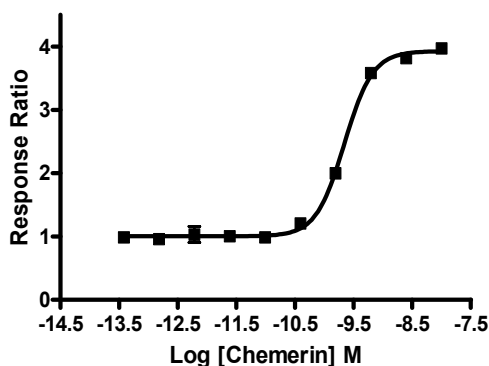


Table of Contents	Page
1. Description .....	1
2. Overview of Tango™ GPCR Cell-Based Assays .....	2
3. Materials Supplied .....	3
4. Materials Required .....	4
5. Detailed Cell Handling Procedures .....	5
5.1 Thawing Method.....	5
5.2 Propagation Method .....	5
5.3 Freezing Method .....	5
6. Media Requirements .....	6
7. Assay Procedure .....	7
7.1 Quick Assay Reference Guides.....	7
7.2 Detailed Assay Protocol .....	8
7.3 Substrate Preparation, Loading and Incubation .....	9
7.4 Detection .....	9
8. Data Analysis.....	10
8.1 Background Subtraction and Ratio Calculation .....	10
8.2 Visual Observation of Intracellular Beta-lactamase Activity Using LiveBLazer™ -FRET B/G Substrate (CCF4-AM)....	10
9. References .....	11
10. Purchaser Notification .....	12

## 1. Description

Tango™ GPR1-*bla* U2OS cells contain the human G protein-coupled receptor 1 linked to a TEV protease site and a Gal4-VP16 transcription factor stably integrated into the Tango™ GPCR-*bla* U2OS parental cell line. This parental cell line stably expresses a beta-arrestin/TEV protease fusion protein and the beta-lactamase reporter gene under the control of a UAS response element. The Tango™ GPR1-*bla* U2OS cells have been functionally validated for a response to Chemerin.



EC <sub>50</sub>	119pM
Z'-factor at EC <sub>100</sub>	0.70

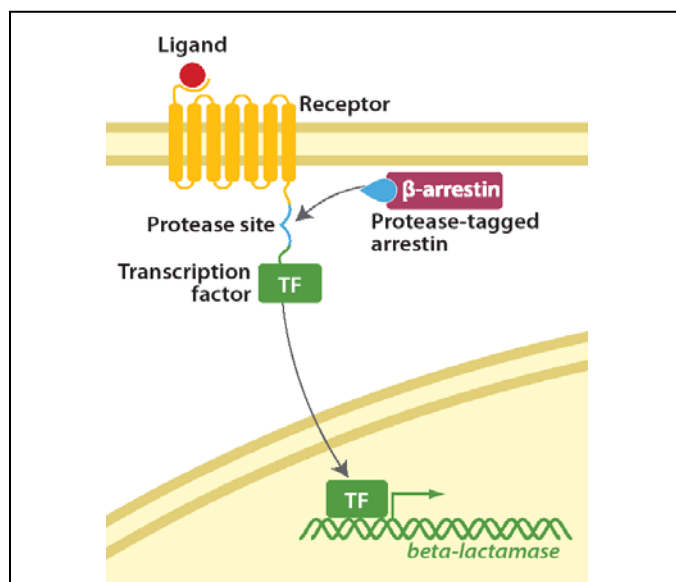
Dose response of Tango™ GPR1-*bla* U2OS cells to Chemerin.

## 2. Overview of Tango™ GPCR Cell-Based Assays

The Tango™ GPCR Assay technology combines the benefits of the Tango™ assay platform with the highly accurate, sensitive, and easy-to-use GeneBLAzer® beta-lactamase reporter system. The Tango™ assay platform is based upon ligand binding to G-Protein Coupled Receptors (GPCRs) that triggers desensitization, a process mediated by the recruitment of intracellular arrestin proteins to the activated receptor. As a result, the ligand-induced activation of GPCRs may be assayed by monitoring the interaction of arrestin with the test GPCR. A major advantage of this approach is that it does not depend on knowledge of the G-protein signaling specificity of the target receptor.

