	Abbreche	en

- 9. To start the measurement, enter the desired Plate ID to identify the measurement. If you want to edit the wells to be measured, click on Settings and select the desired wells (see point 3). When you are ready, click Start. The plate tray will open; insert the plate and click OK to start the measurement.
- 10. When the measurement has finished, click Export to export the data for further calculation, if necessary. Example raw data values are displayed below.

Page 8 of 14

 ThermoFisher
 Microplate Reader Documentat

 S C I E N T I F I C
 Sature Cuide on the Berthold J

Setup Guide on the Berthold Technologies Tristar² S LB 942 Microplate Reader

Donor signal (495 nm)				
[Acceptor]	5000 nM	2500 nM	1250 nM	625 nM	200 nM
Α	72080	114730	138331	154378	166714
В	84620	110952	167138	164687	172279
С	84353	119646	163230	158120	179057
D	93596	119719	133115	178583	181955
E	81901	116300	159747	170857	175962
F	78187	121427	151868	173419	176770
G	90046	109680	153657	173173	183518
н	75351	125453	155489	171273	163278
I	76437	118729	149655	167632	176547
J	72668	109586	140843	165270	177934
Acceptor sign	al (520 nm)				
[Acceptor]	5000 nM	2500 nM	1250 nM	625 nM	200 nM
Α	59800	47679	31795	20861	14986
В	58916	44836	35194	21540	14895
С	58518	45205	33611	20813	16441
D	66299	41950	29785	22536	15696
E	59961	40919	34110	22562	15093
F	54247	46138	32433	22146	16135
G	59240	40088	32845	21755	15922
н	58345	47410	31741	21392	14084
I	58952	46061	31982	21586	14125
J	57458	45350	31469	21430	14488

- 11. Plots of ratios corresponding to these raw data are displayed below.
- A. Ratio Data

B. Normalized Data



12. These values were obtained using the procedure detailed in the next section. Additional representative data from the Berthold Technologies Tristar² S are available at the end of the section.

Version No.: March 2022

Page 9 of 14

Setup Guide on the Berthold Technologies Tristar² S LB 942 Microplate Reader

Test Your Plate Reader Set-up Before Using LanthaScreen® Tb Assays

<u>Purpose</u>

This LanthaScreen[®] Tb Microplate Reader Test provides a method for verifying that a fluorescent plate reader is able to detect a change in time-resolved fluorescence energy transfer (TR-FRET) signal, confirming proper instrument set-up and a suitable response. The method is independent of any biological reaction or equilibrium and uses reagents that are on-hand for the LanthaScreen[®] assay.

<u>At a Glance</u>

Step 1: This document can be found at <u>www.thermofisher.com/instrumentsetup</u>.

Step 2: Prepare individual dilutions of the TR-FRET acceptor (fluorescein-labeled substrate, e.g. PV3610 or other).

2X = 10,000 nM, 5,000 nM, 2,500 nM, 1,250 nM and 400 nM.

Note: This application is NOT suited for LanthaScreen[®] GFP-tagged physiological substrates or Nuclear Receptor Fluormones. Instead, use Fluorescein-poly GT (e.g. PV3610) as the acceptor.

Note: To avoid propagating dilution errors, we do NOT recommend using serial dilutions. See page 10.

Step 3: Prepare a dilution of the TR-FRET donor (Tb-Antibody, e.g. PV3550 or other)

$2X = 125 \text{ nM Tb}^{3+}$ chelate.

Note: Concentration is based on the molarity of the Tb chelate (found on the Certificate of Analysis), NOT the molarity of the antibody, to account for normal variation in antibody labeling. See pages 11 - 12 for calculations and method.

Step 4: Prepare plate and read.

Step 5: Contact Technical Support with your results. E-mail us directly at <u>drugdiscoverytech@thermofisher.com</u> or in the US call 1-800-955-6288 ext. 40266. We will determine Z'-factors by comparing each concentration of acceptor to the 200 nM acceptor data. Example results and data analysis are available on page 14.

Introduction

This LanthaScreen[®] Tb Microplate Reader Test uses diffusion-enhanced TR-FRET to generate a detectable TR-FRET signal. At high donor or acceptor concentrations, donor and acceptor diffuse to within a suitable distance from one another to allow TR-FRET to take place, resulting in a signal. The response in diffusion-enhanced TR-FRET is easy to control because it is directly proportional to the concentrations of donor and acceptor in solution and is not related to a binding event.

In this method, acceptor concentration varies while the donor concentration remains fixed. As the concentration of acceptor increases, the diffusion-enhanced TR-FRET signal increases. The signal from the acceptor concentrations are compared to the signal from the lowest acceptor concentration to simulate assay windows from high to low to help you assess whether your instrument is properly set-up and capable of detecting TR-FRET signals in the LanthaScreen[®] Assays.

