

BacMam ERK2 [pThr185/pTyr187] Cellular Assay User Guide

Catalog no. A12888

Shipping: Varies Storage: Varies

Protocol part no. A12888.pps

Revision date: 10 January 2011

Table of Contents

Kit Contents and Handling	. 2
Overview LanthaScreen® Cellular Assay BacMam Technology	. 2
Before Starting Materials Required but Not Provided CHO Cell Culture Reagents (optional) U-2 OS Cell Culture Reagents (optional) Guidelines for Working with BacMam Reagent Guidelines for Optimizing LanthaScreen® Assays	. 3 . 4 . 4 . 4 . 5
Quick Reference Protocols for CHO Cells GPCR Agonist Assay—CHO Cells GPCR Antagonist Assay—CHO Cells	. 6
Quick Reference Protocols for U-2 OS Cells Agonist Assay—U-2 OS Cells Antagonist Assay—U-2 OS Cells	. 8
Detailed Protocol for GPCR-mediated ERK2 phosphorylation in CHO Cells	10
Detailed Protocol for ERK2 Phosphorylation in U-2 OS cells	12
LanthaScreen® Detection	14
Appendix	
Troubleshooting Guide	16
Purchaser Notification	17

Component	SKU#	Amount	Storage	Handling
BacMam ERK2 Reagent	A12889	2 × 25 mL	4°C	 Do not freeze Avoid extended exposure to ambient room light Use sterile technique Aliquot into sterile containers to minimize handling, if necessary
LanthaScreen® Tb-anti-ERK2 [pThr185/pTyr187] Antibody	PV5833*	10 µg	-20°C	Aliquot if necessary to avoid multiple freeze/thaw cycles
6X LanthaScreen [®] Cellular Assay Lysis Buffer	A12891	6 mL	4°C	On the day of assay, supplement with inhibitor cocktails (not provided; see page 3) and antibody
BacMam Enhancer Solution (1000X)	PV5835	150 μL	-20°C	Aliquot if necessary to avoid multiple freeze/thaw cycles

Kit Contents and Handling

*To order additional antibody, use catalog no. PV5269 (25 µg amount).

Overview

BacMam Cellular Assays use the BacMam gene delivery system in conjunction with LanthaScreen[®] Cellular Assays to measure post-translational modifications of a target substrate, as described below. The combination of the two technologies provides a fast, convenient, and robust method for interrogating specific signal transduction events in a cell background of choice.

LanthaScreen® Cellular Assay

LanthaScreen[®] Cellular Assays are high-throughput screening (HTS) compatible immunoassays used to interrogate target-specific post-translational modifications in a cell-based format. Target proteins are expressed as fusions with green fluorescent protein (GFP) in living cells, and modification-specific antibodies labeled with terbium (Tb) are used to detect stimulus-induced post-translational modifications in a time-resolved fluorescence resonance energy transfer (TR-FRET) format.

The use of GFP as a FRET acceptor circumvents the need to use complex antigen-capturing reagents, thereby providing a high-throughput alternative to commonly used analytical methods such as Western blot and ELISA.

For more information, visit www.invitrogen.com/lanthascreen.

BacMam Technology

BacMam technology uses a modified baculovirus to efficiently deliver and express genes (in this case, GFP-substrate fusion gene) in mammalian cells. The virus is non-replicating in mammalian cells, rendering them safe as research reagents.

This technology has several advantages over traditional transient methods for heterologous gene expression:

- High transduction efficiency across a broad range of cell types, including primary and stem cells
- Little-to-no observable cytopathic effects
- Reproducible and titratable target gene expression
- Compatibility with simultaneous delivery of multiple genes.

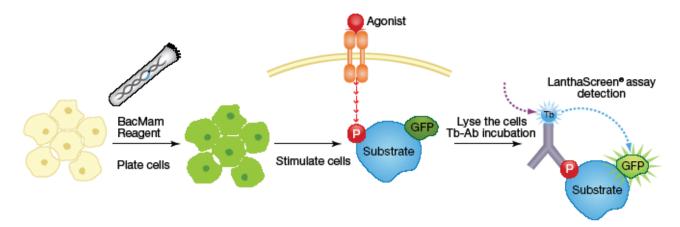


Figure 1. Illustration of Representative Assay Workflow. Cells are treated with BacMam reagent encoding a GFP-fusion protein and plated in 384-well format; 24 hours post-transduction, the cells are stimulated to induce post-translational modification of the GFP-substrate (e.g., phosphorylation as shown). Cells are then lysed in the presence of a terbium-anti-modification-specific antibody prior to the LanthaScreen[®] assay readout.

See Kost, TA et. al *Drug Disc. Today* **2007**, *12*, 396-403 for more information on BacMam gene expression in cells. For more information on BacMam, visit <u>www.invitrogen.com/bacmam</u>.

Intended Use

For research use only. Not intended for human or animal diagnostic or therapeutic uses.

Before Starting

Materials Required but Not Provided

Materials	Recommended Source	Part #
Positive Control Growth Factor We recommend: EGF (if using U-2 OS cells)	Invitrogen	PHG0314
Positive Control GPCR agonist We recommend: 5-HT (if using CHO cells expressing 5-HT1A)	Sigma	H9523
Cell line of interest	various	various
DMSO	Fluka	41647
Protease Inhibitor Cocktail	Sigma	P8340
Phosphatase Inhibitor Cocktail	Sigma	P0044
White tissue culture-treated, 384-well assay plates	Corning	3570
Fluorescence plate reader with top-read and TR-FRET capability	Visit <u>www.invitrogen.com/i</u>	<u>nstrumentsetup</u>
Optional: Clear-bottom, tissues culture treated, 384-well assay plates for visualization of GFP-fusion	Corning	3712

CHO Cell Culture Reagents (option	onal)
-----------------------------------	-------

