

## Lenti-Luciferase-P2A-Neo Plasmid

Catalog number: 12553  
Unit size: 1 ug

| Component                        | Storage           | Amount (Cat No. 12553) |
|----------------------------------|-------------------|------------------------|
| Lenti-Luciferase-P2A-Neo Plasmid | Freeze (< -15 °C) | 1 ug                   |

### OVERVIEW

Lenti-Luciferase-P2A-Neo Plasmid is a lentiviral expression vector designed for reporter-based studies in mammalian cells. The plasmid encodes firefly luciferase followed by a neomycin resistance gene separated by a P2A self-cleaving peptide. During translation, the P2A sequence mediates ribosomal skipping, allowing production of two separate proteins from a single mRNA transcript. Luciferase expression provides a luminescent reporter signal that can be measured using standard luciferase assay reagents. Following lentiviral packaging and transduction, cells expressing the construct can be selected using neomycin (G418). This system allows establishment of stable cell populations while enabling quantitative monitoring of luciferase activity. Lentiviral vectors are commonly used for efficient gene delivery and stable integration in dividing and non-dividing mammalian cells.

### AT A GLANCE

#### Specifications:

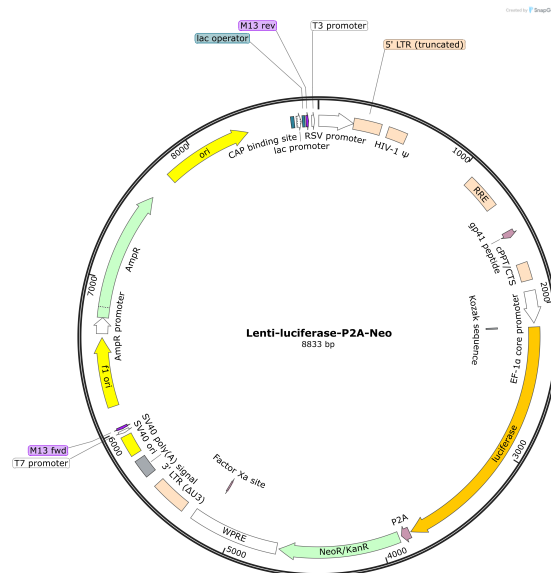
Concentration: 0.02 µg/µL in TE  
Volume: 50 µL  
Backbone size: 8833 bp  
Vector type: Cloning vector  
Promoter: EF-1α  
Bacterial Resistance(s): Ampicillin  
Copy number: High Copy

### SAMPLE EXPERIMENTAL PROTOCOL

1. Plate mammalian cells in appropriate culture medium and allow them to reach the recommended confluency (typically 60–80%) prior to transfection or transduction.
2. Introduce the plasmid into cells using a suitable transfection reagent according to the manufacturer's instructions for the chosen method.  
AAT Bioquest offers a range of Transfectamine reagents that can be used for plasmid delivery into mammalian cells.
3. Incubate cells under standard culture conditions to allow reporter expression. For constructs containing a selectable marker (e.g., neomycin resistance), apply the appropriate antibiotic to enrich for successfully transfected or transduced cells.
4. Measure luciferase activity using a compatible luciferase assay reagent and a luminometer or plate reader capable of detecting bioluminescence.

**Note:** Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.

### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Schematic map of the Lenti-Luciferase-P2A-Neo plasmid (8833 bp).

### DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.