

## Data Sheet

### **Expression Negative Control Lentivirus (Puromycin)**

Catalog #: 79902-P

#### **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 (Puromycin) expresses the gene for puromycin N-acetyl-transferase, which confers resistance to puromycin.

#### **Application**

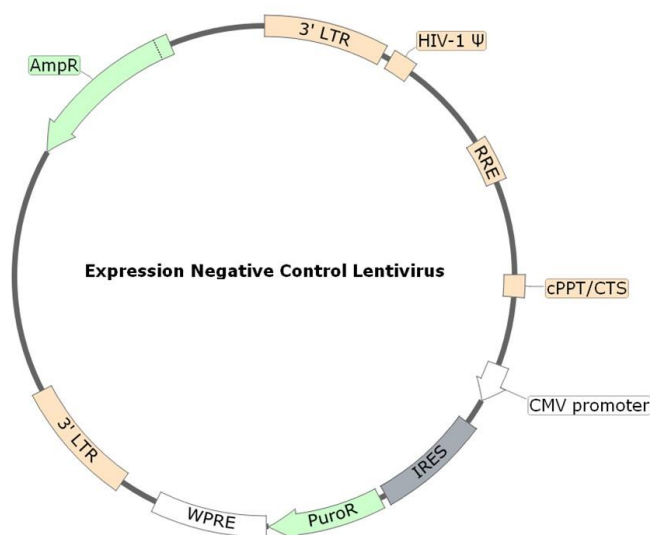
1. Useful as controls for the target expression lentivirus having puromycin selection marker.
2. Generation of stable cell line with puromycin selection without expressing a specific protein under the CMV promoter, to validate the specificity of any target expression effects.

#### **Formulation**

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

#### **Titer**

Two vials (500  $\mu$ l x 2) of Expression Negative Control Lentivirus at a titer  $1 \times 10^7$  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](mailto:info@bpsbioscience.com)

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

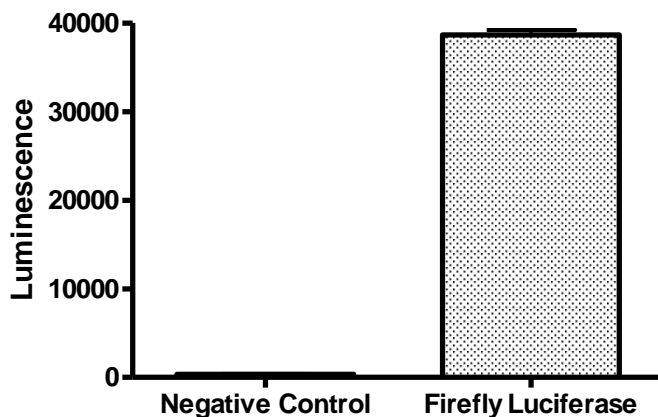
### Figure 1. Schematic of the lenti-vector used to generate the Expression Negative Control Lentivirus

#### 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-P) or expression negative control lentivirus (BPS Bioscience #79902-P) 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](mailto:info@bpsbioscience.com)

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

## Related Products

| <b><u>Product</u></b>                             | <b><u>Cat. #</u></b> | <b><u>Size</u></b> |
|---|----------------------|--------------------|
| 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          |
| FcGR11A 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](mailto:info@bpsbioscience.com)

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