Media/Reagents	Recommended Source	Part #		
GeneBLAzer [®] 5HT1A Gα15-NFAT- <i>bla</i> CHO-K1 Cells (CHO cells expressing 5-HT1A Receptor)	Invitrogen	K1712		
Dulbecco's Modified Eagle Medium (D-MEM) (1X), liquid (with GlutaMAX [™] -I)				
Opti-MEM [®] I Reduced-Serum Medium (with HEPES and L-glutamine, without Phenol Red)	Invitrogen 11058.			
Fetal Bovine Serum (dialyzed)	Invitrogen	26400-036		
Nonessential amino acids (NEAA)	Invitrogen	11140-050		
Sodium Pyruvate	Invitrogen	11360-070		
Penicillin/Streptomycin (antibiotic)	Invitrogen	15140-122		
Dulbecco's Phosphate-buffered saline (D-PBS) without Ca ²⁺ and Mg ²⁺	Invitrogen	14190-136		
Dulbecco's Phosphate-buffered saline (D-PBS) with Ca ²⁺ and Mg ²⁺	Invitrogen	14040-133		
HEPES Buffer Solution (1 M)	Invitrogen	15630-080		
Trypsin/EDTA	Invitrogen	25300-062		

U-2 OS Cell Culture Reagents (optional)

Media/Reagents	Recommended Source	Part #
U-2 OS cells	ATCC®	HTB-96 [™]
Opti-MEM [®] I Reduced-Serum Medium (with HEPES and L- glutamine, without Phenol Red)	Invitrogen	11058-021
McCoy's 5A Medium (modified) (1X), liquid	Invitrogen	16600-108
Fetal Bovine Serum (dialyzed)	Invitrogen	26400-036
Nonessential amino acids (NEAA)	Invitrogen	11140-050
Sodium Pyruvate	Invitrogen	11360-070
Penicillin/Streptomycin (antibiotic)	Invitrogen	15140-122
Dulbecco's Phosphate-buffered saline (PBS) without Ca ²⁺ and Mg ²⁺	Invitrogen	14190-136
HEPES Buffer Solution (1 M)	Invitrogen	15630-080
Trypsin/EDTA	Invitrogen	25300-062

Guidelines for Working with BacMam Reagent

- For first time users of BacMam reagent, we recommend including a control cell line that can be transduced exceptionally well, such as U-2 OS (ATCC[®] catalog number HTB-96[™]).
- Most cell types can be transduced efficiently using the protocol described for U-2 OS cells. However, some challenging cell types require the alternative protocol described for CHO cells.
- BacMam Enhancer Solution is not required for this BacMam reagent if using U-2 OS cells.
- We recommend the following steps for optimizing the transduction conditions for your cell background of interest in a small scale in order to minimize reagent consumption:
 - 1. Prepare a virus titration to determine the optimal percentage of virus to use.
 - 2. For untested cell backgrounds, perform a transduction in the presence and absence of BacMam Enhancer Solution included in the kit and then analyze the expression of GFP-ERK2.

B. CHO-5HT1A (5-HT stimulation)

Titration of BacMam ERK2 Reagent

A. U-2 OS (EGF stimulation)

We recommend performing a titration of the BacMam ERK2 Reagent to determine the optimal percentage of virus for the transduction in your cell background of interest. Select the lowest percentage of BacMam reagent that yields the largest assay window (response ratio). See the examples below.

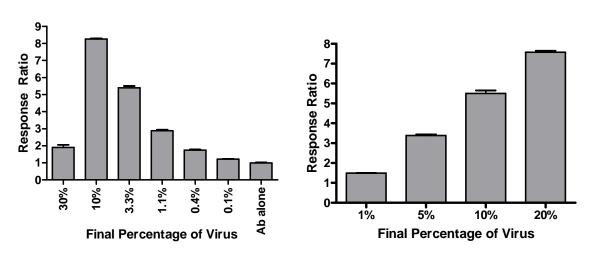


Figure 2. Detection of ERK2 phosphorylation at Thr185/Tyr187 in (**A**) U-2 OS cells transduced with various concentrations of BacMam ERK2 Reagent and stimulated with 100 ng/mL EGF; (**B**) CHO-5HT1A cells transduced with various concentrations of BacMam ERK2 and stimulated with 0.3 μ M 5-HT.

Guidelines for Optimizing LanthaScreen® Assays

First-time LanthaScreen® users

LanthaScreen[®] assays require the detection of terbium TR-FRET. For more information about your specific instrument and to purchase filters, visit <u>www.invitrogen.com/instrumentsetup</u>. For a protocol describing how to test whether your instrument is able to detect a terbium/GFP TR-FRET signal, please contact Drug Discovery Technical Support at <u>drugdiscoverytech@invitrogen.com</u> or 760-603-7200 (extension 40266).

Important assay parameters for optimization

- Confluence of cells at harvest for assay set-up may impact results, such as the assay window. In general, cells should be harvested at the maximum density at which they are still healthy.
- Cell plating density (i.e., cell number per well in the assay plate)
- Ligand stimulation time
- Assay equilibration time

Validation packet

Visit <u>www.invitrogen.com</u> and search for A12888 to download the validation packet for the BacMam ERK2 Cellular Assay. The validation packet is located under the "How to Use" tab on the product page, and contains more information about ERK2 applications.

Quick Reference Protocols for CHO Cells

The following agonist and antagonist GPCR-mediated ERK2 phosphorylation protocols have been developed for CHO cells over-expressing the serotonin receptor HTR1A. Conditions such as plating density, stimulant, or stimulation time may need to be optimized for CHO cell lines expressing different GPCRs. Phosphorylation of ERK2 at [Thr185/Tyr187] can also be induced in CHO cells with either 10% serum or 100 nM PMA.

