

# LanthaScreen<sup>®</sup> LRRK2 [pSer935] Cellular Assay Kit

Catalog no. A14685 or A14686 Protocol part no. A14685PIS Rev. date: 14 June 2012

Storage: Varies (see Kit Contents)Pub. no. MAN0006945Shipping: Varies

## **Kit Contents and Handling**

Component	Catalog no.	Amount	Storage	Handling
BacMam LRRK2-GFP Reagent (WT or G2019S)	A14687 or A14688	$5 \times 1 \text{ mL}$	4°C	DO NOT FREEZE Use sterile technique Protect from light
Tb-anti-LRRK2 pSer935 Antibody	A14689	5 µg	-20°C	Protect from light
LanthaScreen <sup>®</sup> 6X Cellular Assay Lysis Buffer	A12891	6 mL	4°C	On day of assay, add protease inhibitor*, phosphatase inhibitor*, and antibody
Instrument Control Terbium TR-FRET Kit	A14138	1 kit	4°C	Protect from light

\*See Materials Required but Not Provided on page 2.

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

## **Overview**

The LanthaScreen<sup>®</sup> LRRK2 [pSer935] Cellular Assay Kit combines the flexibility of BacMam gene delivery system with the robustness and power of the LanthaScreen<sup>®</sup> TR-FRET technology. It allows researchers to interrogate phosphorylation at serine 935 on Leucine-Rich Repeat Kinase-2 (LRRK2) protein in a variety of cell backgrounds to identify LRRK2 kinase inhibitors in a high throughput format. This kit provides enough material to perform the assay in a 384-well plate; larger sizes of all reagents are available through our catalog.

Figure 1 Assay Workflow. Day 1: Cells are transduced with BacMam reagent. Day 2: Cells are harvested and transferred to assay plate. Day 3: Following the treatment with or without LRRK2 inhibitor, cells are lysed in the presence of a Tb-labeled anti-LRRK2 pSer935 antibody and the levels of phosphorylation are measured using a Tb TR-FRET-compatible plate reader.

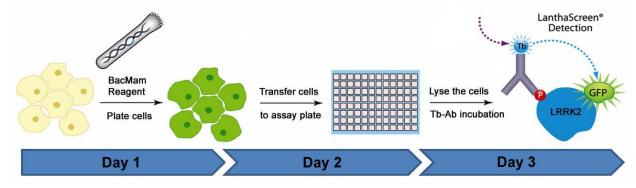
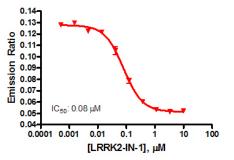


Figure 2 Representative data collected using U-2 OS cells transduced with BacMam LRRK2 G2019S.



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

## **Before Starting**

### Materials Required but Not Provided

Materials	<b>Recommended Source</b>	Part no.			
Fluorescence plate reader with Terbium (Tb) TR-FRET (LanthaScreen®) capability.					
Prior to setting up an assay, we strongly recommend checking your plate reader setup for Tb- based TR-FRET (LanthaScreen <sup>®</sup> ) detection using the Instrument Controls included with this kit.	For instrument set-up guidance, refer to www.lifetechnologies.com/instrumentsetup. For technical assistance, contact Drug Discovery Technical Support at drugdiscoverytech@invitrogen.com or 760-603-7200, extension 40266).				
Tb-based TR-FRET (LanthaScreen <sup>®</sup> ) technology requires specific instrument settings that are critical to experimental success. Also note that for this assay we do not recommend using monochromator-based instruments as the sensitivity of these instruments is generally not sufficient to detect the TR-FRET signal.					
<b>Cell Line of Interest</b> (visit <b>www.lifetechnologies.com/bacmam</b> for more information regarding BacMam compatible cell types; we have used this assay successfully with U-2 OS, HEK293T and SH-SY5Y cells)	Various	Various			
BacMam Enhancer Solution (1000X) (Do not use this solution if using U-2 OS cells)	Life Technologies	PV5835			
Protease Inhibitor Cocktail	Sigma	P8340			
Phosphatase Inhibitor Cocktail	Sigma	P0044			
Assay Plate					
• for TR-FRET, use white tissue culture-treated 384-well plates	Corning	3570			
• for visualization of LRRK2-GFP, use clear-bottom, tissue culture-treated plates	Corning	3712			

## **Assay Protocol**

## Day 1: BacMam LRRK2-GFP Transduction

The first critical experiment is a titration of the BacMam virus in your cell line of interest. For easy-to-transduce cells, such as U-2 OS, HEK293, HeLa, and human mammary epithelial cells, we recommend following the standard transduction protocol. For more information, also see the protocol described for BacMam GFP control (Cat. no. B10383) available at www.lifetechnologies.com.

