



6405 Mira Mesa Blvd., Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

## **Data Sheet**

### **IL-8 Promoter Luciferase Reporter Lentivirus**

**Catalog #: 79827**

#### **Product Description**

The IL-8 Promoter Luciferase Reporter Lentivirus are replication incompetent, HIV-based, VSV-G pseudotyped lentiviral particles that are ready to be transduced into almost all types of mammalian cells, including primary and non-dividing cells. The particles contain a firefly luciferase gene driven by the human IL-8 promoter (Figure 1). After transduction, activation of the IL-8 signaling pathway in the target cells can be monitored by measuring the luciferase activity.

#### **Application**

- Screen for activators or inhibitors of IL-8 signaling pathway in transduced target cells
- Generation of IL-8 promoter luciferase reporter stable cell line

#### **Formulation**

The lentiviruses were produced from HEK293T cells in medium containing 90% DMEM + 10% FBS.

#### **Titer**

Two vials (500  $\mu$ l x 2) of IL-8 promoter luciferase reporter lentivirus at a titer  $>5 \times 10^6$  TU/ml. The titer will vary with each lot; the exact value is provided with each shipment.

#### **Storage**

Lentiviruses are shipped with dry ice. For long term storage, it is recommended to store the virus at  $-80^{\circ}\text{C}$ . Avoid repeated freeze-thaw cycles. Titers can drop significantly with each freeze-thaw cycle.

#### **Biosafety**

The lentiviruses are produced with the third generation SIN (self-inactivation) lentivector which ensures self-inactivation of the lentiviral construct after transduction and integration into the genomic DNA of the target cells. None of the HIV genes (gag, pol, rev) will be expressed in the transduced cells, as they are expressed from packaging plasmids lacking the packing signal.

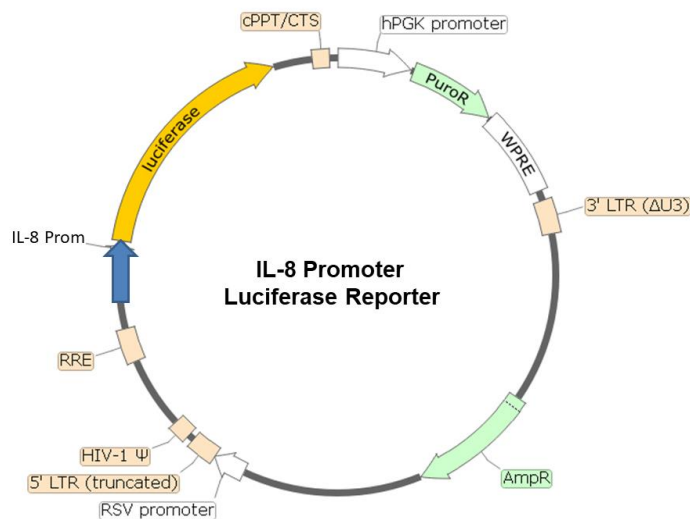
Although the pseudotyped lentiviruses are replication-incompetent, they require the use of a Biosafety Level 2 facility. BPS recommends following all local federal, state, and institutional regulations and using all appropriate safety precautions.

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



**Figure 1. Schematic of the lenti-vector used to generate the IL-8 promoter luciferase reporter lentivirus**

**Materials Required but Not Supplied**

- Jurkat cells (ATCC #TIB-152)
- Jurkat growth medium (Thaw Medium 2, BPS Bioscience, #60184)
- PMA (LC Laboratories, #P1680)
- Ionomycin (Sigma, #I3909)
- Polybrene (Millipore, #TR-1003-G)
- 96-well tissue culture treated white clear-bottom assay plate (Corning, #3610)
- One-Step luciferase assay system (BPS Bioscience, #60690)
- Luminometer

**Assay protocol**

The following protocol is a general guideline for transducing Jurkat cells using IL-8 promoter luciferase reporter lentivirus. The optimal transduction conditions (e.g. MOI, concentration of polybrene, time of assay development) should be optimized according to the cell type and the assay requirements. In most cell types, the expression of the reporter gene can be measured approximately 72 hours after transduction. For cell types with low transduction efficacy, it may be necessary to select the cells stably expressing the reporter gene with puromycin prior to carrying out the reporter assays.

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd., Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

1. Harvest the Jurkat cells by centrifugation and resuspend the cells in fresh growth medium. Dilute the cells to  $1 \times 10^6$  /ml in growth medium. Mix 500  $\mu$ l of the Jurkat cells and 100  $\mu$ l of IL-8 promoter luciferase reporter lentivirus in a 1.5-ml Eppendorf tube. Add polybrene to a final concentration of 8  $\mu$ g/ml. Gently mix and incubate the virus with the Jurkat cells for 20 minutes at room temperature in the tissue culture hood.
2. Centrifuge the virus/cells mixture for 30 minutes at 800 x g at 32°C. Remove the virus containing medium and resuspend the cell pellet in 2 ml of fresh Jurkat growth medium. Transfer the cells into one well in a 6-well plate. Incubate the plate at 37°C with 5% CO<sub>2</sub> for 48-66 hours. The transduced Jurkat cells are ready for assay development.
3. Harvest the cells and resuspend the cells into 1 ml of fresh Jurkat growth medium. Add 90  $\mu$ l of the cells to each well of the 96-well plate. Dilute ionomycin and PMA using Jurkat growth medium, and add 5  $\mu$ l of diluted ionomycin and 5  $\mu$ l of diluted PMA to the stimulated wells. Add 10  $\mu$ l of Jurkat growth medium to the unstimulated control wells (for measuring the uninduced level of IL-8 promoter reporter activity).
4. Incubate at 37°C with 5% CO<sub>2</sub> for 5-6 hours.
5. Prepare the ONE-Step™ Luciferase reagent per recommended protocol. Add 100  $\mu$ l of ONE-Step™ Luciferase Assay reagent per well. Incubate at room temperature for ~15 to 30 minutes and measure luminescence using a luminometer.

