

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

TCF/LEF Luciferase Reporter Lentivirus (Wnt/β-catenin Signaling Pathway) Catalog #: 79787

Product Description

The TCF/LEF 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 under the control of TCF/LEF-responsive element located upstream of the minimal TATA promoter (Figure 1). After transduction, activation of the Wnt/β-catenin signaling pathway in the target cells can be monitored by measuring the luciferase activity.

Applications

- Screen for activators or inhibitors of Wnt signaling pathway in the transduced target cells
- Generation of TCF/LEF luciferase reporter stable cell line

Formulation

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

Titer

Two vials (500 μ I x 2) of TCF/LEF luciferase reporter lentivirus at a titer \geq 1 x 10⁷ TU/mI. 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°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.



6042 Cornerstone Court W, Ste B San Diego, CA 92121 **Tel:** 1.858.202.1401

Fax: 1.858.481.8694

Email: info@bpsbioscience.com

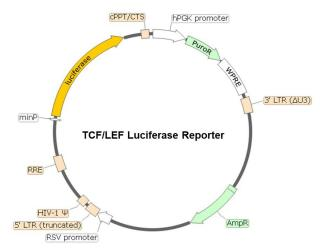


Figure 1. Schematic of the lenti-vector used to generate the TCF/LEF luciferase reporter lentivirus

Materials Required but Not Supplied

- LiCl (Sigma, #L7026)
- Mouse Wnt3a (R&D Systems 1324-WN)
- HEK293 growth medium or use Thaw Medium 9 (BPS Bioscience, #79665): MEM with 10% FBS, 0.1 mM nonessential amino acids, 1 mM sodium pyruvate
- 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 HEK293 cells using TCF/LEF 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.

1. Day 1: Harvest HEK293 cells from culture and seed cells at a density of 5,000-10,000 cells per well into white opaque 96-well microplate in 50 μl of HEK growth medium. Incubate cells at 37°C with 5% CO₂ overnight.



6042 Cornerstone Court W, Ste B San Diego, CA 92121 **Tel:** 1.858.202.1401

Fax: 1.858.481.8694
Email: info@bpsbioscience.com

2. Day 2: To each well add 10 μl of TCF/LEF luciferase reporter lentivirus. Add polybrene to each well at a final concentration of 5 μg/ml. Gently swirl the plate to mix. Incubate the plate at 37°C with 5% CO₂ for 18-24 hours.

Alternatively, seeding cells and the transduction can be performed at the same day.

- 3. Day 3: Remove the medium containing the lentivirus from the wells. Add 100 µl of fresh growth medium to each well.

 If neither the polybrene nor the lentivirus adversely affects the target cells, it is not necessary to change the medium on Day 3. The target cells can be incubated with the virus for 48-72 hours before changing the medium.
- 4. Day 4: Treat transduced cells with LiCl (final concentration 10 mM) in 90 μl fresh growth medium. Incubate the cells at 37°C with 5% CO₂ for 16 hours.
- 5. Day 5: Prepare diluted mouse Wnt3a in HEK growth medium. Add 10 μl diluted Wnt3a to stimulated cells; add 10 μl growth medium to unstimulated control cells; add 100 μl growth medium to cell-free control wells (for determining background luminescence). Set up each treatment in at least triplicate. Incubate at 37°C with 5% CO₂ for 5-6 hours.
- 6. Prepare the ONE-Step™ Luciferase reagent per recommended protocol. Add 100 µl 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 TCF/LEF luciferase reporter stable cell line, on day 4 remove HEK growth medium and replaced 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:
 - 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.
 - 3) Firefly Luciferase (Fluc) Lentivirus (BPS Bioscience, #79692-G, #79692-H): 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.



6042 Cornerstone Court W, Ste B San Diego, CA 92121

Tel: 1.858.202.1401 Fax: 1.858.481.8694 Email: info@bpsbioscience.com

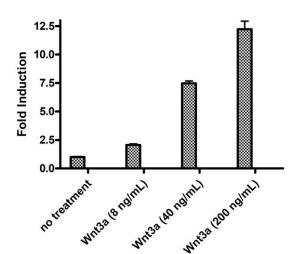


Figure 2. TCF/LEF luciferase reporter activity stimulated by mouse Wnt3a in HEK293 cells. Approximately 10,000 HEK293 cells/well were transduced with 100,000 TU/well TCF/LEF luciferase reporter lentivirus. After 48 hours of transduction, the cells were treated with LiCl (10 mM) for 16 hours. The cells were then treated with mouse Wnt3a for 5 hours. The results are shown as fold induction of luciferase reporter expression. Fold induction was determined by comparing values against the mean value for control cells without Wnt3a treatment.

Related Products

<u> </u>	
Cat. #	<u>Size</u>
79564	500 µl x2
79580	500 µl x2
79579	500 µl x2
79744	500 µl x2
79745	500 µl x2
79806	500 µl x2
79578	500 µl x2
79565	500 µl x2
79692-G	500 µl x2
79692-H	500 µl x2
60500	500 rxns
60501	2 vials
60690-1	10 ml
60690-2	100 ml
60683	10 ml
79665	100 ml
	79564 79580 79579 79744 79745 79806 79578 79565 79692-G 79692-H 60500 60501 60690-1 60690-2 60683

References

- 1. Tian S., et al., Blood. 1994; 84(6):1760-1764.
- 2. Zhong, Z., et al., Science. 1994; **264(5155):**95-98.

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