

Data Sheet

Expression Negative Control Lentivirus (G418)

Catalog #: 79902-G

Product Description

The Expression Negative Control 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 controls package the same virus particles as the target expression virus, but they do not express a specific protein under the CMV promoter (Figure 1). The Expression Negative Control Lentivirus (G418) expresses the gene for aminoglycoside 3' phosphotransferase, which confers resistance to kanamycin, neomycin, and geneticin (G418).

Application

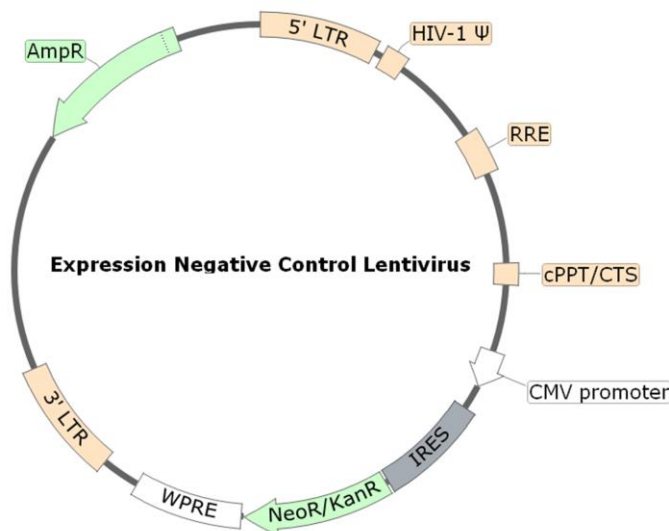
1. Useful as controls for the target expression lentivirus having G418 selection marker.
2. Generation of stable cell line with G418 selection without expressing a specific protein under the CMV promoter, to validate the specificity of any target expression effects.

Formulation

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

Titer

Two vials (500 μ l x 2) of Expression Negative Control Lentivirus at a titer 1×10^7 TU/ml. The titer will vary with each lot; the exact value is provided with each shipment.



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

Figure 1. Schematic of the lenti-vector used to generate the Expression Negative Control Lentivirus

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

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.

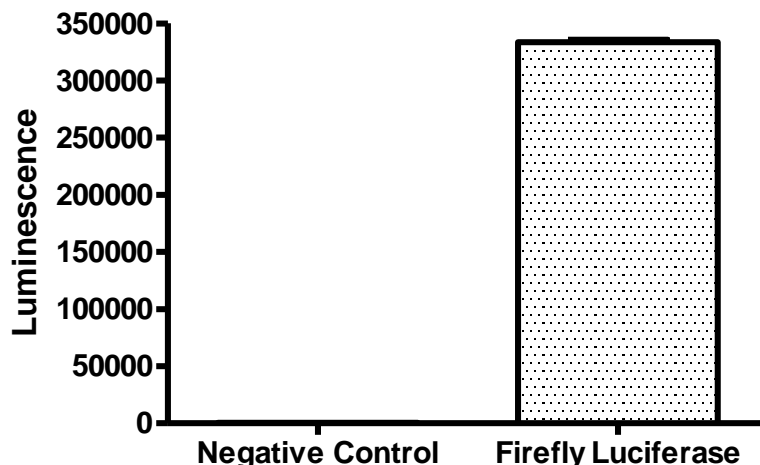


Figure 2. Luciferase activity in Jurkat cells transduced with firefly luciferase lentivirus.

Approximately 20,000 Jurkat cells/well were transduced with 200,000 TU/well firefly luciferase lentivirus (BPS Bioscience #79692-G) or expression negative control lentivirus (BPS Bioscience #79902-G) using spinoculation. After 66 hours of transduction, medium was changed to Thaw Medium 2 (BPS Bioscience #60184). The luciferase assay was performed using the ONE-Step™ Luciferase assay system (BPS Bioscience, #60690), following the recommended protocol in the user manual. 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: 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</u>	<u>Cat. #</u>	<u>Size</u>
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
NFκB 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
ISRE Luciferase Reporter Lentivirus	79824	500 µl x2
IL-2 Promoter Luciferase Reporter Lentivirus	79825	500 µl x2
IL-8 Promoter Luciferase Reporter Lentivirus	79827	500 µl x2
AP-1 Luciferase Reporter Lentivirus	79823	500 µl x2
SBE Luciferase Reporter Lentivirus	79806	500 µl x2
TEAD Luciferase Reporter Lentivirus	79833	500 µl x2
ARE Luciferase Reporter Lentivirus	79869	500 µl x2
Negative Control Lentivirus	79578	500 µl x2
Renilla Luciferase (Rluc) Lentivirus	79565	500 µl x2
FcGR11IA Lentivirus	79876	500 µl x2
FcGR11B Lentivirus	79877	500 µl x2
FcER1G Lentivirus	79878	500 µl x2
Expression negative Control Lentivirus	79902	500 µl x2
TCR Activator Lentivirus	79894	500 µl x2
Non-Secreted Gaussia Luciferase Lentivirus	79893	500 µl x2
Thaw Medium 2	60184	100 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: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com