		Cell-Free Control Wells	Unstimulated Control Wells	Stimulated Control Wells	Test Compound Wells	
BacMam Transduction	Step 1 Grow and Transduce Cells	 Grow cells in Growth Medium* to 90–100% confluence (~1–2 × 10⁵ cells/cm²). Perform five 3-fold serial dilutions of BacMam reagent in PBS with Ca²⁺ and Mg²⁺. Remove media from cells and add the serial diluted BacMam virus. Incubate at room temperature protected from light for 3–4 hrs. 				
BacMan	Step 2 Add Enhancer and Incubate Cells	Remove virus and add Incubate		plus 1X BacMam Er % CO ₂ for 20–24 hor		
	Step 3 Harvest Cells	Harvest cells, wash once 0.75 × 10 ⁶ cells/mL		m** and resuspend i mal cell density for		
	Step 4 Plate Cells	20 μL/well Assay20 μL transduced cells/well (about 15,000 cells/welMedia onlyquick spin of the plate				
	Step 5 Serum-starve and Incubate Cells	Incubate the plate at 37°C/5% CO ₂ for 4 hours				
Assay	Step 6 Prepare Complete 6X Lysis Buffer	To 1 mL 6X Lysis Buffer, add 30 μL of 100x protease inhibitor, 30 μL 100x phosphatase inhibitor , and Tb-anti-ERK2 [pThr185/pTyr187] Antibody to 12 nM .				
LanthaScreen® Assay	Step 7 Add Agonist	Assay Modia 5-HT (900 nM) in Test Compou			10 μL/well of 3X Test Compound in Assay Media	
anthaS	Step 8 Stimulate Cells	Incubate the plate at $37^{\circ}C/5\%$ CO ₂ for 6–8 minutes				
ت	Step8 Add Lysis Buffer (including Tb-Ab)	Add 6 μ L/ well of Complete 6X Lysis Buffer to each well				
	Step 9 Cell Lysis/Assay Equilibration	Incubate plate for ~3 hours at room temperature in the dark				
	Step 10 Read Plate and Analyze Data	See LanthaScreen [®] bandwidth); Emissio Media (Invitrogen 10569)	n filters: 490 nm (10 bandwie	lth)	d 520 nm (25 nm	

* Growth Media: D-MEM Media (Invitrogen 10569) with 10% dFBS, 10 mM HEPES, 0.1 mM NEAA, and 100 U/mL Penicillin/100 μ g/mL Streptomycin

** Assay Media: Opti-MEM[®] I (Invitrogen 11058) with 0.1% cdFBS (or dialyzed FBS), 0.1 mM NEAA, 1 mM Sodium Pyruvate, and 100 U/mL Penicillin/100 μg/mL Streptomycin

GPCR Antagonist Assay—CHO Cells

		Cell-Free Control Wells	Unstimulated Control Wells	Stimulated Control Wells	Control Compound wells	Test Compound Wells
BacMam Transduction	Step 1 Grow and Transduce Cells	 Grow cells in Growth Medium* to 90–100% confluence (~1–2 × 10⁵ cells/cm²). Perform five 3-fold serial dilutions of BacMam reagent in PBS with Ca²⁺ and Mg²⁺. Remove media from cells and add the serial diluted BacMam virus. Incubate at room temperature protected from light for 3–4 hrs. 				
Ba Tran	Step 2 Add Enhancer and Incubate Cells		Remove virus and add Growth Medium* plus 1X BacMam Enhancer Solution. Incubate the plate at 37°C/5% CO2 for 20–24 hours.			
	Step 3 Harvest Cells	Harvest cells, was 0.75×10^6 cells/mL				
	Step 4 Plate Cells	20 µL/well Assay Media only	20 µL transduced spin of the plate	d cells /well (abo	ut 15,000 cells/v	vell), quick
	Step 5 Serum-starve and Incubate Cells	Incubate the plate at $37^{\circ}C/5\%$ CO ₂ for 4 hours				
	Step 6 Prepare Complete Lysis Buffer	To 1 mL 6X Lysis Buffer, add 30 μL of 100x protease inhibitor, 30 μL 100x phosphatase inhibitor , and Tb-anti-ERK2 [pThr185/pTyr187] Antibody to 12 nM .				
LanthaScreen® Assay	Step 7 Add Antagonist	5 μL/well of5 μL/well of5 μL/well of5 μL/well of 0.5% DMSO in Assay Medium5 μL/well of5X controlin AssayAssayAssayMediumMedium				5X test compound in Assay
laScre	Step 8 Incubate	Incubate the plate at 37°C/5% CO ₂ for 30–45 minutes				
Lanth	Step 9 Add Agonist	5 μL/well Ass	ay Medium		Cagonist (for 5-H n Assay Mediun	
	Step10 Stimulate Cells	Incubate the plate at $37^{\circ}C/5\%$ CO ₂ for 6–8 minutes				
	Step 11 Add Lysis Buffer (including Tb-Ab)	Add 6 µL/ well of Complete 6X Lysis Buffer to each well				
	Step 12 Cell Lysis/Assay Equilibration	Incubate plate for ~3 hours at room temperature in the dark				
	Step 13 Read Plate and Analyze Data Growth Media: D-ME	See LanthaScreen [®] Detection, page 14—Excitation filter: 337 nm (30 nm bandwidth); Emission filters: 490 nm (10 nm bandwidth) and 520 nm (25 nm bandwidth)				andwidth)

* Growth Media: D-MEM Media (Invitrogen 10569) with 10% dFBS, 10 mM HEPES, 0.1 mM NEAA, and 100 U/mL Penicillin/100 μg/mL Streptomycin

** Assay Media: Opti-MEM[®] I (Invitrogen 11058) with 0.1% cdFBS (or dialyzed FBS), 0.1 mM NEAA, 1 mM Sodium Pyruvate, and 100 U/mL Penicillin/ 100 µg/mL Streptomycin

Quick Reference Protocols for U-2 OS Cells

In the following agonist and antagonist ERK2 phosphorylation protocols have been developed for U-2 OS cells. They may be applied to many other cell types such as HEK293, HeLa, and A549. The cell harvesting and plating densities, growth medium, and assay medium must be optimized for your particular cell line(s).