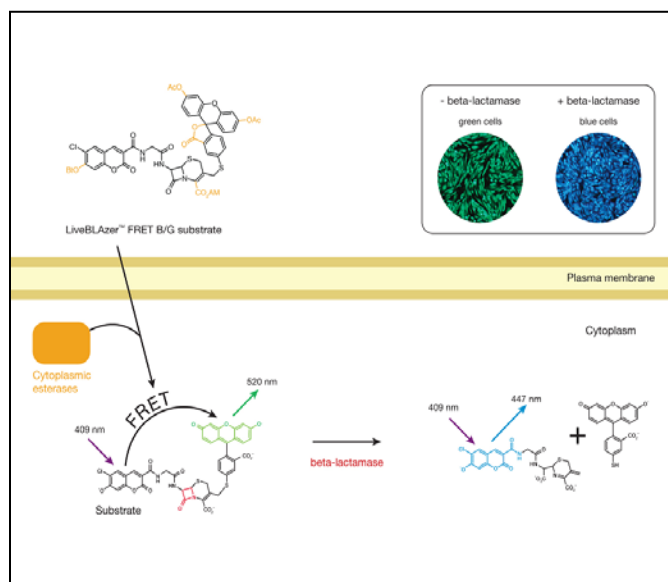
The design of the Tango™ GPCR assay is shown in Figure 1. The target GPCR is fused at its intracellular C-terminus to an exogenous transcription factor. Interposed between the receptor and the transcription factor is a specific cleavage sequence for a non-native protease. This chimeric receptor protein is expressed in a cell line containing the *bla* reporter gene responsive to the transcription factor. The cell line also expresses an arrestin-protease fusion protein that recognizes and cleaves the site between the receptor and transcription factor. The assay is performed by adding a ligand to the growing cells for a defined period and measuring the activity of the reporter gene. Activation of the reporter gene provides a quantifiable measurement of the degree of interaction between the target receptor and the protease-tagged arrestin partner. Additionally, it is unaffected by other signaling pathways in the cell, thus providing a highly selective readout of target receptor activation.

Figure 1.



**Figure 1—Tango™ GPCR cell-based beta-arrestin signaling.** Upon ligand binding and receptor activation, a protease-tagged beta-arrestin molecule is recruited to the GPCR that has been modified at the C-terminus to include a transcription factor linked by a protease cleavage site. The protease cleaves the transcription factor from the GPCR, which translocates to the nucleus and activates the expression of beta-lactamase.

Figure 2.



**Figure 2—Sensitive fluorescent detection of beta-lactamase reporter gene response using GeneBLAzer® technology.** After loading the cell permeable LiveBLAzer™ FRET B/G substrate, cellular fluorescence is measured. In the absence of beta-lactamase expression, cells generate green fluorescence. In the presence of beta-lactamase expression, the substrate is cleaved and the cells generate blue fluorescence.

The Tango™ assay technology uses a mammalian-optimized beta-lactamase (*bla*) reporter gene combined with a FRET-enabled substrate to provide reliable and sensitive detection in cells (1) (Figure 2). Cells are loaded with an engineered fluorescent substrate containing two fluorophores, coumarin and fluorescein. In the absence of *bla* expression, the substrate molecule remains intact. In this state, excitation of the coumarin results in fluorescence resonance energy transfer to the fluorescein moiety and emission of green fluorescent light. However, in the presence of *bla* expression, the substrate is cleaved separating the fluorophores and disrupting energy transfer. Excitation of the coumarin in the presence of *bla* enzyme activity results in a blue fluorescence signal.

The resulting coumarin:fluorescein ratio provides a normalized reporter response that can minimize experimental noise that masks subtle underlying biological responses of interest. The Tango™ assay technology has been proven in high-throughput screening (HTS) campaigns for a range of target classes, including G-protein coupled receptors (GPCRs) (2, 3), nuclear receptors (4-6) and kinase signaling pathways (7).

