# LanthaScreen<sup>™</sup> TR-FRET Farnesoid X Receptor Coactivator Assay

Catalog no. PV4833

Shipping Condition: Dry Ice

**Initial Storage: Varies** 

Protocol part no. PV4833.pps

Rev. date: 10-Oct-2011

# **TABLE OF CONTENTS**

invitrogen<sup>®</sup>

<ul> <li>2. Materials Required but Not Supplied</li></ul>	1.	. Kit Contents1				
3.       Introduction         3.1       Principle of FRET and TR-FRET         3.2       Assay Overview         4.       Instrument Settings         5.       Guidelines and Recommendations         5.1       Reagent Handling         5.2       Ligand Dilutions         5.3       Solvent Tolerance         5.4       Note on Reagent Order of Addition         5.5       Incubation Conditions         6.       Agonist Assay         6.1       Agonist Assay         6.2       Plate Layout         6.3       Agonist Assay         6.4       Agonist Assay         6.5       Agonist Assay         6.6       Agonist Assay         7.       Assay Pharmacology	2.		2			
3.1       Principle of FRET and TR-FRET			2			
3.2       Assay Overview         4.       Instrument Settings         5.       Guidelines and Recommendations         5.1       Reagent Handling         5.2       Ligand Dilutions         5.3       Solvent Tolerance         5.4       Note on Reagent Order of Addition         5.5       Incubation Conditions         6.       Agonist Assay         6.1       Agonist Assay         6.2       Plate Layout.         6.3       Agonist Assay         6.4       Agonist Assay         6.5       Agonist Assay         7.       Assay Pharmacology			2			
<ul> <li>5. Guidelines and Recommendations</li> <li>5.1 Reagent Handling</li> <li>5.2 Ligand Dilutions</li> <li>5.3 Solvent Tolerance</li> <li>5.4 Note on Reagent Order of Addition</li> <li>5.5 Incubation Conditions</li> </ul> 6. Agonist Assay 6.1 Agonist Assay Procedure <ul> <li>6.2 Plate Layout.</li> <li>6.3 Agonist Assay</li> <li>6.4 Agonist Assay</li> <li>7. Assay Pharmacology</li> </ul>	3.2		2			
<ul> <li>5. Guidelines and Recommendations</li> <li>5.1 Reagent Handling</li> <li>5.2 Ligand Dilutions</li> <li>5.3 Solvent Tolerance</li> <li>5.4 Note on Reagent Order of Addition</li> <li>5.5 Incubation Conditions</li> </ul> 6. Agonist Assay 6.1 Agonist Assay Procedure <ul> <li>6.2 Plate Layout.</li> <li>6.3 Agonist Assay</li> <li>6.4 Agonist Assay</li> <li>7. Assay Pharmacology</li> </ul>	4.	Instrument Settings	3			
5.1       Reagent Handling         5.2       Ligand Dilutions         5.3       Solvent Tolerance         5.4       Note on Reagent Order of Addition         5.5       Incubation Conditions         6.       Agonist Assay         6.1       Agonist Assay         6.2       Plate Layout         6.3       Agonist Assay         6.4       Agonist Assay         6.5       Agonist Assay         7.       Assay Pharmacology	5.	-	4			
5.2       Ligand Dilutions         5.3       Solvent Tolerance         5.4       Note on Reagent Order of Addition         5.5       Incubation Conditions         6.       Agonist Assay         6.1       Agonist Assay         6.2       Plate Layout.         6.3       Agonist Assay         6.4       Agonist Assay         6.5       Agonist Assay         7.       Assay Pharmacology	•••		4			
5.4       Note on Reagent Order of Addition         5.5       Incubation Conditions         6.       Agonist Assay         6.1       Agonist Assay         6.2       Plate Layout.         6.3       Agonist Assay         6.4       Agonist Assay         6.5       Agonist Assay         7.       Assay Pharmacology	5.2		4			
5.5       Incubation Conditions         6.       Agonist Assay         6.1       Agonist Assay—Procedure.         6.2       Plate Layout.         6.3       Agonist Assay—Data Analysis         6.4       Agonist Assay—Reagent Volumes and Controls         6.5       Agonist Assay—Representative Data         7.       Assay Pharmacology	5.3					
<ul> <li>6. Agonist Assay</li></ul>						
<ul> <li>6.1 Agonist Assay—Procedure</li></ul>	5.5	Incubation Conditions	5			
<ul> <li>6.1 Agonist Assay—Procedure</li></ul>	6.	Agonist Assay	ò			
<ul> <li>6.3 Agonist Assay—Data Analysis</li> <li>6.4 Agonist Assay—Reagent Volumes and Controls</li> <li>6.5 Agonist Assay—Representative Data</li> <li>7. Assay Pharmacology</li> </ul>	6.1	Agonist Assay—Procedure				
<ul> <li>6.4 Agonist Assay—Reagent Volumes and Controls</li> <li>6.5 Agonist Assay—Representative Data</li> <li>7. Assay Pharmacology</li> </ul>	6.2	Plate Layout	7			
<ul> <li>6.5 Agonist Assay—Representative Data</li> <li>7. Assay Pharmacology</li> </ul>	6.3	Agonist Assay—Data Analysis	7			
7. Assay Pharmacology						
	6.5	Agonist Assay—Representative Data	3			
8. References	7.	Assay Pharmacology	)			
	8.	References	9			
9. Notice to Purchaser1	9.	Notice to Purchaser10	)			

