

## Certificate of Analysis

### pGL4.51[*luc2*/CMV/Neo] Vector:

**Part No.**  
E132A

**Size**  
20µg

Part# 9PIE132  
Revised 10/16



Instructions for use of this product can be found in the pGL4 Luciferase Reporter Vectors Technical Manual #TM259, available online at:  
[www.promega.com/protocols](http://www.promega.com/protocols)

**Description:** The pGL4.51[*luc2*/CMV/Neo] Vector<sup>(a,b,c)</sup> (Cat.# E1320) encodes the luciferase reporter gene *luc2* (*Photinus pyralis*), which has been codon optimized for mammalian expression. This vector is also engineered with fewer consensus regulatory sequences for reduced backgrounds and a decreased risk of anomalous transcription.

This vector contains the following features:

- *luc2* reporter gene for expression in mammalian cells
- CMV promoter for high translational expression
- SV40 late poly(A) signal sequence is positioned downstream of *luc2* to provide efficient transcription termination and mRNA polyadenylation
- Binding region for RVprimer 3 and RVprimer 4
- Synthetic poly(A) signal/transcription start site
- Synthetic Neomycin-resistance gene for mammalian cell selection of the plasmid
- Plasmid replication origin
- *Amp<sup>r</sup>* gene for bacterial selection for vector amplification

For more information, see the *pGL4 Luciferase Reporter Vectors Technical Manual* #TM259, available online at:  
[www.promega.com/protocols](http://www.promega.com/protocols)

**Concentration:** 1µg/µl.

**GenBank® Accession Number:** EU921841.

**Storage Buffer:** The pGL4.51[*luc2*/CMV/Neo] Vector is supplied in 10mM Tris-HCl (pH 7.4), 1mM EDTA.

**Storage Conditions:** See the Product Information Label for storage temperature recommendations. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. These fluctuations can greatly alter product stability. See the label for expiration date.

#### Usage Note:

Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.

## Quality Control Assays

### Contaminant Assays

**Contaminating Nucleic Acids:** RNA, single-stranded DNA and chromosomal DNA are not evident in a specified sample of this vector as determined by agarose gel electrophoresis.

**Nuclease Assay:** Following incubation of 1µg of this vector in Restriction Enzyme Buffer at 37°C for 16–24 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

**Physical Purity:**  $A_{260}/A_{280} \geq 1.80$ ;  $A_{260}/A_{250} \geq 1.05$ .

### Functional Assays

**Identity Assay:** The vector has been sequenced completely and has 100% identity with the published sequence available at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Restriction Digestion:** The functional purity of this vector DNA is verified by successful incubation with a variety of restriction enzymes at 37°C for one hour. Samples are examined by agarose gel electrophoresis, comparing cut and uncut vector DNA with marker DNA.



AF9PIE132 1016E132



**Promega**

#### Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	<a href="http://www.promega.com">www.promega.com</a>

#### PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 2008–2016 Promega Corporation. All Rights Reserved.

GeneBank is a registered trademark of US Department of Health and Human Services.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Signed by:

R. Wheeler, Quality Assurance

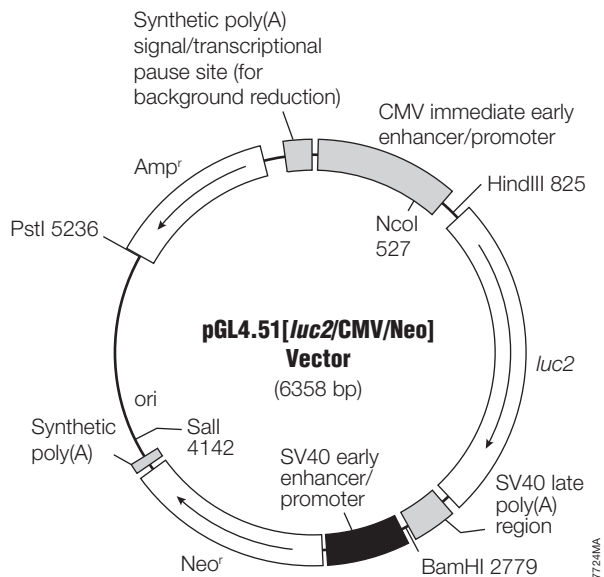
Part# 9PIE132  
Printed in USA Revised 10/16.



# Promega

## Features list and map for the pGL4.51[*luc2*/CMV/Neo] Vector

CMV immediate early enhancer/promoter	14–755
<i>luc2</i>	859–2511
SV40 late poly(A) region	2546–2767
SV40 early enhancer/promoter	2815–3233
Synthetic neomycin phosphotransferase coding region (Neo <sup>r</sup> )	3258–4055
Synthetic poly(A)	4077–4125
Reporter vector primer 4 binding region	4357–4365
Replication origin	4449
Synthetic beta-lactamase (Amp <sup>r</sup> ) coding region	5240–6100
Synthetic poly(A) signal/transcriptional pause region	6205–6358
Reporter vector primer 3 binding region	6307–6326



©BY USE OF THIS PRODUCT, RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE LABEL LICENSE. If the researcher is not willing to accept the terms of this label license, and the product is unused, Promega will accept return of the unused product and provide the researcher with a full refund.

Researchers may use this product for research use only, no commercial use is allowed. "Commercial use" means any and all uses of this product and derivatives by a party for money or other consideration and may include but is not limited to use in: (1) product manufacture; and (2) to provide a service, information or data; and/or resale of the product or its derivatives, whether or not such product or derivatives are resold for use in research. Researchers shall have no right to modify or otherwise create variations of the nucleotide sequence of the luciferase gene except that researchers may: (1) create fused gene sequences provided that the coding sequence of the resulting luciferase gene has no more than four deoxynucleotides missing at the affected terminus compared to the intact luciferase gene sequence, and (2) insert and remove nucleic acid sequences in splicing research predicated on the inactivation or reconstitution of the luminescence of the encoded luciferase. No other use or transfer of this product or derivatives is authorized without the prior express written consent of Promega. In addition, researchers must either: (1) use luminescent assay reagents purchased from Promega for all determinations of luminescence activity of this product and its derivatives; or (2) contact Promega to obtain a license for use of the product and its derivatives. Researchers may transfer derivatives to others for research use provided that at the time of transfer a copy of this label license is given to the recipients and recipients agree to be bound by the terms of this label license. With respect to any uses outside this label license, including any diagnostic, therapeutic or prophylactic uses, please contact Promega for supply and licensing information. PROMEGA MAKES NO REPRESENTATIONS OR WARRANTIES OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING FOR MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE WITH REGARDS TO THE PRODUCT. The terms of this label license shall be governed under the laws of the State of Wisconsin, USA. This label license relates to Promega patents and/or patent applications on improvements to the luciferase gene.

©U.S. Pat. No. 8,008,006 and European Pat. No. 1341808.

©U.S. Pat. No. 7,728,118.

Part# 9PIE132  
Printed in USA Revised 10/16.