The cells are incubated with virus at the time of plating onto the assay plate. These protocols include a 20–24 hour incubation period.

Agonist Assay—U-2 OS Cells

		Cell-Free Control Wells	Unstimulated Control Wells	Stimulated Control Wells	Test Compound Wells	
BacMam Transduction	Step 1 Grow, Harvest Cells and Transduce	 Grow cells in Growth Medium* to 90-100% confluence (~0.8–1.2 × 10⁵ cells/cm²). Harvest and wash cells and resuspend in Assay Medium** at 8.4 × 10⁵ cells/mL. Perform five 3-fold serial dilutions of BacMam reagent in Assay Medium. Add 0.4 mL undiluted or serial diluted BacMam reagent to 1 mL cells to generate a virus titration range of 0.1% to 30% (v/v) final concentration. 				
icMam Tr	Step 2 Plate Cells/virus mixture	Add 20 μL/well Assay Media only		lls and BacMam mix s/well), and quick sj		
Ba	Step 3 Incubate Cells	Incubate	the plate at 37°C/5%	6 CO2 for 20–24 hour	s	
	Step 4 Prepare Complete 6X Lysis Buffer	To 1 mL 6X Lysis Buffer, add 30 μL of 100x protease inhibitor, 30 μL 100x phosphatase inhibitor , and Tb-anti-ERK2 [pThr185/pTyr187] Antibody to 12 nM				
® Assay	Step 5 Add media or ligand	Add 10 µL/well of Assay Media with 0.3% DMSO Add 10 µL/well of Assay Media with 0.3% DMSO ECE use			Add 10 μL/well of 3X Test Compound in Assay Media (0.3% DMSO)	
creen	Step 6 Stimulate Cells	Incubate the plate at 37°C/5% CO ₂ for 6–8 minutes				
LanthaScreen® Assay	Step 7 Add Lysis Buffer (including Tb-Ab)	Add 6 μ L/well of Complete 6X Lysis Buffer to each well				
	Step 8 Cell Lysis/Assay Equilibration	Incubate plate for ~3 hours at room temperature in the dark				
	Step 9 Read Plate and Analyze Data	See LanthaScreen® Detection , page 14—Excitation filter: 337 nm (30 nm bandwidth); Emission filters: 490 nm (10 nm bandwidth) and 520 nm (25 nm bandwidth)				

* Growth Media for U-2 OS Cells: McCoy's 5A Media with 10% dFBS, 10 mM HEPES, 0.1 mM NEAA, 1 mM Sodium Pyruvate, and 100 U/mL Penicillin/ 100 µg/mL Streptomycin

** Assay Media: Opti-MEM[®] I (Invitrogen 11058) with 0.1% cdFBS (or dialyzed FBS), 0.1 mM NEAA, 1 mM Sodium Pyruvate, and 100 U/mL Penicillin/ 100 μg/mL Streptomycin

Antagonist Assay—U-2 OS Cells

		Cell-Free Control Wells	Unstimulated Control Wells	Stimulated Control Wells	Control Compound wells	Test Compound Wells
BacMam Transduction	Step 1 Grow, Harvest Cells and Transduce	 Grow cells in Growth Medium* to 70–90% confluency (~0.8–1.2 × 10⁵ cells/cm²). Harvest, wash cells and resuspend in Assay Medium** at 8.4 × 10⁵ cells/mL. Perform five- 3 fold serial dilutions of BacMam reagent in Assay Medium. Add 0.4 mL undiluted or serial diluted BacMam reagent to 1 mL cells to generate a virus titration range of 0.1% to 30% (v/v) final concentration. 				
acMam Tr	Step 2 Plate Cells/virus mixture	Add 20 µL/well Assay Media only			acMam mixture p and quick spin th	
B	Step 3 Incubate Cells	Ir	ncubate the plate a	tt 37°C/5% CO2	for 20–24 hours	
	Step 4 Add media, control, or test compounds	Add 5 µL/well of 0.5% DMSO in Assay Media 5X Iressa (2 5 µM) in Compound			Add 5 µL/well of 5X Test Compound in Assay Media	
	Step 5 Compound pretreatment	Incubate the plate at 37°C/5% CO ₂ for 30–60 minutes				
Assay	Step 6 Prepare Complete 6X Lysis Buffer	To 1 mL 6X Lysis Buffer, add 30 μL of 100x protease inhibitor, 30 μL 100x phosphatase inhibitor , and Tb-anti-ERK2 [pThr185/pTyr187] Antibody to 12 nM				
LanthaScreen® Assay	Step 7 Add media or ligand	Add 5 μL/wellAdd 5 μL/well of 6X Agonist (for EGF, 600 ng/mL) in AssayAssay MediaMedia				;/mL) in Assay
antha	Step 8 Stimulate Cells	Incubate the plate at 37°C/5% CO ₂ for 6–8 minutes				
	Step 9 Add Lysis Buffer (including Tb-Ab)	Add 6 µL/well of Complete 6X Lysis Buffer to each well				
	Step 10 Cell Lysis/Assay Equilibration	Incubate plate for ~3 hours at room temperature in the dark				
	Step 11 Read Plate and Analyze Data	See LanthaScreen [®] Detection, page 14—Excitation filter: 337 nm (30 nm bandwidth); Emission filters: 490 nm (10 nm bandwidth) and 520 nm (25 nm bandwidth)				

* **Growth Media for U-2 OS Cells:** McCoy's 5A Media with 10% dFBS, 10 mM HEPES, 0.1 mM NEAA, 1 mM Sodium Pyruvate, and 100 U/mL Penicillin/ 100 µg/mL Streptomycin

** Assay Media: Opti-MEM[®] I (Invitrogen 11058) with 0.1% cdFBS (or dialyzed FBS), 0.1 mM NEAA, 1 mM Sodium Pyruvate, and 100 U/mL Penicillin/ 100 μg/mL Streptomycin

Detailed Protocol for GPCR-mediated ERK2 phosphorylation in CHO Cells

The following detailed protocol provides separate workflows for agonist and antagonist assays. Cells are allowed to adhere to the tissue-culture flasks before transduction with BacMam reagent. For easy-to-transduce cells, an alternative detailed transduction protocol is provided starting on page 14.