We designed the LanthaScreen[®] Tb technical note to use components and reagents that are generally used in most LanthaScreen[®] Assays. Please note that the LanthaScreen[®] GFP-tagged physiological substrates and Nuclear Receptor Fluormone tracers are not suited for this method. Instead, use Fluorescein-poly GT, PV3610, as the acceptor.



Materials Required

Component	Storage	Part Number	Example Reagents
LanthaScreen [®] Tb-labeled antibody (donor)	-20°C	Various	PV3550
LanthaScreen [®] fluorescein-labeled substrate (acceptor)	-20°C	Various	PV3610
TR-FRET Dilution Buffer or any Nuclear Receptor Co- regulator Buffer	Various	PV3574 or Various	PV3574

96-well polypropylene microplate or 1.5 mL microcentrifuge tubes

384-well plate (typically a white, low-volume Corning 4513 or black, low-volume Corning 4514)

Plate seals

Suitable single and multichannel pipettors

Plate reader capable of reading TR-FRET

Handling

To reread the plate on another day, seal and store the plate at room temperature for up to 5 days. To reread the plate, centrifuge the plate at 300 xg for 1 minute, remove seal and read.

Important: Prior to use, centrifuge the antibody at approximately 10,000 xg for 5 minutes, and carefully pipette the volume needed for the assay from the supernatant. This centrifugation pellets aggregates present that can interfere with the signal.

Procedure

Step 1: Set up your instrument using the information in this document.

Step 2: Prepare the Acceptor (LanthaScreen[®] fluorescein substrate or peptide)

Acceptor concentrations (2X) are individually prepared from a 30 μ M stock to prevent propagation of error that can occur with serial dilutions. We suggest preparing 10 replicates for calculation of a Z'-factor. To accommodate replicates that use 10 μ L per well, prepare 120 μ L of each concentration. Prepare each concentration in micro-centrifuge tubes or a 96-well polypropylene plate and then transfer it to a 384-well plate.

Note: This application is NOT suited for LanthaScreen[®] GFP-tagged physiological substrates or Nuclear Receptor Fluormones. Instead, use Fluorescein-poly GT (Cat no. PV3610) or other Fluorescein-peptides as the acceptor.

Fluorescein Substrate/Peptide	Cat #	Concentration as Sold	Dilution to prepare a 30 µM solution
Fluorescein-Poly GT	PV3610	30 µM	No dilution needed
Fluorescein-Poly GAT	PV3611	30 µM	No dilution needed
Fluorescein peptides for kinases	Various	Various	Add 6 μ L of 1 mg/mL peptide stock to 94 μ L of TR-FRET Dilution Buffer (peptide at 1 mg/mL with a MW ~ 2kDa = ~500 μ M)
Fluorescein co-regulator peptides for NRs	Various	100 μM	Add 30 μ L of 100 μ M peptide stock to 70 μ L of TR-FRET Co-regulator Buffer. Do not add DTT.

1. Prepare 30 µM acceptor stock solution:

2. Prepare 120 µL of each 2X acceptor concentration from the 30 µM stock:

96-well plate or tubes	A1	B1	C1	D1	E1
2X Acceptor Concentration	10,000 nM	5,000 nM	2,500 nM	1,250 nM	400 nM
Final 1X Acceptor Concentration	5,000 nM	2,500 nM	1,250 nM	625 nM	200 nM
Volume TR-FRET Dilution Buffer or NR Coregulator Buffer	80.0 μL	100.0 μL	110.0 μL	115 μL	118.4 μL
Volume 30 µM Acceptor (prepared above)	40.0 µL	20.0 µL	10.0 µL	5.0 μL	1.6 μL

Step 3: Prepare the Donor (Tb-chelate labeled antibody)

Prepare a 2X stock of Tb-chelate at 125 nM that will result in a final assay concentration of 62.5 nM. This method relies on the concentration of Tb-chelate, NOT the concentration of antibody. The lot-to-lot variation in the number of Tb-chelates covalently bound to antibody can be accounted for by referring to the Tb-chelate-to-antibody ratio listed on the lot-specific Certificate of Analysis for your antibody. Multiply this ratio by the antibody concentration to calculate the Tb-chelate concentration.