#### **KIT CONTENTS** 1.

The LanthaScreen<sup>™</sup> TR-FRET Farnesoid X Receptor Coactivator Assay contains the following:

Component	Composition	Amount	Storage Temp.	Individual Catalog no.
Fluorescein-SRC2-2 coactivator peptide (Voegel <i>et al.</i> , 1996)	100 μM in 50 mM HEPES, pH 7.5 Sequence: LKEKHKILHRLLQDSSSPV	100 µl	-20°C	PV4586
FXR LBD, GST	FXR ligand-binding domain in a buffer (pH 8.0) containing protein stabilizing reagents and glycerol.	10 µg	-80°C	PV4834
LanthaScreen <sup>™</sup> Tb-anti- GST antibody	Terbium labeled anti-GST antibody in HEPES buffered saline (137 mM NaCl, 2.7 mM KCl, 10 mM HEPES pH 7.5). See Certificate of Analysis for lot specific concentration	25 μg	-20°C	PV3550
Coregulator Buffer G	Proprietary buffer (pH 7.4)	25 ml	4°C	PV4553
DTT, 1 M	In water	1 ml	-20°C or -80°C	P2325

For Technical Support on this and other Drug Discovery Products, dial 760-603-7200, option 3, extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to <u>www.invitrogen.com/lanthascreen</u>

# 2. MATERIALS REQUIRED BUT NOT SUPPLIED

The following materials are required but not supplied in the kit:

- A fluorescence plate reader with excitation at 340 nm and with the appropriate filter sets installed for detecting the fluorescent emission signals of terbium at 495 nm and fluorescein at 520 nm (see Section 4).
- Pipetting devices for 1–1000-µl volumes, suitable repeater pipettors, or multi-channel pipettors.
- Black, 384-well assay plates. We recommend black Corning<sup>®</sup> 384-well, low-volume, round-bottom (non-binding surface) assay plates; Corning #3677. Other plate types may give satisfactory results as well.
- 96-well polypropylene plate that can accommodate a 400-μl volume per well. We recommend Nalgene Nunc #249944.
- A known FXR agonist, such as GW4064, to serve as a positive control. We recommend Tocris catalog no. 2473.
- DMSO

# 3. INTRODUCTION

Invitrogen's LanthaScreen<sup>™</sup> Farnesoid X receptor TR-FRET Coactivator Assay provides a sensitive and robust method for high-throughput screening of potential FXR ligands as agonists of coactivator recruitment. The kit uses a terbium-labeled anti-GST antibody, a fluorescein-labeled coactivator peptide, and an FXR ligand-binding domain that is tagged with glutathione-S-transferase (GST) in a homogenous mix-and-read assay format. This kit contains enough reagents for 800 assays of 20 µl each.

# 3.1 Principle of FRET and TR-FRET

For screening libraries of compounds, time-resolved FRET (TR-FRET) is a recognized method for overcoming interference from compound autofluorescence or light scatter from precipitated compounds. The premise of a TR-FRET assay is the same as that of a standard FRET assay: when a suitable pair of fluorophores is brought within close proximity of one another, excitation of the first fluorophore (the donor) can result in energy transfer to the second fluorophore (the acceptor). This energy transfer is detected by an increase in the fluorescence emission of the acceptor and a decrease in the fluorescence emission of the donor. In HTS assays, FRET is often expressed as a ratio of the intensities of the acceptor and donor fluorophores. The ratiometric nature of such a value corrects for differences in assay volumes between wells and corrects for quenching effects due to colored compounds.