Day 1. BacMam Transduction

- 1. Begin with cells grown to near-complete confluence under normal growth conditions (e.g., CHO cells should be grown to 90–100% confluence, about 1–2 × 10⁵cells/cm²). **Confluence of cells may impact results, such as the assay window (i.e., less than 90% confluence may result in a significantly lower assay window).**
- 2. Perform five 3-fold serial dilutions of BacMam reagent in D-PBS with Ca²⁺ and Mg^{2+.}
- 3. Remove media from cells and add the serial diluted BacMam virus. A typical final concentration of BacMam reagent is 0.12-30% (v/v).
- 4. Incubate at room temperature for 3–4 hours in the dark.
- 5. Remove the BacMam reagent from the cells. Add growth medium containing 1X BacMam Enhancer Solution.
- 6. Incubate the cells in a $37^{\circ}C/5\%$ CO₂ incubator for 20–24 hours.

Day 2. LanthaScreen® Cellular Assay

- 7. Harvest the cells and wash once with assay medium, then resuspend the cells in assay medium at 0.75×10^6 cells/mL. You will need to determine the optimal cell density for your cell line of interest.
- 8. Plate 20 μ L/well cells in assay medium onto a 384-well assay plate (about 15,000 cells/well). Plate 20 μ L/well assay medium in cell-free control wells.
- 9. **Optional:** Transfer some cells to the wells of a clear-bottom 384-well plate for image analysis of GFP expression.
- 10. Incubate the cells in a $37^{\circ}C/5\%$ CO₂ incubator for ~4 hours for serum-starvation. Depending on the cell type, you may need to incubate longer.
- 11. **Optional:** Analyze GFP expression levels in the clear-bottom 384-well plate by fluorescence microscopy using standard FITC filter sets.
- During incubation, prepare Complete 6X Lysis Buffer by adding both protease inhibitor and phosphatase inhibitor cocktails to the provided 6X Lysis Buffer, at a 1:33 dilution of 100X stock (e.g., 30 μL of 100X stock inhibitors per 1,000 μL of 6X Lysis Buffer) and add LanthaScreen[®] Tb-anti-ERK2 [pThr185/pTyr187] Antibody to 12 nM Mix gently by inversion. Store on ice until use.
- 13. Proceed with the following agonist or antagonist assay setup.

Agonist Assay Setup:

- a. Prepare a stock solution of 0.3% DMSO in assay medium. **Note:** If you are using a solvent other than DMSO for the agonist, change the solvent used in the control wells accordingly. Be careful to keep the amount of solvent consistent in all wells.
- b. Prepare 3X control agonist in assay medium with 0.3% DMSO (e.g., if treating CHO-5HT1A cells with 5-HT, prepare 900 nM 5-HT in assay medium for a 3X concentration). Run a dose response curve to determine the EC₁₀₀ for your control agonist solution.
- c. Prepare 3X test compound in assay medium (if the test compound is dissolved in DMSO, make sure the DMSO concentration for the 3X solution is 0.3%).

- d. Add 10 μL of assay medium with 0.3% DMSO to each unstimulated control and cell-free control well.
- e. Add 10 µL of the 3X control agonist (5-HT) in assay medium to each stimulated control well.
- f. Add 10 µL of the 3X test compound in assay medium to each test compound well.
- g. Proceed to the next numbered step.

Antagonist Assay Setup:

- a. Prepare a stock solution of 0.5% DMSO in assay medium. **Note:** If you are using a solvent other than DMSO for the agonist, change the solvent used in the control wells accordingly. Be careful to keep the amount of solvent consistent in all wells.
- b. Prepare a 5X control antagonist in assay medium. Run a dose response curve to determine the IC_{100} for your control inhibitor solution.
- c. Prepare a 5X test compound in assay medium (if the test compound is dissolved in DMSO, make sure the DMSO concentration for the 5X solution is 0.5%).
- d. Add 5 μL of the assay medium with 0.5% DMSO to each unstimulated, stimulated, and cell-free control well.
- e. Add 5 µL of the 5X control antagonist in assay medium to each control compound well.
- f. Add 5 µL of the 5X test compound in assay medium to each test compound well.
- g. Incubate the plate in a humidified $37^{\circ}C/5\%$ CO₂ incubator for 30–60 minutes.
- a. Prepare a 6X stock of agonist in assay medium (e.g., if using CHO-5HT1A cells, prepare 180 nM 5-HT in assay medium for a 6X concentration). Run a dose response curve to determine the EC_{80} for your agonist solution.
- h. Add 5 µL of the assay medium to each unstimulated and cell-free control well.
- i. Add 5 µL of 6X stock of agonist (5-HT) to each stimulated, control, and test compound well
- 14. Cover the plate and stimulate the cells by incubating in a humidified $37^{\circ}C/5\%$ CO₂ incubator for 6–8 minutes.
- *Tip:* The stimulation time must be optimized for each cell type and agonist. "Stimulation time" refers to the time from agonist addition to the addition of Complete 6X Lysis Buffer.
- 15. After stimulation, immediately add 6 µL of Complete 6X Lysis Buffer to each well. Cover the plate.
- 16. Incubate the covered plate at room temperature in the dark for 3 hours or other desired equilibration time. The equilibration time can be optimized for your cell line of interest.
- *Note:* Assay plates may be stored at 4°C overnight prior to reading. Let the plate warm to room temperature prior to reading.
- 17. Proceed to LanthaScreen[®] Detection, page 14.