Example chelate concentrations

Antibody Concentration	Antibody Molarity	Chelate: Antibody Ratio	Chelate Concentration
0.5 mg/mL	3.3 μM	11	$36.3 \ \mu M = 36,300 \ nM$
0.25 mg/mL	1.7 μM	8	$13.6 \mu\text{M} = 13,600 \text{nM}$

Example Calculation: Prepare 1,000 µL of Tb-chelate:

Tb-antibody = $0.5 \text{ mg/mL} (3.3 \mu \text{M})$ with a chelate:antibody ratio of 11

Chelate: Stock = $3.3 \ \mu M \ x \ 11 = 36.3 \ \mu M = 36,300 \ nM.$

 $1X = 62.5 \text{ nM}; \quad 2X = 125 \text{ nM}$

Т	he	9	'n	nc	F	Ĩ.	sł	1	er	
S	С	Т	Е	Ν	т	Т	F	Т	С	

LanthaScreen® Tb CompatibleVersion No.:Page 12 of 14Microplate Reader DocumentationMarch 2022

Setup Guide on the Berthold Technologies Tristar² S LB 942 Microplate Reader

Formula	V ₁	X	C ₁	=	V_2	X	C ₂
			[Stock]				[2X]
Th-Chelate	V ₁	Х	36,300 nM	=	1,000 µL	Х	125 nM
10-Chelate	V ₁ = 3	.4 μL					

Add 3.4 µL of 36,300 nM stock to 996.6 µL TR-FRET dilution buffer or NR coregulator buffer.

Step 4: Add Reagents to the 384-well plate and read

1. Donor

Transfer 10 μ L of 2X Tb-chelate to rows A through J and columns 1 through 5 of the 384-well assay plate. Since you need only a single concentration, you can transfer this solution with a multichannel pipettor from a basin to all 50 wells. We recommend preparing the 1 mL solution in a 1.5 mL micro-centrifuge tube before transferring into the basin.

2. Acceptor

Note: To eliminate carryover, we recommend changing pipette tips for each concentration of acceptor. *Note:* After adding 2X acceptor, mix the reagents by pipetting up and down.

Transfer 10 μ L of the indicated concentration of 2X acceptor to the rows A-J of the corresponding column of the 384- well plate.

2X Acceptor	Column
10,000 nM	1
5,000 nM	2
2,500 nM	3
1,250 nM	4
400 nM	5

3. Read the plate

This step does not require any equilibration time.

Step 5: Contact Technical Support

Send us your results by e-mailing us directly at <u>drugdiscoverytech@thermofisher.com</u> or in the US call 1-800-955-6288 ext. 40266.

Version No.: March 2022

Page 13 of 14

Setup Guide on the Berthold Technologies Tristar² S LB 942 Microplate Reader

B. Normalized Data

We will help you evaluate your results by performing the following data analysis:

- 1. Obtain the emission ratios by dividing the acceptor signal (520 nm) by the donor signal (495 nm, exact wavelength varies with instrument) for each well.
- Calculate the average ratio for each column (1 through 5). These values can be plotted against the final 1X concentrations (5,000 nM, 2,500 nM, 1,250 nM, 625 nM, and 200 nM) of acceptor (see graph A). Dilution curves form diffusion-enhanced TR-FRET do not plateau and, therefore, do not fit the normal sigmoidal shape produced by binding curves.
- 3. Using the data from column 5 (200 nM acceptor) as the bottom of the "assay window," divide the average ratios from the other columns b the average ratio from column 5 to obtain a range of simulated "assay window" sizes. See the example data below. This "normalized" data can be plotted against the acceptor concentration as shown below in graph B.
- 4. Calculate the Z'-factor for each "assay window." Very general guidance is that you should observe a satisfactory Z'-factor (>0.5) for at least the "small window" that compares columns 3 to 5 (1,250 nM to 200 nM). In our hands and on certain instruments, the data in columns 4 and 5 produces suitable Z'-factors (>0.5) with a simulated assay window of less than 2.



Thermo Físher

SCIENTIFIC

Columns Compared	Description
1 to 5	Largest window
2 to 5	Intermediate window
3 to 5	Small window
4 to 5	Smallest window, less than 2-fold

Version No.:

March 2022

[Acceptor]	5,000 nM	2,500 nM	1,250 nM	625 nM	200 nM
Row A	0.830	0.416	0.230	0.135	0.090
Row B	0.696	0.404	0.211	0.131	0.086
Row C	0.694	0.378	0.206	0.132	0.092
Row D	0.708	0.350	0.224	0.126	0.086
Row E	0.732	0.352	0.214	0.132	0.086
Row F	0.694	0.380	0.214	0.128	0.091
Row G	0.658	0.365	0.214	0.126	0.087
Row H	0.774	0.378	0.204	0.125	0.086
Row I	0.771	0.388	0.214	0.129	0.080
Row J	0.791	0.414	0.223	0.130	0.081

Example data: Ratiometric data obtained on a Berthold Technologies Tristar² LB 942 microplate reader.

Data Analysis:

[Acceptor]	5,000 nM	2,500 nM	1,250 nM	625 nM	200 nM
Average Ratio	0.735	0.382	0.215	0.129	0.087
St dev	0.054	0.023	0.008	0.003	0.004
% CV	7.4	6.1	3.8	2.5	4.4
Assay Window	8.49	4.42	2.49	1.49	Reference
Z'-factor	0.73	0.73	0.72	0.50	

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.