In contrast to standard FRET assays, TR-FRET assays use a long-lifetime lanthanide chelate as the donor species. Lanthanide chelates are unique in that their excited-state lifetime (the average time that the molecule spends in the excited state after accepting a photon) can be on the order of a millisecond or longer. This is in sharp contrast to the lifetime of common fluorophores used in standard FRET assays, which are typically in the nanosecond range. Because interference from autofluorescent compounds or scattered light is also on the nanosecond timescale, these factors can negatively impact standard FRET assays. To overcome these interferences, TR-FRET assays are performed by measuring FRET after a suitable delay, typically 50 to 100 microseconds after excitation by a flashlamp excitation source in a microtiter plate reader. This delay not only overcomes interference from background fluorescence or light scatter, but also avoids interference from direct excitation due to the non-instantaneous nature of the flashlamp excitation source.

The most common lanthanides used in TR-FRET assays for HTS are terbium and europium. Terbium offers unique advantages over europium when used as the donor species in a TR-FRET assay. In contrast to europium-based systems that employ APC as the acceptor, terbium-based TR-FRET assays can use common fluorophores such as fluorescein as the acceptor. Because it is straightforward (and inexpensive) to label a molecule such as a peptide with fluorescein, directly labeled molecules may be used in terbium-based TR-FRET assays, rather than biotinylated molecules that must then be indirectly labeled via streptavidin-mediated recruitment of APC. The use of directly labeled molecules in a terbium-based TR-FRET assay reduces costs, improves kinetics, avoids problems due to steric interactions involving large APC conjugates, and simplifies assay development since there are fewer independent variables requiring optimization in a directly labeled system.

### 3.2 Assay Overview

Binding of agonist to the nuclear receptor (Figure 1) causes a conformational change around helix 12 in the ligand binding domain, resulting in higher affinity for the coactivator peptide. When the terbium label on the anti-GST antibody is excited at 340 nm, energy is transferred to the fluorescein label on the coactivator peptide and detected as emission at 520 nm.

When running the LanthaScreen<sup>TM</sup> TR-FRET FXR Coactivator Assay, FXR -LBD is added to ligand test compounds followed by addition of a mixture of the fluorescein-coactivator peptide and terbium anti-GST antibody. After an incubation period at room temperature, the TR-FRET ratio of 520:495 is calculated and can be used to determine the EC<sub>50</sub> from a dose response curve of the compound. Based on the

For Technical Support on this and other Drug Discovery Products, dial 760-603-7200, option 3, extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to www.invitrogen.com/lanthascreen biology of the FXR-coactivator peptide interaction, this ligand  $EC_{50}$  is a composite value representing the amount of ligand required to bind to receptor, effect a conformational change, and recruit coactivator peptide (see Figure 1).



**Figure 1. Principle of the nuclear receptor (NR) agonist dependent coactivator peptide recruitment assay:** Tb-anti-GST antibody indirectly labels the nuclear receptor by binding to the GST tag. Binding of the agonist to the NR causes a conformational change that results in an increase in the affinity of the NR for a coactivator peptide. The close proximity of the fluorescently labeled coactivator peptide to the terbium-labeled antibody causes an increase in the TR-FRET signal.

# 4. INSTRUMENT SETTINGS

General Settings:		
Excitation	Excitation 340 nm filter (30 nm bandwidth)	
Emission 520 nm filter (25 nm bandwidth)		
Emission	490 or 495 nm filter (10 nm bandwidth)	
Delay Time	100 μs	
Integration Time	200 µs	

The excitation and emission spectra of terbium and fluorescein are shown in Figure 2. As with other TR-FRET systems, the terbium donor is excited using a 340-nm excitation filter with a 30-nm bandwidth. However, the exact specifications of the excitation filter are not critical, and filters with similar specifications will work well. In general, excitation filters that work with europium-based TR-FRET systems will perform well with the LanthaScreen<sup>TM</sup> terbium chelates.

### **Spectra of Terbium and Fluorescein**



Figure 2: Excitation and Emission spectra of fluorescein and terbium.