Detailed Protocol for ERK2 Phosphorylation in U-2 OS cells

The following detailed protocol provides separate workflows for agonist and antagonist assays. Cells are incubated with virus at the time of plating onto the assay plate. This protocol includes a 20–24-hour incubation.

Day 1. BacMam Transduction

- 1. Begin with cells grown to near-complete confluence under normal growth conditions (e.g., U-2 OS cells should be grown to 100% confluence). Confluence of cells may impact results, such as the assay window.
- *Note:* For many cell types, such as U-2 OS, a cell seeding density of ~30,000 cells/cm² for 3 days with a harvest density of ~ $0.8-1.2 \times 10^5$ cells/cm² is optimal. In general, cells should be transduced at the maximum density at which the cells are still healthy.
- 2. Harvest and wash the cells and resuspend in low-serum assay medium, using the appropriate conditions for your particular cell line (for U-2 OS cells, resuspend at 7×10^5 cells/mL).
- *Note:* For the BacMam ERK2 assay, it is important to wash away any residual serum following harvest to ensure optimal starve.
- 3. Add BacMam ERK2 Reagent to the cells. A typical final concentration of BacMam reagent is 1-30% (v/v). Mix gently by inversion.
- *Note:* We recommend testing a range of v/v dilutions of BacMam reagent. For U-2 OS cells, we recommend preparing five three-fold serial dilutions of the BacMam reagent in assay medium. Then mix 1 mL cells at 8.4×10^5 cells/mL with 0.4 mL of each BacMam reagent dilution. This should yield a final cell concentration of 6×10^5 cells/mL and final reagent concentrations of ~30%, 10%, 3%, 1%, 0.3%, and 0.1% (v/v).
- 4. Transfer 20 μL/well of cells/BacMam reagent mixture to a white 384-well assay plate. The number of cells per well may need to be optimized (e.g., seed U-2 OS cells at ~12,000 cells/well).
- 5. **Optional:** Transfer 20 µL/well of cells to at least one well of a clear-bottom 384-well plate for image analysis of GFP expression.
- 6. Incubate plates for 20–24 hours in a humidified incubator at 37°C/5% CO₂.

Day 2. LanthaScreen® Cellular Assay

- 7. **Optional:** Twenty-four hours post-transduction, analyze GFP expression levels in the clear-bottom 384-well plate by fluorescence microscopy using standard FITC filter sets.
- During incubation, prepare Complete 6X Lysis Buffer by adding both protease inhibitor and phosphatase inhibitor cocktails to the provided 6X Lysis Buffer, at a 1:33 dilution of 100X stock (e.g., 30 μL of each 100X stock inhibitor per 1,000 μL of 6X Lysis Buffer) and add LanthaScreen[®] Tb-anti-ERK2 [pThr185/pTyr187] Antibody to 12 nM Mix gently by inversion. Store on ice until use.
- 9. Proceed with the following agonist or antagonist assay setup.

Agonist Assay Setup:

- a. Prepare a stock solution of 0.3% DMSO in assay medium. **Note:** If you are using a solvent other than DMSO for the agonist, change the solvent used in the control wells accordingly. Be careful to keep the amount of solvent consistent in all wells.
- b. Prepare 3X control agonist in assay medium with 0.3% DMSO (e.g., if treating U-2 OS cells with EGF, prepare 600 ng/mL in Assay Media for a 3X concentration). Run a dose response curve to determine the EC₁₀₀ for your control agonist solution.
- c. Prepare 3X test compound in assay medium (if the test compound is dissolved in DMSO, make sure the DMSO concentration for the 3X solution is 0.3%).

- d. Add 10 μL of assay medium with 0.3% DMSO to each unstimulated control and cell-free control well.
- e. Add 10 µL of the 3X control agonist (EGF) in assay medium to each stimulated control well.
- f. Add 10 µL of the 3X test compound in assay medium to each test compound well.
- g. Proceed to the next numbered step.

Antagonist Assay Setup:

- a. Prepare a stock solution of 0.5% DMSO in assay medium. **Note:** If you are using a solvent other than DMSO for the agonist, change the solvent used in the control wells accordingly. Be careful to keep the amount of solvent consistent in all wells.
- b. Prepare a 5X control compound Iressa in assay medium (e.g., prepare a 10 mM stock solution of Iressa in DMSO, then dilute to 2.5 μ M in Assay Media for a 5X concentration). Run a dose response curve to determine the IC₁₀₀ for your control inhibitor solution.
- c. Prepare a 5X test compound in assay medium (if the test compound is dissolved in DMSO, make sure the DMSO concentration for the 5X solution is 0.5%).
- d. Add 5 µL of the assay medium with 0.5% DMSO to each unstimulated, stimulated and cell-free control well.
- e. Add 5 µL of the 5X control inhibitor Iressa in assay medium to each control compound well.
- f. Add 5 µL of the 5X test compound in assay medium to each test compound well.
- g. Incubate the plate in a humidified $37^{\circ}C/5\%$ CO₂ incubator for 30–60 minutes.
- h. Prepare a 6X stock of agonist (EGF) in assay medium (e.g., if using U-2 OS cells, prepare 600 ng/mL in Assay Media for a 6X concentration). Run a dose response curve to determine the EC₈₀ for your agonist solution.
- i. Add 5 µL of the assay medium to each unstimulated and cell-free control well.
- j. Add 5 µL of 6X stock of agonist (EGF) to each stimulated, control and test compound well
- 10. Cover the plate and stimulate the cells by incubating in a humidified $37^{\circ}C/5\%$ CO₂ incubator for 6 minutes.
- *Tip:* The stimulation time must be optimized for each cell type and agonist. "Stimulation time" refers to the time from agonist addition to the addition of Complete 6X Lysis Buffer.
- 11. After stimulation, immediately add 6 µL of Complete 6X Lysis Buffer to each well. Cover the plate.
- 12. Incubate the covered plate at room temperature in the dark for 3 hours or other desired equilibration time. The equilibration time can be optimized for your cell line of interest.
- *Note:* Assay plates may be stored at 4°C overnight prior to reading. Let the plate warm to room temperature prior to reading.
- 13. Proceed to LanthaScreen[®] Detection.

