# **Certificate of Analysis**

## pGL4.13[*luc2*/SV40] Vector:

**Part No. Size** E668A 20µg



**Instructions for use** of this product can be found in the pGL4 Vectors Technical Manual #TM259, available online at: **www.promega.com/protocols** 

**Description:** The pGL4.13[*luc2*/SV40] Vector<sup>(a,b,c)</sup> encodes the luciferase reporter gene *luc2* (*Photinus pyralis*) and is designed for high expression and reduced anomalous transcription. The pGL4 Vectors are engineered with fewer consensus regulatory sequences and a synthetic gene, which has been codon optimized for mammalian expression.

The pGL4.13[/uc2/ SV40] Vector contains the /uc2 reporter gene and the SV40 early enhancer/promoter for use as an expression control or a co-reporter vector.

Concentration: 1µg/µl.

GenBank® Accession Number: AY738225.

Storage Buffer: The pGL4.13[/uc2/SV40] Vector is supplied in 10mM Tris-HCI (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 expiration date on the product information label.

#### **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 specified quantities of the vector as determined by agarose gel electrophoresis.

**Nuclease Assay:** Following incubation of 1µg of the 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:**  $A_{260}/A_{280} \ge 1.80$ ,  $A_{260}/A_{250} \ge 1.05$ .

### **Functional Assays**

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

**Restriction Digestion:** The functional purity of the vector DNA is verified by successful digestion with restriction enzymes at the optimal temperature for one hour. Samples are examined by agarose gel electrophoresis, comparing cut and uncut vector DNA with marker DNA.

# Part# 9PIE668 Revised 10/16



AE9PIE668 1016E668



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	www.promega.com

#### 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 armful 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 WARANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT IMITATION, 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 but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tort (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tort (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tort (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tort (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.

GenBank is a registered trademark of the U.S. 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.

Part# 9PIE668 Printed in USA. Revised 10/16.



R. Wheeler, Quality Assurance



## pGL4.13[luc2/SV40] Vector Features List and Map

SV40 early enhancer/promoter	51-469
luc2 reporter gene	499-2151
SV40 late poly(A) region	2186-2407
Reporter Vector primer 4 (RVprimer4) binding region	2475-2494
ColE1-derived plasmid replication origin	2732
Synthetic β-lactamase (Amp <sup>r</sup> ) coding region	3523-4383
Synthetic poly(A) signal/transcriptional pause region	4488-4641
Reporter Vector primer 3 (RVprimer3) binding region	4590-4609

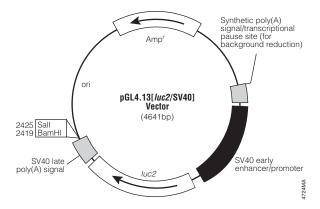


Figure 1. pGL4.13[luc2/SV40] Vector circle map.

Sequence information and restriction enzyme tables for the pGL4 Vectors are available

online at: www.promega.com/vectors

Further information on the use of pGL4 Vectors is available in Technical Manual #TM259,

which is available online at:  $\begin{tabular}{ll} www.promega.com/protocols \\ \end{tabular}$ 

(a)BY USE OF THIS PRODUCT, RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE LABEL LICENSE.

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, and (2) insert and remove nucleic acid sequences in splicing research. No other use or transfer of this product or derivatives is authorized. Researchers must either (1) use luminescence activity of this product and its derivatives; or (2) contact Promega to obtain a license for use of the product. For any uses outside this label license, contact Promega for supply and licensing information. This product is for research use only; no commercial use is allowed. For a full copy of this label license, including the definition of "commercial use," go to: www.promega.com/LULL

(b)U.S. Pat. No. 8,008,006 and European Pat. No. 1341808.

(c)Patents Pending.

Part# 9PIE668 Printed in USA. Revised 10/16.