For Technical Support for this or other Invitrogen Drug Discovery Solutions Products, dial 760-603-7200 extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to www.invitrogen.com/lanthascreen As shown in the figure, the terbium emission spectrum is characterized by four sharp emission peaks, with silent regions between each peak. The first terbium emission peak (centered between approximately 485 and 505 nm) overlaps with the maximum excitation peak of fluorescein. Energy transfer to fluorescein is then measured in the silent region between the first two terbium emission peaks. Because it is important to measure energy transfer to fluorescein without interference from terbium, a filter centered at 520 nm with a 25 nm bandwidth is used for this purpose. The specifications of this filter are more critical than those of the excitation filter. In general, standard "fluorescein" filters may not be used, because such filters also pass light associated with the terbium spectra as well. The emission of fluorescein due to FRET is referenced (or "ratioed") to the emission of the first terbium peak, using a filter that isolates this peak. This is typically accomplished with a filter centered at 490 or 495 nm, with a 10 nm bandwidth. In general, a 490 nm filter will reduce the amount of fluorescein emission that "bleeds through" into this measurement, although instrument dichroic mirror choices (such as those on the Tecan Ultra instrument) may necessitate the use of a 495 nm filter. The effect on the quality of the resulting measurements is minimal in either case. Filters suitable for LanthaScreen<sup>TM</sup> assays are available from Chroma (www.chroma.com) as filter set PV001, or from other vendors. A LanthaScreen for the BMG PheraStar is available direct from BMG Instruments.

Aside from filter choices, instrument settings are similar to the settings used with europium-based technologies. In general, the guidelines provided by the instrument manufacturer can be used as a starting point for optimization. A delay time of 100  $\mu$ s followed by a 200- $\mu$ s integration time is typical for a LanthaScreen<sup>TM</sup> assay. The number of flashes or measurements per well is highly instrument dependent and should be set as advised by your instrument manufacturer. In general, LanthaScreen<sup>TM</sup> assays can be run on any filter-based instrument capable of time-resolved FRET, such as the Tecan Ultra, BMGLabTech PHERAStar, Molecular Devices Analyst, or PerkinElmer Envision. LanthaScreen<sup>TM</sup> assays have also been performed successfully on the Tecan Safire<sup>2</sup> and Molecular Devices M5 monochromator-based instrument. Contact Invitrogen Technical Services for instrument-specific setup guidelines.

# 5. GUIDELINES AND RECOMMENDATIONS

# 5.1 Reagent Handling

#### FXR-LBD

Store FXR-LBD at -80°C. Thaw on ice before use and perform all dilutions while on ice. Never vortex the FXR-LBD stock or dilutions.

#### Coregulator Buffer G

Store Coregulator buffer G at 4°C. Mix well before first use, as the buffer is viscous.

# LanthaScreen<sup>™</sup> Tb anti-GST Antibody

Store Tb anti-GST antibody at -20°C and thaw before use.

### Fluorescein-SRC2-2 Coactivator Peptide

Store fluorescein-SRC2-2 coactivator peptide at -20°C and thaw before use.

# 5.2 Ligand Dilutions

The procedure described below involves diluting the ligand to 100X in DMSO followed by transfer into complete assay buffer, resulting in a 2X ligand, 2% DMSO dilution. This may be done in a DMSO-tolerant assay plate or polypropylene tubes. (Dilution in DMSO increases solubility of ligands and reduces sticking to plastics.) After addition of all reagents to the assay, the final concentration will be 1X ligand and 1% DMSO.

*Note:* Handling of some ligands can be problematic due to their tendency to stick to various plastics. These ligands may show varying degrees of stickiness, causing differences in the actual concentration of the ligand, depending on the type of pipette tips and plates used in the set up of the assay. Therefore, you may observe different  $EC_{50}$  values than reported here.

### 5.3 Solvent Tolerance

The assay was validated in the presence of 1% DMSO. However, the assay has been performed with up to 8% DMSO, 8% ethanol, and 8% methanol (in addition to the 1% DMSO present from the ligand dilution) with good results.