#### Important Notes:

1. To generate the IL-8 promoter luciferase reporter stable cell line, remove the growth medium 48 hours after transduction and replace it with fresh growth medium containing the appropriate amount of puromycin for antibiotic selection of transduced cells.
2. The following Lenti Reporter Controls are also available from BPS Bioscience to meet your experimental needs:
  - 1) Reporter Negative Control Lentivirus (BPS Bioscience, #79578): Ready-to-transduce lentiviral particles expressing firefly luciferase under the control of a minimal promoter. The negative control is important to establish the specificity of any treatments and to determine the background reporter activity.
  - 2) Renilla Luciferase (Rluc) Lentivirus (BPS Bioscience, #79565): Ready-to-transduce lentiviral particles expressing Renilla luciferase under the CMV promoter. The RLuc lentivirus can serve as an internal control to overcome sample-to-sample variability when performing dual-luciferase reporter assays.

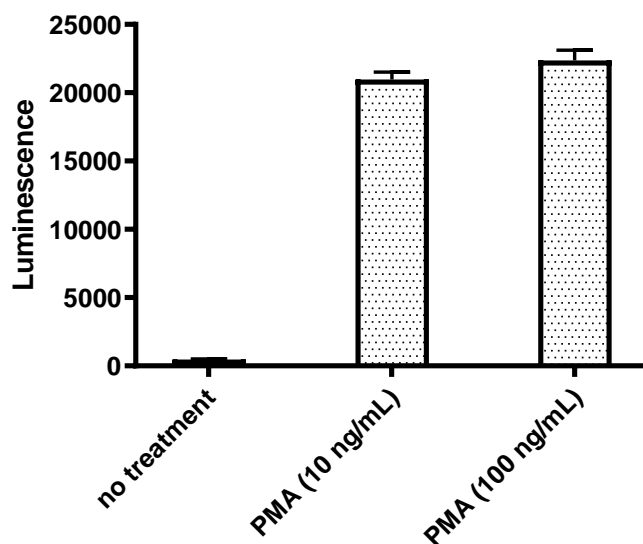
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

- 3) Firefly Luciferase (Fluc) Lentivirus (BPS Bioscience, #79692-G, #79692-H, #79692-P): Ready-to-transduce lentiviral particles expressing firefly luciferase under the CMV promoter. The Fluc lentivirus can serve as a positive control for transduction optimization studies.



**Figure 2. IL-8 promoter luciferase reporter activity stimulated by Ionomycin/PMA in Jurkat cells.** Appropriate 50,000 Jurkat cells were transduced with 100,000 TU IL-8 promoter luciferase reporter lentiviruses. After 48 hours of transduction, medium was changed to fresh Jurkat growth medium. Cells were stimulated with PMA in the presence of 0.5  $\mu$ M ionomycin for 5 hours. The results are shown as the raw luminescence reading.

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd., Ste 100

San Diego, CA 92121

**Tel:** 1.858.202.1401

**Fax:** 1.858.481.8694

**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## Related Products

<b><u>Product</u></b>	<b><u>Cat. #</u></b>	<b><u>Size</u></b>
NF- $\kappa$ B Luciferase Reporter Lentivirus	79564	500 $\mu$ l x2
CRE/CREB Luciferase Reporter Lentivirus	79580	500 $\mu$ l x2
NFAT Luciferase Reporter Lentivirus	79579	500 $\mu$ l x2
STAT3 Luciferase Reporter Lentivirus	79744	500 $\mu$ l x2
STAT5 Luciferase Reporter Lentivirus	79745	500 $\mu$ l x2
TCF/LEF Luciferase Reporter Lentivirus	79787	500 $\mu$ l x2
ISRE Luciferase Reporter Lentivirus	79824	500 $\mu$ l x2
IL-2 Promoter Luciferase Reporter Lentivirus	79825	500 $\mu$ l x2
Reporter Negative Control Lentivirus	79578	500 $\mu$ l x2
Renilla Luciferase (Rluc) Lentivirus	79565	500 $\mu$ l x2
Firefly Luciferase (Fluc) Lentivirus (G418)	79692-G	500 $\mu$ l x2
Firefly Luciferase (Fluc) Lentivirus (Hygromycin)	79692-H	500 $\mu$ l x2
Firefly Luciferase (Fluc) Lentivirus (Puromycin)	79692-P	500 $\mu$ l x2
GFP Reporter Lentivirus	79703	1 ml x 2
Anti-CD3 Agonist Antibody	71274-1	50 $\mu$ g
Anti-CD28 Agonist Antibody	100182	50 $\mu$ g
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml
Dual Luciferase (Firefly-Renilla) Assay System	60683	10 ml

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)