



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data sheet
LT β R:LIGHT[Biotinylated] Inhibitor Screening Assay Kit
Catalog #79808
Size: 96 reactions

BACKGROUND: Lymphotoxin β receptor (LT β R) and herpes virus entry mediator (HVEM) are the physiological receptors of LIGHT (tumor necrosis factor superfamily member 14 (TNFSF14) or CD258). The activation of the LT β R/HVEM:LIGHT co-signaling pathways induces the production of pro-inflammatory molecules, like NF- κ B, chemokines and adhesion molecules. LT β R is up regulated in hepatitis and hepatocellular carcinoma, and influences lipid metabolism and atherosclerosis. It also regulates cell growth and can initiate inflammation related carcinogenesis.

DESCRIPTION: The *LT β R:LIGHT[Biotinylated] Inhibitor Screening Assay Kit* is designed for screening and profiling inhibitors of LIGHT binding to LT β R. This kit comes in a convenient 96-well format, with biotin-labeled LIGHT, purified LT β R, streptavidin-labeled HRP, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of biotin-labeled LIGHT by streptavidin-HRP. Only a few simple steps on a microtiter plate are required for the assay. First, LT β R is coated on a 96-well plate. Next, biotinylated LIGHT is incubated with LT β R on the plate. Finally, the plate is treated with streptavidin-HRP followed by addition of an HRP substrate to produce chemiluminescence, which can be measured using a chemiluminescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
71112	Human LTBR, Fc fusion HiP™	10 μ g	-80°C	Avoid multiple freeze/thaw cycles!
100032	LIGHT, Biotin-labeled	10 μ g	-80°C	
79311	3x Immuno Buffer 1	50 ml	-20°C	
79728	Blocking Buffer 2	50 ml	+4°C	
79742	Streptavidin-HRP	15 μ l	+4°C	
79670	ELISA ECL Substrate A (transparent bottle)	6 ml	Room Temp	
	ELISA ECL Substrate B (brown bottle)	6 ml	Room Temp	
79699	96-well white microplate	1	+4°C	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

PBS (Phosphate buffered saline)
Luminometer or fluorescent microplate reader capable of reading chemiluminescence
Adjustable micropipettor and sterile tips

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

APPLICATIONS: This kit is useful for screening for inhibitors of LIGHT binding to LT β R.

STABILITY: Up to 6 months from date of receipt, when stored as recommended.

CONTRAINDICATIONS: Keep final DMSO concentration at or below 1%. Excess DMSO interferes with binding and decreases sensitivity of the assay.

REFERENCES:

Chang, J. H., *et al. J Biomed Sci* 2005, **12(2)**: 363-375.
del Rio, M. L., *et al. MAbs* 2016, **8(3)**: 478-490.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Coating the plate with LT β R :

- 1) Thaw **LT β R** on ice. Upon first thaw, briefly spin tube containing **LT β R** to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining **LT β R** in aliquots at -80°C. Note: **LT β R** is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.
- 2) Dilute **LT β R** to 2 μ g/ml in PBS.
- 3) Add 50 μ l of diluted LT β R solution to each well and incubate overnight at 4°C. Leave a couple of wells empty (uncoated), for use with the "Ligand Control" (see below).
- 4) Dilute **3x Immuno Buffer 1** to **1x Immuno Buffer 1** with water. (Keep a portion of the **3x Immuno Buffer 1** undiluted, for use in Step 1 below).
- 5) Decant to remove supernatant. Wash the plate three times with 100 μ l **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 6) Block wells by adding 100 μ l of **Blocking Buffer 2** to each well. Incubate for 1 hour at room temperature with slow shaking. Remove supernatant as described in step 5.

Step 1:

- 1) Prepare the master mixture: N wells \times (10 μ l **3x Immuno Buffer 1** + 15 μ l distilled water)
- 2) Add 25 μ l of master mixture to each well. Use uncoated wells for the "Ligand Control".
- 3) Prepare 10X concentration of test inhibitor in an aqueous-based solution. If the test inhibitor is dissolved in DMSO, make sure the final DMSO concentration in the assay is \leq 1%. (For example, to test an inhibitor at 10 μ M final concentration, prepare the inhibitor at 1 mM in 100% DMSO. Then dilute 1 mM inhibitor in water to 100 μ M, which contains 10% of DMSO. Use 5 μ l of the 100 μ M inhibitor solution (10% DMSO) in the assay to make a 1% DMSO concentration in the final 50 μ l reaction mixture.)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

- 4) For the inhibitor buffer, prepare the same solution as above, but without the test inhibitor (e.g. 10% DMSO in water). The DMSO concentration should be the same as in the 10X inhibitor solution above.
- 5) Add 5 µl of inhibitor solution to each well designated “Test Inhibitor”. For the “Positive Control”, “Ligand Control” and “Blank”, add 5 µl of the same solution without inhibitor (inhibitor buffer). Incubate at room temperature for one hour with slow shaking.
- 6) Thaw **LIGHT-biotin** on ice. Upon first thaw, briefly spin tube containing the protein to recover full contents of the tube. Aliquot **LIGHT-biotin** into single use aliquots. Immediately store remaining undiluted protein in aliquots at - 80 °C. Note: **LIGHT-biotin** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 7) Dilute **LIGHT-biotin** to 3 ng/µl (approximately 50 nM) in **1x Immuno Buffer 1**. Keep diluted protein on ice until use. Discard any unused diluted protein after use.
- 8) Add 20 µl of **1x Immuno Buffer 1** to the well designated “Blank”.