## 5.4 Note on Reagent Order of Addition

The assay was validated using three additions per well in which FXR-LBD was added to agonist dilutions, followed by the addition of premixed fluorescein-SRC2-2 coactivator peptide and Tb anti-GST antibody (agonist, receptor, peptide/antibody). A pre-mixture of FXR-LBD, fluorescein-SRC2-2 coactivator peptide, and Tb anti-GST antibody may also be added to the agonist dilutions for a total of 2 additions per well, although the assay was not fully validated in this manner. However, it is important to consider the effect of time and temperature on this three- component pre-mixture when developing the assay.

For Technical Support on this and other Drug Discovery Products, dial 760-603-7200, option 3, extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to www.invitrogen.com/lanthascreen

# 5.5 Incubation Conditions

#### Incubation Time

The incubation time can be set by the user. As a guide, results for various time points using GW4064 as the agonist are shown in Table 1 below. The  $EC_{50}$  and Z' are stable from 1 to 24 hours (Table 1).

Incubation Time (hours)	EC <sub>50</sub> GW4064	Z'-Factor
1	55 nM	0.75
2	55 nM	0.83
4	57 nM	0.83
6	60 nM	0.84
24	65 nM	0.81

**Table 1. Effect of Incubation Time on Assay Performance.** Sample data represents mean values from 3 separate experiments (n = 24). EC<sub>50</sub> values were determined by fitting the data to a sigmoidal dose response (variable slope) equation in GraphPad Prism 4.0 (data not shown). Z'-factors were calculated using the method of Zhang *et al.* (Zhang *et al.*, 1999) on the 24 replicates of maximum inverse agonist and no inverse agonist. Z'-factor is an indication of the robustness of the assay, where values  $\geq 0.5$  indicate an excellent assay, while a value of 1 indicates a theoretically ideal assay with no variability.

#### Temperature

We recommend that assays be conducted at room temperature (20–23°C).

For Technical Support for this or other Invitrogen Drug Discovery Solutions Products, dial 760-603-7200 extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to <u>www.invitrogen.com/lanthascreen</u>

# 6. AGONIST ASSAY

The procedure in this section describes a method for determining the  $EC_{50}$  of an agonist (n = 4) using the agonist-induced recruitment of fluorescein-SRC2-2 coactivator peptide to FXR-LBD. The only variable is the agonist concentration. All other assay components (FXR-LBD, peptide, Tb anti-GST antibody) are fixed at concentrations optimized to provide an assay with the lowest  $EC_{50}$  with the control agonist. The recommended final concentrations for optimal assay performance are listed in the following table. If component concentrations are changed, the assay must be reoptimized. The data in Figure 3 was generated using a method adapted from this protocol.

Component	Final Assay Concentration
Fluorescein-SRC2-2	500 nM
Tb anti-GST antibody	5 nM
FXR LBD-GST	5 nM

#### 6.1 Agonist Assay—Procedure

#### Prepare Complete Coregulator buffer I and Agonist Controls

- *Note:* The FXR-LBD should be thawed on ice just prior to use. Equilibrate all other assay components to room temperature.
- 1. Prepare Complete Coregulator buffer G by adding 1 M DTT to Coregulator buffer G for a final concentration of 10 mM DTT. Complete Coregulator buffer G must be prepared fresh daily.

Example: Add 60 µl of 1 M DTT to 5.94 ml of Coregulator buffer G

- *Note:* Keep Complete Coregulator buffer G at room temperature for the preparation of all reagents except for 4X FXR-LBD, which should be prepared with cold buffer. Reserve an appropriate volume of buffer **on ice** for preparation of 4X FXR-LBD.
- Prepare a solution of agonist at 100X of the final desired maximum starting concentration using DMSO.

*Example:* If the final desired maximum starting concentration of agonist is  $10 \,\mu$ M, prepare a solution of  $1000 \,\mu$ M agonist in DMSO.

#### Prepare 2X Agonist Dilution Series

1. Prepare a 12-point 100X dilution series of agonist in a 96-well plate by serially diluting the 100X agonist solution from step 3 three-fold using DMSO.

*Example:* Add 20  $\mu$ l of DMSO to wells A2–A12 in a 96-well polypropylene plate. To well A1, add 30  $\mu$ l of the 100X agonist solution prepared in step 3. Perform a three-fold serial dilution by transferring 10  $\mu$ l of the 100X agonist solution from well A1 to the 20  $\mu$ l of DMSO in well A2. Mix by pipetting up and down. Repeat for wells A2–A12.