### 3. Materials Supplied

Product:	Name	Size	Catalog #
	<b>Tango™ GPR1-<i>bla</i> U2OS cells</b> Includes: <ul style="list-style-type: none"><li>• Tango™ GPR1-<i>bla</i> U2OS cells (K1839A)</li><li>• Protocol</li><li>• Certificate of Analysis (lists the number of cells provided)</li></ul>	1 tube	K1839
<b>Shipping Condition:</b>	Dry ice		
<b>Storage Condition of Cells:</b>	Liquid nitrogen. <b>Immediately upon receipt, cells must be stored in liquid nitrogen or thawed for immediate use. Cells stored at -80°C can quickly lose viability.</b>		
<b>Growth Properties of Cells:</b>	Adherent		
<b>Cell Phenotype:</b>	Epithelial		
<b>Selection Marker(s) for Cells:</b>	Zeocin™ 200 µg/mL, Geneticin® 100 µg/mL, and Hygromycin 50 µg/mL		
<b>Mycoplasma Testing:</b>	Negative		
<b>BioSafety Level:</b>	1		

## 4. Materials Required

Use the table below to determine the additional media and reagents required for use with each kit:

Media/Reagents	Recommended Source	Part #
LiveBLAzer™-FRET B/G Loading Kit: LiveBLAzer™-FRET B/G Substrate (CCF4-AM) DMSO for Solution A Solution B Solution C	Invitrogen	K1427 (70 µg) K1095 (200 µg) K1096 (1 mg) K1030 (5 mg)
Solution D	Invitrogen	K1156 (1 mL) K1157 (25 mL)
Recovery™ Cell Culture Freezing Medium	Invitrogen	12648-010
FreeStyle™ Expression Medium	Invitrogen	12338-018
McCoy's 5A Medium (modified) (1X)	Invitrogen	16600-082
DMSO	Fluka	41647
Fetal bovine serum (FBS), dialyzed, ( <b>DO NOT SUBSTITUTE!</b> )	Invitrogen	26400-036
Non-essential amino acids (NEAA)	Invitrogen	11140-050
Penicillin/Streptomycin (antibiotics)	Invitrogen	15140-122
Sodium Pyruvate	Invitrogen	11360-070
Phosphate-buffered saline without calcium and magnesium [PBS(-)]	Invitrogen	14190-136
HEPES (1 M, pH 7.3)	Invitrogen	15630-080
0.05% Trypsin/EDTA	Invitrogen	25300-054
Chemerin	R&D Systems	2324-CM
Zeocin™	Invitrogen	R250-01
Hygromycin	Invitrogen	10687-010
Geneticin®	Invitrogen	10131-027

The following table lists additional items for use with all kits:

Consumables	Recommended Source	Part #
Black-wall, clear-bottom, 384-well assay plates (with low fluorescence background)	Corning	3712
Compressed air	Various	---
Equipment	Recommended Source	
Fluorescence plate reader with bottom-read capabilities	Various	
Filters if required for plate reader (see <b>Section 7.4.1</b> )	Chroma Technologies	
<b>Optional:</b> Epifluorescence- or fluorescence-equipped microscope, with appropriate filters	Various	
<b>Optional:</b> Microplate centrifuge	Various	

## 5. Detailed Cell Handling Procedures

---

**Note:** Refer to **Section 6, Media Requirements** for specific media recipes.

### 5.1 Thawing Method

**Note:** Cells are shipped to you on dry ice and as such may require a short period of time prior to full recovery and normal growth.

1. The number of cells provided is listed on the Certificate of Analysis. For  $2 \times 10^6$  cells, add 14 mL of Thawing Medium to a T75 flask; for  $8 \times 10^6$  cells, add 30 mL of Thawing Medium to a T225 flask.
2. Place the flask in a humidified 37°C/5% CO<sub>2</sub> incubator for 15 minutes to allow medium to equilibrate to the proper pH and temperature.
3. Remove the vial of cells to be thawed from liquid nitrogen and rapidly thaw by placing at 37°C in a water bath with gentle agitation for 1–2 minutes. Do not submerge vial in water.
4. Decontaminate the vial by wiping with 70% ethanol before opening in a Class II biological safety cabinet.
5. Transfer the vial contents drop-wise into 10 mL of Thawing Medium in a sterile 15-mL conical tube.
6. Centrifuge cells at  $200 \times g$  for 5 minutes.
7. Aspirate supernatant and resuspend the cell pellet in 1 mL of fresh Thawing Medium.
8. Transfer the resuspended cells to the tissue culture flask containing pre-equilibrated Thawing Medium (from Step 2) and place the flask into the humidified 37°C/5% CO<sub>2</sub> incubator.
9. At first passage, switch to Growth Medium.