	Blank	Ligand Control	Positive Control	Test Inhibitor
3x Immuno Buffer 1	10 µl	10 µl	10 µl	10 µl
Distilled water	15 µl	15 µl	15 µl	15 µl
Test Inhibitor	-	-	-	5 µl
Inhibitor buffer (no inhibitor)	5 µl	5 µl	5 µl	-
1x Immuno Buffer 1	20 µl	-	-	-
LIGHT-biotin (3 ng/µl)	-	20 µl	20 µl	20 µl
Total	50 µl	50 µl	50 µl	50 µl

- 9) Initiate reaction by adding 20 µl of diluted **LIGHT-biotin** (see Step 1-7) to wells labeled “Positive Control”, “Ligand Control” and “Test Inhibitor”. Incubate at room temperature for one hour with slow shaking.
- 10) Decant to remove supernatant. Wash the plate 3 times with 100 µl/well **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 11) Block wells by adding 100 µl of **Blocking Buffer 2** to each well. Incubate for 10 minutes at room temperature. Remove supernatant as in Step 1-10.

Step 2:

- 1) Dilute **Streptavidin-HRP** 1000-fold with **Blocking Buffer 2**.
- 2) Add 100 µl to each well. Incubate for 1 hour at room temperature with slow shaking.
- 3) Wash plate three times with **1x Immuno Buffer 1**. Tap plate onto clean paper towel to remove liquid.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

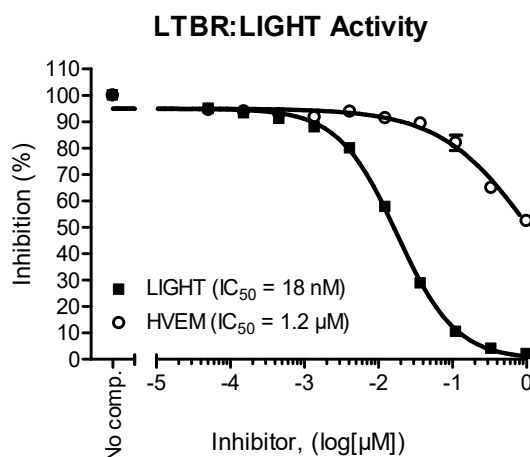
- 4) Block wells by adding 100 μ l of **Blocking Buffer 2** to each well. Incubate for 10 minutes at room temperature. Decant to remove supernatant. Tap plate onto clean paper towels to remove liquid.
- 5) Just before use, mix on ice 50 μ l **ELISA ECL Substrate A** and 50 μ l **ELISA ECL Substrate B**, then add 100 μ l to each well. Discard any unused chemiluminescent reagent after use.
- 6) Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence. "Blank" value is subtracted from all readings.

Reading Chemiluminescence:

Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry.

To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second; delay after plate movement is 100 milliseconds. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the "hole" position on the filter wheel; Optics position: Top; Read type: endpoint. Sensitivity may be adjusted based on the luminescence of a control assay without enzyme (typically we set this value as 100).

Example of Assay Results:



LT β R:LIGHT binding activity, measured using the using the *LT β R:LIGHT[Biotinylated] Inhibitor Screening Assay Kit*, BPS Bioscience #79808. Inhibition of LT β R:LIGHT binding using unlabeled LIGHT (closed squares, BPS Bioscience #71266) and HVEM (open circles, BPS Bioscience #71142) in the *LT β R:LIGHT[Biotinylated] Screening Assay Kit*. Luminescence was measured using a Bio-Tek fluorescent microplate reader. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
 San Diego, CA 92121
 Tel: 1.858.202.1401
 Fax: 1.858.481.8694
 Email: info@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog#</u>	<u>Size</u>
Human LTBR, Fc fusion HiP™	71112	100 µg
Mouse LTBR, Fc fusion HiP™	71122	100 µg
LIGHT, His-Tag (Human)	71266	100 µg
LIGHT, His-Tag (Mouse)	79068	100 µg
LIGHT-CHO Recombinant Cell Line	79262	2 vials
LIGHT:HVEM [Biotinylated] Inhibitor Screening Assay Kit	79684	96 reactions

TROUBLESHOOTING GUIDE

Problem	Possible cause	Solution
Luminescence signal of positive control reaction is weak	LTβR or LIGHT-biotin has lost activity	Proteins lose activity upon repeated freeze/thaw cycles. Use fresh LIGHT-biotin, (#100032) and fresh LTβR (#71112). Store proteins in single-use aliquots. Increase time of protein incubation. Increase protein concentration.
	Incorrect settings on instruments	Refer to instrument instructions for settings to increase sensitivity of light detection.
	Chemiluminescent reagents mixed too soon	Chemiluminescent solution should be used within 15 minutes of mixing. Ensure both reagents are properly mixed.
Luminescent signal is erratic or varies widely among wells	Inaccurate pipetting/technique	Run duplicates of all reactions. Use a multichannel pipettor. Use master mixes to minimize errors.
	Bubbles in wells	Pipette slowly to avoid bubble formation. Tap plate lightly to disperse bubbles; be careful not to splash between wells.
Background (signal to noise ratio) is high	Insufficient washes	Increase number of washes. Increase wash volume. Increase Tween-20 concentration to 0.1% in PBST
	Sample solvent is inhibiting the enzyme	Run negative control assay including solvent. Maintain DMSO level at <1% Increase time of protein incubation.
	Results are outside the linear range of the assay	Use different concentrations of LIGHT-biotin (BPS Bioscience #100032) to create a standard curve

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com