2. Dilute each 100X agonist serial dilution from step 1 to 2X using Complete Coregulator buffer G.

*Example:* Transfer 5 µl of each of the 100X agonist serial dilutions from row A of the 96-well plate (wells A1–A12) to row B (wells B1–B12). Add 245 µl of Complete Coregulator buffer G to each well in row B of the 96-well plate. Mix by pipetting up and down.

3. Transfer 10  $\mu$ l of each of the 2X agonist serial dilutions to duplicate columns of rows A and B of a 384-well assay plate as shown in the plate layout in **Section 6.2**. (n = 4).

*Example:* Columns 1 and 2 of rows A and B of the 384-well assay plate will receive 10-µl aliquots from well B1 of the 96-well assay plate, wells A and B in columns 3 and 4 of the 384-well assay plate will receive 10-µl aliquots from well B2 of the 96-well assay plate, and so on.

#### Prepare 4X FXR-LBD

1. Prepare 4X FXR-LBD using **cold** Complete Coregulator buffer G from step 1. The recommended molar concentration of FXR for this kit is listed on the Certificate of Analysis. *Never vortex the FXR-LBD stock or dilutions*. Mix by pipetting or gentle inversion. Keep this solution on ice until needed for use in the assay.

*Example:* If the FXR-LBD has a stock concentration of 20000 nM and the recommended concentration for this kit is 5 nM, prepare a 4X solution at 20 nM by adding 5  $\mu$ l of FXR-LBD stock to 4995  $\mu$ l of **cold** Complete Coregulator buffer G.

2. Add 5 µl of 4X FXR-LBD to rows A–D, columns 1–24 of the 384-well assay plate.

#### Prepare 4X Fluorescein-SRC2-2/4X Tb anti-GST Antibody

1. Prepare a solution containing 2.0  $\mu$ M fluorescein-SRC2-2 (4X) and 20 nM Tb anti-GST antibody (4X) using Complete Coregulator buffer G at room temperature. The stock concentration of fluorescein-SRC2-2 is 100  $\mu$ M and the concentration of Tb anti-GST antibody is indicated on both the vial label and the Certificate of Analysis (1 mg/ml = ~6.7  $\mu$ M antibody).

*Example:* Add 40 µl of 100 µM fluorescein-SRC2-2 and 6.0 µl of 6.7 µM Tb anti-GST antibody to 1954 µl of Complete Coregulator buffer G.

Add 5 μl of 4X peptide/4X antibody solution to rows A–D, columns 1–24 of the 384-well assay plate (see plate layout in Section 6.2).

#### Plate Incubation and Reads

- 1. Gently mix the 384-well plate on a plate shaker and incubate at room temperature protected from light. The plate may be sealed with a cover to minimize evaporation.
- Read the plate between 1 and 24 hours at wavelengths of 520 nm and 495 nm, using the instrument settings described in Section
   If using a plate seal, spin the sealed plate in a centrifuge with an appropriate balance to spin down any condensation on the bottom of the seal, shake the plate gently, and read. Then proceed to data analysis as described in the next section.

# 6.2 Plate Layout



# 6.3 Agonist Assay—Data Analysis

Calculate the TR-FRET ratio by dividing the emission signal at 520 nm by the emission signal at 495 nm. Generate a binding curve by plotting the emission ratio vs. the log [ligand]. To determine the  $EC_{50}$  value, fit the data using an equation for a sigmoidal dose response (varying slope), as provided by GraphPad<sup>TM</sup> Prism 4.0 or another comparable graphing program.

The "maximum agonist" and "no agonist" control data can be used to calculate Z'-factor based on the equation of Zhang et al., 1999).

*Note:* The ligand EC<sub>50</sub> determined in the assay is a composite of multiple equilibria, including ligand binding to receptor and peptide binding to ligand/receptor complex.

For Technical Support for this or other Invitrogen Drug Discovery Solutions Products, dial 760-603-7200 extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to <u>www.invitrogen.com/lanthascreen</u>

Page 8 of 10

## 6.4 Agonist Assay—Reagent Volumes and Controls

The following table summarizes the reagent volumes, order of addition, and potential controls for developing an agonist mode assay. The protocol in **Section 6.1** may be adapted to screen test compounds using the following summary table as a guide.