### 5.2 Propagation Method

1. Passage or feed cells at least twice a week. Maintain cells between 25% and 95% confluence. Do not allow cells to reach confluence.
2. To passage cells, aspirate medium, rinse once in PBS, add 0.05% Trypsin/EDTA (3 mL for a T75 flask, 5 mL for a T175 flask, and 7 mL for T225 flask) and swirl to coat the cells evenly. Cells usually detach after ~2–5 minutes exposure to 0.05% Trypsin/EDTA. Add an equal volume of Growth Medium to inactivate 0.05% Trypsin/EDTA.
3. Verify under a microscope that cells have detached and clumps have completely dispersed.
4. Centrifuge cells at  $200 \times g$  for 5 minutes and resuspend in Growth Medium.

### 5.3 Freezing Method

1. Harvest the cells as described in **Section 5.2** (above), Step 2. After detachment, count the cells, centrifuge cells at  $200 \times g$  for 5 minutes, and resuspend in 4°C Freeze Medium to a density of  $8 \times 10^6$  cells/mL.
2. Dispense 1.0-mL aliquots into cryogenic vials.
3. Place in an insulated container for slow cooling and store overnight at –80°C.
4. Transfer to liquid nitrogen the next day for storage.

## 6. Media Requirements

**Note:** Unless otherwise stated, have all media and solutions at least at room temperature (we recommend 37°C for optimal performance) before adding to cells.

**Note:** Make **NO MEDIA SUBSTITUTIONS**, as these cell lines have been specifically validated for optimal assay performance with these media. For dividing cells, we recommend that you create and store an aliquot for back up.

Component	Assay Medium	Growth Medium	Thawing Medium	Freeze Medium
FreeStyle™ Expression Medium	100%	—	—	—
McCoy's 5A Medium	—	90%	90%	—
Dialyzed FBS (Do not substitute!)	—	10%	10%	—
NEAA	—	0.1 mM	0.1 mM	—
HEPES (pH 7.3)	—	25 mM	25 mM	—
Sodium Pyruvate	—	1 mM	1 mM	—
Penicillin (antibiotic)	—	100 U/mL	100 U/mL	—
Streptomycin (antibiotic)	—	100 µg/mL	100 µg/mL	—
Recovery™ Cell Culture Freezing Medium	—	—	—	100%
Zeocin™	—	200 µg/mL	—	—
Hygromycin	—	50 µg/mL	—	—
Geneticin®	—	100 µg/mL	—	—

## 7. Assay Procedure

The following instructions outline the recommended procedure for determining activity of compounds as modulators of GPR1 using LiveBLAzer™-FRET B/G Substrate as the readout. If alternative substrates are used (*e.g.*, ToxBLAzer™ DualScreen or LyticBLAzer™ Loading kits), follow the loading protocol provided with the product.

### 7.1 Quick Assay Reference Guides

For a more detailed assay protocol, see **Section 7.2**.

#### Agonist Assay Quick Reference Guide

	Unstimulated Wells	Stimulated Wells	Cell-free Wells	Test Compound Wells
<b>Step 1</b> Plate cells, incubate	32 $\mu$ L cells in Assay Medium (10,000 cells/well)	32 $\mu$ L cells in Assay Medium (10,000 cells/well)	32 $\mu$ L Assay Medium (no cells)	32 $\mu$ L cells in Assay Medium (10,000 cells/well)
Incubate cells for 48 hrs. at 37°C / 5%CO <sub>2</sub>				
<b>Step 2</b> Add Agonist or Test Compounds	8 $\mu$ L Assay Medium with 0.5% DMSO	8 $\mu$ L 5X agonist in Assay Medium with 0.5% DMSO	8 $\mu$ L Assay Medium with 0.5% DMSO	8 $\mu$ L 5X Test Compounds in 0.5% DMSO
<b>Step 3</b> Incubate cells	Incubate in a humidified 37°C/5% CO <sub>2</sub> incubator for 24 hours			
<b>Step 4</b> Prepare 6X Substrate Mix	6 $\mu$ L of 1 mM LiveBLAzer™-FRET B/G (CCF4-AM) Substrate + 60 $\mu$ L of solution B, mix. Add 904 $\mu$ L of Solution C, mix. Add 30 $\mu$ L of Solution D, mix.			
<b>Step 5</b> Add Substrate Mixture	8 $\mu$ L per well			
<b>Step 6</b> Incubate Substrate Mix + cells	2 hours at room temperature in the dark			
<b>Step 7</b> Detect activity	See <b>Section 7.4</b>			
<b>Step 8</b> Analyze data	See <b>Section 8</b>			