*Note:* Expression levels of GFP Control and LRRK2-GFP do not correlate; therefore, optimal concentration of the GFP control virus may not apply to the BacMam LRRK2-GFP virus. To determine the optimal concentration of BacMam LRRK2-GFP reagents for your cell line of interest, you must perform a virus titration experiment.

We recommend testing a range of BacMam reagent dilutions (v/v) to determine the optimal percentage of virus for your transduction in the presence or absence of 0.5X BacMam Enhancer Solution (Cat. no. PV5835). As a starting point, we recommend using 50%, 30%, 20%, 10%, and 2% (v/v) of BacMam reagent for initial optimization.

1. Harvest and resuspend cells in growth medium. Add BacMam LRRK2-GFP to cell suspension at the optimal concentration. For U-2 OS we recommend 20% (v/v). Mix gently by inversion. (For some cell types except U-2 OS, 0.5X BacMam Enhancer solution can be added at this step to increase expression level.) Transfer cells/BacMam reagent mixture to appropriate cell-culture plates (such as 6-well plates). Incubate plates in a 37°C incubator with a humidified atmosphere of 5% CO<sub>2</sub> for 20–24 hours.

# For SH-SY5Y cells and other difficult-to-transduce cells, such as human astrocytes and primary neurons, we recommend the following transduction protocol.

- Start with SH-SY5Y cells already plated in a 6-well plate at 1.5 to 2 million cells/well in growth medium. Remove growth medium and wash cells once with PBS with Ca<sup>2+</sup> and Mg<sup>2+</sup> (Cat. no. 14040117). Add 1.5 mL of PBS with Ca<sup>2+</sup> and Mg<sup>2+</sup> to each well, and then add 0.5 mL of BacMam LRRK2-GFP reagent to 25% (v/v) final. Incubate the plate at room temperature in the dark with gentle rocking for 3–4 hours. Remove virus/PBS solution and add 2 mL/well of growth medium containing 0.3X BacMam Enhancer Solution. Incubate plate in a 37°C incubator with a humidified atmosphere of 5% CO<sub>2</sub> for 20–24 hours.
- *Note*: Growth medium for SH-SY5Y consists of DMEM/F12 (Cat. no. 10565) containing 10% dialyzed FBS (Cat. no. 26400) and 1X penicillin/streptomycin (Cat. no. 15140).

## Day 2: Cell Plating onto a 384-well assay Plate

1. After transduction, harvest and resuspended cells in Assay Medium at the following density (U-2 OS:  $0.5 \times 10^6$  cells/mL; HEK293:  $0.75 \times 10^6$  cells/mL; SH-SY5Y:  $1 \times 10^6$  cells/mL). Plate cells onto a 384-well assay plate (Corning, Cat. no. 3570) in 20 µL/well volume.

Optional: Plate cells into wells of a clear bottom plate (Corning, Cat. no. 3712) for GFP visualization.

- *Note:* Assay Medium consists of Opti-MEM I (Cat. no. 11058021) with 0.1% dialyzed FBS (Cat. no. 26400), 0.1 mM NEAA (Cat. no. 11140-050, 1 mM Sodium Pyruvate (Cat. no. 11360-070), and 100 U/mL Penicillin/100 μg/mL Streptomycin (Cat. no. 15140-122)
- 2. Centrifuge plate briefly at  $300 \times g$ . Incubate plates in a 37°C incubator with a humidified atmosphere of 5% CO<sub>2</sub> for 20–24 hours.