Assay	Reagent Additions	Purpose
Test Compound	<ol> <li>10 μl 2X Test Compound (or dilution series)</li> <li>5 μl 4X FXR-LBD</li> <li>5 μl 4X FI-SRC2-2/Tb anti-GST Ab</li> </ol>	Assess coactivator recruitment upon binding of test compound.
Positive Control	<ol> <li>10 μl 2X agonist (or dilution series)</li> <li>5 μl 4X FXR-LBD</li> <li>5 μl 4X FI-SRC2-2 /Tb anti-GST Ab</li> </ol>	Assess coactivator recruitment upon binding of a known FXR agonist.
Negative Control	<ol> <li>10 μl 2X Compound solvent</li> <li>5 μl 4X FXR-LBD</li> <li>5 μl 4X FI-SRC2-2 /Tb anti-GST Ab</li> </ol>	Provides indication of basal signal or possible ligand-independent coactivator recruitment and accounts for possible interference from a compound's vehicle solvent.
No FXR-LBD Control	<ol> <li>10 μl 2X compound solvent</li> <li>5 μl Complete Coregulator buffer G</li> <li>5 μl 4X Fl-SRC2-2 /Tb anti-GST Ab</li> </ol>	Provides absolute bottom baseline for assay, not accounting for ligand-independent coactivator recruitment.

# 6.5 Agonist Assay—Representative Data



**Figure 3. Representative experiment of LanthaScreen**<sup>TM</sup> **TR-FRET FXR-Coactivator assay.** Assay composition: serial dilution of agonist GW4064, 5 nM FXR-LBD, 500 nM Fluorescein-SRC2-2, and 5 nM Tb anti-GST antibody. Results for 1-hour, 2-hour, 4-hour, 6-hour and 24-hour incubations are shown with the corresponding  $EC_{50}$  value. The curves were generated using a sigmoidal dose response (variable slope) equation in GraphPad<sup>TM</sup> Prism<sup>®</sup> 4.0.

*Note:* The ligand EC<sub>50</sub> determined in the assay is a composite of multiple equilibria, including ligand binding to receptor and peptide binding to ligand/receptor complex.

For Technical Support on this and other Drug Discovery Products, dial 760-603-7200, option 3, extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to <u>www.invitrogen.com/lanthascreen</u>

### 7. ASSAY PHARMACOLOGY



- Fexaramine, 63 nM
- GW4064, 41 nM
- CDCA, 3.2 μM
- Cholic acid, ND
- × UDCA, 40 μM



- *Note:* Binding of different ligands may result in different conformations in the nuclear receptor and thus different affinities for the coregulator peptide. A lower peptide affinity will result in a decreased TR-FRET signal and a lower plateau in the dose response curve.
- *Note:* The ligand EC<sub>50</sub> determined in the assay is a composite of multiple equilibria, including ligand binding to receptor and peptide binding to ligand/receptor complex.

# 8. **REFERENCES**

- Voegel, J. J., Heine, M. J., Zechel, C., Chambon, P., and Gronemeyer, H. (1996) TIF2, a 160 kDa transcriptional mediator for the ligand-dependent activation function AF-2 of nuclear receptors. *Embo J*, 15, 3667-3675
- Zhang, J. H., Chung, T. D., and Oldenburg, K. R. (1999) A simple statistical parameter for use in evaluation and validation of high-throughput screening assays. J. Biomol. Screen., 4, 67-73

For Technical Support for this or other Invitrogen Drug Discovery Solutions Products, dial 760-603-7200 extension 40266 For information on frequently asked questions regarding the LanthaScreen<sup>™</sup> technology, please go to <u>www.invitrogen.com/lanthascreen</u>

# 9. NOTICE TO PURCHASER

#### Limited Use Label License No. 176: Lanthanide Chelates

This product is the subject of U.S. Patent No. 5,622,821, 5,639,615 and 5,656,433 and foreign equivalents licensed exclusively to Invitrogen Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use, or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

The performance of this product is guaranteed for six months from the date of purchase if stored and handled properly.

This product and/or the use of this product is covered under one or more of the following US patents:

5,622,821, 5,639,615, 5,656,433, and 4,822,733, as well as pending patent applications owned by or licensed by Invitrogen Corporation.

Prism<sup>®</sup> is a registered trademark of GraphPad<sup>™</sup> Software, Inc., San Diego, CA.

Page 10 of 10

©2007 Invitrogen Corporation. All rights reserved. Reproduction without permission forbidden.