LanthaScreen® Detection

Instruments and Filters

Detection can be performed on a variety of plate readers, including the PE Envision. The data presented on page 4 were generated using a BMG PHERAstar plate reader using the LanthaScreen[®] filter block available from BMG.

Visit <u>www.invitrogen.com/instrumentsetup</u> or contact Invitrogen Discovery Sciences technical support (<u>drugdiscoverytech@invitrogen.com</u> or 760-603-7200 (extension 40266)) for more information on performing LanthaScreen[®] Cellular Assays on your particular instrument.

Note: We do not recommend using monochromator-based instruments without adjustable bandwidth, as the sensitivity of these instruments is not sufficient to adequately detect the TR-FRET signal.

Reading the Assay Plate and Data Analysis

All measurements should be taken at room temperature from the top of the wells.

- 1. Let the assay plate warm to room temperature before reading, if necessary.
- 2. Set the fluorescence plate reader to top/time-resolved read mode.
- 3. Allow the lamp in the plate reader to warm up for at least 10 minutes before making measurements.
- 4. Use the filter selections described below. To obtain an assay window, filter bandwidths are critical and cannot be approximated.

	Settings to Measure Donor (Terbium) Signal	Settings to Measure Acceptor (TR-FRET to GFP) Signal		
Excitation filter:	337 nm (30 nm bandwidth)			
Emission filter:	490 nm (10 nm bandwidth)	520 nm (25 nm bandwidth)		
Dichroic Mirror:	Variable			
Delay Time:	100 µs			
Integration Time:	200 µs			

- 5. Calculate the acceptor/donor Emission Ratio (TR-FRET Ratio, 520 nm/490 nm) for each well, by dividing the acceptor emission values by the donor emission values.
- 6. **Optional:** Convert the data to a response ratio by dividing each emission ratio value by the value from unstimulated cells (cells not receiving agonist).

Appendix

Alternative Transduction Protocol (Tested for NIH3T3 and HCT116 cells)

In this protocol, cells are incubated with virus at the time of adhering to the tissue-culture plate, usually 24 hours prior to re-plating onto the assay plate.

Day 1

- 1. Begin with cell cultures grown to complete confluence under normal growth conditions. Confluence of cells may impact results, such as the assay window.
- 2. Trypsinize to harvest adherent cells as recommended by the cell-line manufacturer.

Invitrogen • BacMam ERK2 [pThr185/pTyr187] Cellular Assay User Guide Page 15 of 18

- Resuspend cells in growth medium. Seed cells such that the monolayers will be approximately 50– 80% confluent once attached and spread. Avoid plating the cells such that 80% confluence is exceeded 24 hours post-transduction.
- *Note:* For first time users, it may be useful to work with 6-well plates to minimize reagent consumption during the transduction optimization process.
- *Tip:* For most cell lines (with a doubling time of approximately 24 hours), a seeding density of $2-4 \times 10^4$ cells/cm² is optimal for BacMam transduction. It may be desirable to optimize the cell density for specific cell backgrounds.
- 4. Immediately after seeding the cells, add the desired amount of BacMam reagent to the cells. For initial optimization, we recommend testing 30%, 10%, 3%, and 1% v/v dilutions of BacMam reagent.
- 5. Add BacMam Enhancer Solution at a 1X final concentration if required. **Note:** For untested cell backgrounds, we recommend performing the transduction in the presence and absence of Enhancer Solution and then analyzing the expression of the BacMam target.
- 6. Place cells in a humidified $37^{\circ}C/5\%$ CO₂ incubator for 16–24 hours to allow for the transduction and expression of the GFP fusion protein.

Day 2

- 7. **Optional:** Analyze GFP expression levels by fluorescence microscopy using standard FITC filter sets.
- 8. Harvest the transduced cells and be careful not to over-trypsinize the cells as this can result in poor viability and a decreased assay window.
- 9. Resuspend the harvested cells in growth medium with serum to inactivate the trypsin. Centrifuge the cells at 200 × g for 5 minutes. Aspirate the growth medium and resuspend the cell pellet in assay medium (usually low serum).
- 10. Centrifuge the cells at $200 \times g$ for 5 minutes. Aspirate the assay medium and resuspend the cell pellet in assay medium at the desired density.
- *Tip:* The number of cells per well will affect the assay window and can be optimized. We recommend starting with 10,000–20,000 cells per well seeded in 20 μ L of assay medium. Therefore, resuspend cells to 0.5–1 × 10⁶ cells/mL.
- 11. Plate 20 μ L of transduced cells in assay medium into white tissue culture-treated 384-well plates and incubate the plates in a 37°C/5% CO₂ incubator (or appropriate) for 16–20 hours.
- *Note:* Serum starvation in low-serum or serum-free media is required for most cell types analyzed with this assay. We recommend starting with a 16–20 hour serum starvation and optimizing as needed.