#### Antagonist Assay Quick Reference Guide

	Unstimulated Wells	Stimulated Wells	Antagonist Control Wells	Cell-free Wells	Test Compound Wells
<b>Step 1</b> Plate cells, incubate	32 $\mu$ L cells in Assay Medium (10,000 cells/well)	32 $\mu$ L cells in Assay Medium (10,000 cells/well)	32 $\mu$ L cells in Assay Medium (10,000 cells/well)	32 $\mu$ L Assay Medium (no cells)	32 $\mu$ L cells in Assay Medium (10,000 cells/well)
Incubate cells for 48 hrs. at 37°C / 5%CO <sub>2</sub>					
<b>Step 2</b> Add Antagonist or Test Compounds, incubate	4 $\mu$ L Assay Medium with 0.5% DMSO	4 $\mu$ L Assay Medium with 0.5% DMSO	4 $\mu$ L 10X antagonist in Assay Medium with 0.5% DMSO	4 $\mu$ L Assay Medium with 0.5% DMSO	4 $\mu$ L 10X Test Compounds in Assay Medium with 0.5% DMSO
Incubate plate with Antagonist for 30 minutes before proceeding					
<b>Step 3</b> Add Agonist	4 $\mu$ L Assay Medium with 0.5% DMSO	4 $\mu$ L 10X agonist in Assay Medium with 0.5% DMSO	4 $\mu$ L 10X agonist in Assay Medium with 0.5% DMSO	4 $\mu$ L 10X agonist in Assay Medium with 0.5% DMSO	4 $\mu$ L 10X agonist in Assay Medium with 0.5% DMSO
<b>Step 4</b> Incubate cells	Incubate in a humidified 37°C/5% CO <sub>2</sub> incubator for 24 hours				
<b>Step 5</b> Prepare 6X Substrate Mix	6 $\mu$ L of 1 mM LiveBLAzer™-FRET B/G (CCF4-AM) Substrate + 60 $\mu$ L of solution B, mix. Add 904 $\mu$ L of Solution C, mix. Add 30 $\mu$ L of Solution D, mix.				
<b>Step 6</b> Add Substrate Mixture	8 $\mu$ L per well				
<b>Step 7</b> Incubate Mixture	2 hours at room temperature in the dark				
<b>Step 8</b> Detect activity	See <b>Section 7.4</b>				
<b>Step 9</b> Analyze data	See <b>Section 8</b>				

**For Technical Support for this or other Drug Discovery Products, dial 760-603-7200, option 3, extension 40266**

## 7.2 Detailed Assay Protocol

Plate layouts and experimental outlines will vary; in screening mode, we recommend using at least three wells for each control: Unstimulated Control, Stimulated Control, and Cell-free Control.

**Note:** Some solvents may affect assay performance. Assess the effects of solvent before screening. The cell stimulation procedure described below is carried out in the presence of 0.1% DMSO to simulate the effect that a Test Compound's solvent might have on the assay. If you use other solvents and/or solvent concentrations, optimize the following assay conditions appropriately.

### 7.2.1 Precautions

- Work on a dust-free, clean surface. Always handle the 384-well, black-wall, clear-bottom assay plate by the sides; do not touch the clear bottom of the assay plate.
- If pipetting manually, you may need to centrifuge the plate briefly at room temperature (for 1 minute at  $14 \times g$ ) after additions to ensure all assay components are on the bottom of the wells.

### 7.2.2 Plating Cells

1. Harvest cells and resuspend in Assay Medium to a density of 312,500 cells/mL.
2. Add 32  $\mu$ L per well of the Assay Medium to the Cell-free Control wells. Add 32  $\mu$ L per well of the cell suspension to the Test Compound wells, the Unstimulated Control wells, and Stimulated Control wells. Incubate cells at 37°C/ 5% CO<sub>2</sub> for 48 hours. Proceed to **Section 7.2.3** for an Agonist assay or **Section 7.2.4** for an Antagonist assay.

