

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

Negative Control Luciferase Lentivirus Catalog #: 79578

Product Description

The Negative Control Lentivirus (Firefly Luciferase) are replication incompetent, HIV-based, VSV-G pseudo typed 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 a minimal TATA promoter, without any additional transcriptional response elements (Figure 1).

Application

Useful to determine the background reporter activity and establish the specificity of any treatment effects for the Lenti reporter system

Formulation

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

Titer

Two vials (500 μ l x 2) of negative control lentivirus at a titer 1 x 10⁷ 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°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

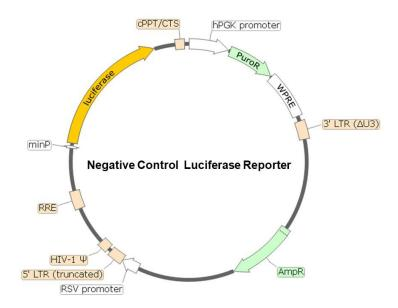


Figure 1. Schematic of the lenti-vector used to generate the negative control lentivirus

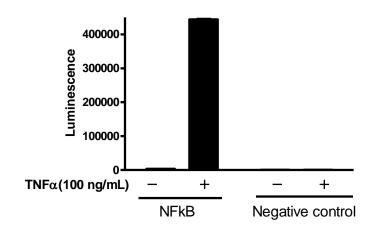


Figure 2. NF-κB luciferase reporter activity stimulated by TNF α in HEK293 cells. 10,000 HEK293 cells/well were transduced with 100,000 TU/well NF-κB luciferase reporter lentivirus or 100,000 TU/well negative control virus. After 48 hours of transduction, medium was changed to HEK growth medium. After 66 hours of transduction, cells were treated with 100 ng/mL of TNF α for ~6 hours. The results are shown as the raw luminescence reading.



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</u>	Cat. #	<u>Size</u>
NF-κB Luciferase Reporter Lentivirus	79564	500 µl x2
CRE Luciferase Reporter Lentivirus	79580	500 µl x2
NFAT Luciferase Reporter Lentivirus	79579	500 µl x2
STAT3 Luciferase Reporter Lentivirus	79744	500 µl x2
STAT5 Luciferase Reporter Lentivirus	79745	500 µl x2
TCF/LEF Luciferase Reporter Lentivirus	79787	500 µl x2
Negative Control Lentivirus	79578	500 µl x2
Firefly Luciferase (Fluc) Lentivirus (G418)	79692-G	500 µl x2
Firefly Luciferase (Fluc) Lentivirus (Hygromycin)	79692-H	500 µl x2
Firefly Luciferase (Fluc) Lentivirus (Puromycin)	79692-P	500 µl x2
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
· · · · · · · · · · · · · · · · · · ·		-

References

- Pessara, U., Koch, N. (1990) Tumor necrosis factor alpha regulates expression of the major histocompatibility complex class II-associated invariant chain by binding of an NFκB-like factor to a promoter element. *Mol Cell Biol.* 10(8):4146-4154.
- 2. Baeuerle, P.A. (1998) Pro-inflammatory signaling: last pieces in the NF-κB puzzle? *Curr Biol.* **8(1):**R19-R22.