## Day 3: Inhibitor Treatment, Cell Lysis and TR-FRET Detection

- 1. Add 5  $\mu$ L/well of compounds or DMSO at 5X concentration in Assay Medium. Centrifuge plate briefly at 300 × g. Incubate plates in a 37°C incubator with a humidified atmosphere of 5% CO<sub>2</sub> for 60–90 minutes.
- 2. Prepare Complete 6X Lysis Buffer by adding protease inhibitor (Sigma, Cat. no. P8340) and phosphatase inhibitor (Sigma, Cat. no. P0044) cocktails to LanthaScreen<sup>®</sup> 6X Lysis Buffer (Cat. no. A12891), at a 1:33 dilution of 100X stock (e.g., 30 µL of 100X stock inhibitors per 1,000 µL of LanthaScreen<sup>®</sup> 6X Lysis Buffer) and add LanthaScreen<sup>®</sup> Tb-anti-LRRK2 [pSer935] Antibody to the 6X Lysis Buffer at a concentration of 12 nM. Mix gently by inversion. Store on ice until use.
- 3. Add 5  $\mu$ L/well of Complete LanthaScreen<sup>®</sup> 6X Lysis Buffer. Centrifuge plate briefly at 300 × g. Incubate at room temperature in the dark for 2–3 hours.
- 4. Read plate using instrument capable of detecting TR-FRET. See below for details on terbium TR-FRET detection or call our technical support team for assistance.

## Testing Terbium TR-FRET Detection Using the Instrument Controls

To test your instrument set-up for performing terbium-based TR-FRET cellular assays:

- 1. Add 60 µl/well of HIGH Control to empty assay plate wells for 96-well format (or 20 µL/well for 384-well format) with a minimum of 3 replicates.
- 2. Add 60 µl/well of the LOW Control to empty assay plate wells for 96-well format (or 20 µL/well for 384-well format) with a minimum of 3 replicates.
- 3. Read plate as described in the next section.
- 4. For each well, calculate the TR-FRET Emission Ratio (e.g., 520 nm/495 nm) by dividing the acceptor emission value (e.g., 520 nm) by the donor emission value (e.g., 495 nm).
- 5. Average the Emission Ratios for the HIGH Control and separately average the Emission Ratios for the LOW Control.
- 6. Determine the HIGH to LOW fold-change by dividing the average Emission Ratio for the HIGH by the average Emission Ratio for the LOW.
- *Note:* The HIGH/LOW fold-change should be 2–4, depending on the plate reader used. Values below 2 may indicate the instrument is not setup properly and/or lacks sufficient sensitivity for Tb-based TR-FRET.

## **Reading the Assay Plate**

- 1. Let assay plate warm to room temperature before reading, if necessary.
- 2. Set fluorescence plate reader to top/time-resolved read mode (allow plate reader lamp to warm up for at least 10 minutes before making measurements).
- 3. Remove lid/seal and read plate using the LanthaScreen<sup>®</sup> Tb TR-FRET instrument-specific filter selection guidelines provided at www.lifetechnologies.com/instrumentsetup.
- *Note:* Filter bandwidths are critical and cannot be approximated. For instruments using a flash lamp light source, we strongly recommend using 200 flashes for the acceptor channel (e.g., 520 nm) and 100 flashes for the terbium channel (e.g., 495 nm). We also recommend performing plate height (or focal height) optimization for best results.

### Data Analysis

- 1. For each well, calculate the TR-FRET Emission Ratio (e.g., 520 nm/495 nm) by dividing the acceptor emission value (e.g., 520 nm) obtained for that well by the donor emission value (e.g., 495 nm) obtained for that well. Do **not** average the 520 nm or 495 nm readings and then take the ratio.
- *Note:* A common practice employed with TR-FRET data is to multiply the Emission Ratio values by 10,000. This is done to convert Emission Ratio values from decimals to integers for easier graphing and data visualization.

In general, we do not recommend performing background subtraction with TR-FRET data since care must be taken to not inadvertently distort the results. However, in certain cases it may be desirable to perform background subtraction by subtracting the Emission Ratio value obtained for the cell-free control from the Emission Ratios of untreated and treated cell wells.

2. *Optional*: Convert TR-FRET Emission Ratio data into normalized Assay Window values, by dividing each Emission Ratio value determined for an untreated (or DMSO control) cells by the value calculated from the treated cells.

## Additional Assay Performance Data and Troubleshooting Guide

Due to the variety of cell types in which this product may be used, further assay optimization including titration of BacMam Reagent, cell plating density, DMSO tolerance, test compound concentrations and treatment times may be required and is recommended for getting best results in your cell-background of interest. For additional assay performance data and troubleshooting guide, visit **www.lifetechnologies.com** and search for A14687 or A14688 to download the **Application Note** for this assay or call the technical support hotline. Protocol and Application Notes are located under the "How to Use" tab on the product page.

## **Limited Product Warranty**

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

### Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

## **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 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 above patents based upon the manufacture, use, or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by 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 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. 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 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,

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

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 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 or for a Commercial Purpose. If the purchaser is not willing to accept the limitations of this limited use statement, 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 outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

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

© 2012 Life Technologies Corporation. All rights reserved. Reproduction forbidden without permission.