### 7.2.3 Agonist Assay Plate Setup

**Note:** This subsection provides directions for performing an Agonist assay. See **Section 7.2.4** for directions for performing an Antagonist assay.

1. Prepare a stock solution of 0.5% DMSO in Assay Medium.
2. Prepare a 5X stock of Test Compounds in Assay Medium with 0.5% DMSO.
3. Prepare a 5X stock of agonist in Assay Medium with 0.5% DMSO. We recommend running a dose response curve to determine the optimal concentration of the agonist solution.
4. Add 8  $\mu$ L of the stock solution of 0.5% DMSO in Assay Medium to the Unstimulated Control and Cell-free Control wells.
5. Add 8  $\mu$ L of the 5X stock solution of agonist to the Stimulated Control wells.
6. Add 8  $\mu$ L of the 5X stock of Test Compounds to the Test Compound wells.
7. Incubate the Agonist assay plate in a humidified 37°C/5% CO<sub>2</sub> incubator for 24 hours. Then proceed to **Section 7.3** for Substrate Loading and Incubation.

### 7.2.4 Antagonist Assay Plate Setup

**Note:** This subsection provides directions for performing an Antagonist assay. See **Section 7.2.3** for directions for performing an Agonist assay.

1. Prepare a stock solution of 0.5% DMSO in Assay Medium.
2. Prepare a 10X stock of Test Compounds in Assay Medium with 0.5% DMSO.
3. Prepare a 10X stock of agonist in Assay Medium with 0.5% DMSO. We recommend running a dose response curve to determine the optimal agonist concentration. For antagonist assays, we recommend stimulating cells initially with an agonist concentration in the EC<sub>50</sub>-EC<sub>80</sub> range.
4. Prepare a 10X stock of antagonist in Assay Medium with 0.5% DMSO. We recommend running a dose response curve to determine the optimal inhibition concentration for the Antagonist solution.
5. Add 4  $\mu$ L of the 10X stock of Test Compounds to the Test Compound wells.
6. Add 4  $\mu$ L of the stock solution of 0.5% DMSO to the Stimulated Control wells, the Unstimulated Control wells, and the Cell-free Control wells.
7. Add 4  $\mu$ L of the 10X stock of antagonist in Assay Medium with 0.5% DMSO to the Antagonist Control wells.
8. If desired, incubate the Test Compounds with the cells humidified 37°C/5% CO<sub>2</sub> incubator before proceeding. Typically, a 30-minute incubation is sufficient.
9. Add 4  $\mu$ L of the 10X stock solution of agonist to the Test Compound wells, the Stimulated Control wells, and the Antagonist Control wells.
10. Add 4  $\mu$ L of Assay Medium with 0.5% DMSO to the Unstimulated Control and Cell-free Control wells.



11. Incubate the Antagonist assay plate in a humidified 37°C/5% CO<sub>2</sub> incubator for 24 hours. Then proceed to **Section 7.3** for Substrate Loading and Incubation.

### 7.3 Substrate Preparation, Loading and Incubation

This protocol is designed for loading cells with LiveBLAzer™-FRET B/G Substrate Mixture (CCF4-AM) Substrate Mixture. If you use alternative substrates, follow the loading protocol provided with the substrate.

Prepare LiveBLAzer™-FRET B/G Substrate Mixture (CCF4-AM) Substrate Mixture and load cells in the absence of direct strong lighting. Turn off the light in the hood.

1. Prepare Solution A: 1 mM LiveBLAzer™-FRET B/G Substrate (CCF4-AM) Substrate Mixture in dry DMSO by adding 912 µL of DMSO per mg of dry substrate. Store the aliquots of the stock solution at -20°C until use. The molecular weight of the LiveBLAzer™-FRET B/G Substrate (CCF4-AM) is 1096 g/mol.
2. Prepare 6X Loading Solution:
  - a. Add 6 µL of Solution A to 60 µL of Solution B and vortex.
  - b. Add 904 µL of Solution C to the above solution and vortex.
  - c. Add 30 µL of Solution D to the above solution and vortex.
3. Remove assay plate from the humidified 37°C/5% CO<sub>2</sub> incubator.  
*Note:* Handle the plate gently and do not touch the bottom.
4. Add 8 µL of the 6X Substrate Mixture to each well.
5. Cover the plate to protect it from light and evaporation.
6. Incubate at room temperature for 2 hours.