Day 3

12. Proceed to the LanthaScreen[®] Cellular Assay on page 10 or 12.

Troubleshooting Guide

Observation	Potential Solutions
Weak/no expression of GFP-	Confirm that your fluorescence microscope is configured appropriately for
fusion in the cell line of interest in a clear-bottom 384- well plates.	detection of GFP/FITC.
	Perform a virus titration to find the optimal virus concentration for your cell background.
	Confirm that no contamination of the BacMam reagent has occurred.
	For first-time users, we recommend the standard transduction protocol using U-2 OS cells at 100% confluence. Overly confluent or unhealthy cells will not transduce efficiently.
	If the standard protocol works for U-2 OS cells but not for your cells, please try Alternative Transduction Protocol in the Appendix.
>50% expression of GFP-fusion is observed, but weak/no detectable stimulation of the posttranslational modification of the GFP-fusion in cell line of interest	Confirm that the fluorescence plate reader is configured appropriately for LanthaScreen [®] detection. Filter bandwidth requirements are exact. For more information about your specific instrument and to purchase filters, visit <u>www.invitrogen.com/instrumentsetup</u> . Contact Drug Discovery Technical Support for more information.
	Be sure that the cells are grown to the highest density at which they remain healthy before proceeding to the transduction step, to enhance the sensitivity of the cells to stimulation.
	For first-time users, we recommend following the standard transduction protocol using U-2 OS cells.
	Perform a stimulation time course experiment to find out the optimal stimulation time for your agonist and cell line of interest. (The stimulation time refers to the time from agonist addition to lysis).
	Image the cells in clear-bottom microtiter plates. Ensure that cells are adhered to the bottom of the plate and are not expressing very high levels of GFP. Dimly green cells are desirable. Excessive expression of the GFP-fusion may be deleterious to cell health.
> 60% expression of GFP-fusion, but very low assay window due to high background (stimulated and unstimulated cells generate high emission ratios compared to cell-free/Tb-antibody alone).	Image the cells in clear-bottom microtiter plates. Ensure that cells are adhered to the bottom of the plate and are not expressing very high levels of GFP. Dimly green cells are desirable. Excessive expression of the GFP-fusion may be deleterious to cell health.
Following transduction with the BacMam reagent, cells have a rounded/unspread phenotype and appear to be in poor health.	Under serum-starvation conditions, many cell types will appear rounded when imaged 24 or 48 hours post-transduction. This is common among cell types such as HEK293 and CHO. Under these conditions, cells are capable of modifying the GFP-fusion target, and should remain adhered despite poor cell spreading.
	Cells that appear rounded and detached from the clear-bottom microtiter plate may be unsuitable for GFP-fusion activation. Ensure that cells are not expressing excessive GFP-fusion as indicated above.
Day-to-day fluctuations in assay window are observed.	Be sure to use cells with the same growth conditions (e.g., same harvest density).

Purchaser Notification

Limited Use Label License No. 21: Bac-to-Bac® and Bac-to-Bac® HT

This product is sold under patent license from Monsanto for research purposes only and no license for commercial use is included. Requests for licenses for commercial manufacture or use should be directed to Director, Monsanto Corporate Research, 800 N. Lindbergh, St. Louis, Missouri 63167.

Limited Use Label License No. 127: GFP with Heterologous Promoter

This product is sold under license from Columbia University. Rights to use this product are limited to research use only, and expressly exclude the right to manufacture, use, sell or lease this product for use for measuring the level of toxicity for chemical agents and environmental samples in cells and transgenic animals. No other rights are conveyed. Not for human use or use in diagnostic or therapeutic procedures. Inquiry into the availability of a license to broader rights or the use of this product for commercial purposes should be directed to Columbia Innovation Enterprise, Columbia University, Engineering Terrace-Suite 363, New York, New York 10027.

Limited Use Label License No. 168: Lanthanide Chelates

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, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Limited Use Label License No. 198: Fluorescent Proteins and Stable Cell Lines Expressing Such Proteins

The purchase of this product conveys to the buyer the nontransferable 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). No rights are conveyed to modify or clone the gene encoding GFP contained in this product. The buyer cannot sell or otherwise transfer (a) this product, (b) its components, or (c) materials made by the employment of this product or its components to a third party or otherwise use this product or its components or materials made by the employment of this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the employment 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. Life Technologies 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 none of this product, or 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 Corporation is willing to accept return of the product with a full refund. For information on purchasing a license to use this product for purposes other than those permitted above, contact Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200 or outlicensing@lifetech.com.

Limited Use Label License No. 272: Humanized GFP

This product is sold for research use only. Not for therapeutic or diagnostic use in humans.

Limited Use Label License No. 306: Baculovirus Vectors

This product is for research use only by those researchers in laboratories of academic, government, industrial and/or clinical institutions engaged in the investigation of biological or biochemical processes, or research and development of biological products. This product is not to be used in the manufacture, use or sale of human or animal diagnostic, therapeutic or prophylactic products.

Limited Use Label License No. 308: WPRE Element

This product contains the Woodchuck Post-transcriptional Regulatory Element ("WPRE") which is the subject of intellectual property owned by The Salk Institute for Biological Studies, and licensed to Life Technologies 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) 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; and/or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. In addition, any use of WPRE outside of this product or the product's authorized use requires a separate license from the Salk Institute. Life Technologies will not assert a claim against the buyer of infringement of patents owned by Life Technologies and claiming 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 or for a Commercial Purpose. 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 on purchasing a license to this product for purposes other than research, contact Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008, Phone (760) 603-7200. Fax (760) 602-6500, or The Salk Institute for Biological Studies, 10010 North Torrey Pines Road, La Jolla, CA 92037, Attn.: Office of Technology Management, Phone: (858) 453-4100 extension 1275, Fax: (858) 546-8093.

Limited Use Label License No. 332: BacMam Virus Use

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 solely in accordance with the accompanying product literature or manual. Purchase of this product does not convey a license to expand, amplify, or otherwise propagate the provided viral particles or to otherwise modify or alter the virus by any means.

Limited Use Label License No. 358: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact <u>outlicensing@lifetech.com</u> or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

Trademarks

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

ATCC and HTB-96 are trademarks of ATCC.

© 2011 Life Technologies Corporation. All rights reserved. Reproduction forbidden without permission.