### 7.4 Detection

Make measurements at room temperature from the bottom of the wells, preferably in 384-well, black-wall, clear-bottom assay plates with low fluorescence background. Before reading the plate, remove dust from the bottom with compressed air.

#### 7.4.1 Instrumentation, Filters, and Plates

- Fluorescence plate reader with bottom reading capabilities.
- Recommended filters for fluorescence plate reader:  
 Excitation filter: 409/20 nm  
 Emission filter: 460/40 nm  
 Emission filter: 530/30 nm

#### 7.4.2 Reading an Assay Plate

1. Set the fluorescence plate reader to bottom-read mode with optimal gain and 5 reads.
2. Allow the lamp in the fluorescence plate reader to warm up for at least 10 minutes before making measurements.
3. Use the following filter selections:

	Scan 1	Scan 2
<b>Purpose:</b>	Measure fluorescence in the Blue channel	Measure FRET signal in the Green channel
<b>Excitation filter:</b>	409/20 nm	409/20 nm
<b>Emission filter:</b>	460/40 nm	530/30 nm

## 8. Data Analysis

---

### 8.1 Background Subtraction and Ratio Calculation

We recommend that you subtract the background for both emission channels (460 nm and 530 nm).

1. Use the assay plate layout to identify the location of the Cell-free Control wells. These Control wells are used for background subtraction.
2. Determine the average emission from the Cell-free Control wells at both 460 nm (Average Blue Background) and 530 nm (Average Green Background).
3. Subtract the Average Blue background from all of the Blue emission data.
4. Subtract the Average Green background from all of the Green emission data.
5. Calculate the Blue/Green Emission Ratio for each well, by dividing the background-subtracted Blue emission values by the background-subtracted Green emission values.

### 8.2 Visual Observation of Intracellular Beta-lactamase Activity Using LiveBLazer™-FRET B/G Substrate (CCF4-AM)

**Note:** Microscopic visualization of cells will cause photobleaching. Always read the assay plate in the fluorescence plate reader before performing microscopic visualization.

An inverted microscope equipped for epifluorescence and with either a xenon or mercury excitation lamp may be used to view the LiveBLazer™-FRET B/G Substrate (CCF4-AM) signal in cells. To visually inspect the cells, you will need a long-pass filter passing blue and green fluorescence light, so that your eye can visually identify whether the cells are fluorescing green or blue.

Recommended filter sets for observing beta-lactamase activity are described below and are available from Chroma Technologies (800-824-7662, [www.chroma.com](http://www.chroma.com)).

#### Chroma Set # 41031

Excitation filter:	HQ405/20x (405 ± 10)
Dichroic mirror:	425 DCXR
Emission filter:	HQ435LP (435 long-pass)

Filter sizes vary for specific microscopes and need to be specified when the filters are ordered. For epifluorescence microscopes, a long-pass dichroic mirror is needed to separate excitation and emission light and should be matched to the excitation filter (to maximally block the excitation light around 405 nm, yet allow good transmission of the emitted light).

## 9. References

---

1. Zlokarnik, G., *et al*, **Quantitation of Transcription and Clonal Selection of Single Living Cells with Beta-Lactamase as Reporter**, (1998) *Science*; **279**: p84-88.
2. Kunapuli P, Ransom R, Murphy K, Pettibone D, Kerby J, Grimwood S, Zuck P, Hodder P, Lacson R, Hoffman I, Inglese J, Strulovici B, **Development of an Intact Cell Reporter Gene Beta-lactamase Assay for G Protein-coupled Receptors**, (2003) *Analytical Biochem.*; **314**: p16-29.
3. Xing, H., Pollok, B., *et al*, **A Fluorescent Reporter Assay For The Detection of Ligands Acting Through G1 Protein-coupled Receptors**, (2000) *J. Receptor & Signal Transduction Research*; **20**: p189-210.
4. Qureshi, S., *et al*, **A One-Arm Homologous Recombination Approach for Developing Nuclear Receptor Assays in Somatic Cells**, (2003) *Assay and Drug Dev. Tech.*; **1**: p755-766.
5. Peekhaus, N. *et al*, **A Beta-Lactamase-Dependent Gal4-Estrogen Receptor Transactivation Assay for the Ultra-High Throughput Screening of Estrogen Receptor Agonists in a 3,456-Well Format**, (2003) *Assay and Drug Dev Tech*, **1**: p789-800.
6. Chin, J., *et al*, **Miniaturization of Cell-Based, Beta-Lactamase-Dependent FRET Assays to Ultra-High Throughput Formats to Identify Agonists of Human Liver X Receptors**, (2003) *Assay and Drug Dev. Tech.*; **1**: p777-787.
7. Whitney M, Rockenstein E, Cantin G, Knapp T, Zlokarnik G, Sanders P, Durick K, Craig FF, Negulescu PA., **A Genome-wide Functional Assay of Signal Transduction in Living Mammalian Cells**, (1998) *Nat. Biotechnol.*; **16**: p1329-1333.

## 10. Purchaser Notification

### Limited Use Label License No. 5: Invitrogen Technology

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) not to 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. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product 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, Life Technologies is willing to accept return of the product with a full refund. For information about purchasing a license to use this product or the technology embedded in it for any use other than for research use please contact Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008 ; Phone (760) 603-7200 or e-mail: outlicensing@lifetech.com.

### Limited Use Label License No. 108: Lentiviral Technology

The Lentiviral Technology (based upon the lentikat™ system) is licensed from Cell Genesys, Inc., under U.S. Patent Nos. 5,834,256; 5,858,740; 5,994,136; 6,013,516; 6,051,427; 6,165,782 and 6,218,187 and corresponding patents and applications in other countries for internal research purposes only. Use of this technology for gene therapy applications or bioprocessing other than for non-human research use requires a license from Cell Genesys (Cell Genesys, Inc. 342 Lakeside Drive, Foster City, California 94404). 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), including non-gene therapy research and target validation applications in laboratory animals.

### Limited Use Label License No. 150: GeneBLAzer® Technology

This product and/or its use is the subject of one or more of U.S. Patent Nos. 5,741,657, 5,955,604, 6,291,162, and 6,472,205 and foreign equivalents licensed to Life Technologies Corporation. The right to use this product for internal research specifically excludes the right to use this product to identify, discover, and profile compounds that act as a flavor, fragrance or taste-enhancers and modify a target identified in taste, olfaction, or pheromone detection, which compound does not require FDA approval of an NDA for claims of safety and efficacy. The right to use methods claimed in the foregoing patents with this product for research purposes can only be acquired by the use of GeneBLAzer® substrates purchased from Life Technologies Corporation or its authorized distributors.

### Limited Use Label License No. 317: LentiVector® Technology

This product is licensed under U.S. Pat. Nos. 5,817,491; 5,591,624; 5,716,832; 6,312,682; 6,669,936; 6,235,522; 6,924,123 and foreign equivalents from Oxford BioMedica (UK) Ltd., Oxford, UK, and is provided for use in academic and commercial in vitro and in vivo research for elucidating gene function, and for validating potential gene products and pathways for drug discovery and development, but excludes any use of LentiVector® technology for: creating transgenic birds for the purpose of producing useful or valuable proteins in the eggs of such transgenic birds, the delivery of gene therapies, and for commercial production of therapeutic, diagnostic or other commercial products not intended for research use where such products do not consist of or incorporate a lentiviral vector. Information about licenses for commercial uses excluded under this license is available from Oxford BioMedica (UK), Ltd., Medawar Centre, Oxford Science Park, Oxford OX4 4GA UK enquiries@oxfordbiomedica.co.uk or BioMedica Inc 11622 EI Camino Real #100, San Diego CA 92130- 2049 USA. LentiVector is a registered US and European Community trade mark of Oxford BioMedica plc.

### Use of Genetically Modified Organisms (GMO)

**Information for European Customers** The Tango™ GPR1-*bla* U2OS DA and Tango™ GPR1-*bla* U2OS cell lines are genetically modified with the plasmids pTango™ β-Arr/TEV, pLenti-zeo/UAS-*bla* (note this construct was utilized as a plasmid not as a lentiviral stock) and pTangoGPR1. As a condition of sale, use of this product must be in accordance with all applicable local legislation and guidelines including EC Directive 90/219/EEC on the contained use of genetically modified organisms.

Zeocin™ is a trademark of CAYLA Corporation.

© 2009, 2010 Life Technologies Corporation. All rights reserved. Reproduction forbidden without permission.