Certificate of Analysis

pGL4.17[/uc2/Neo] Vector:

Part No. E672A



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/

Description: The pGL4.17[luc2/Neo] Vector(a-d) encodes the luciferase reporter gene luc2 (Photinus pyralis) and is designed for high expression and reduced anomalous transcription. This vector also contains a mammalian selectable marker for neomycin resistance in which the number of transcription factor-binding sites has been reduced and mammalian codon usage optimized. This vector is also engineered with fewer consensus regulatory sequences for reduced background and a decreased risk of anomolous transcription and has a synthetic reporter gene, which is codon-optimized for mammalian

The pGL4.17[/uc2/Neo] Vector is a basic vector with no promoter. However, the vector contains a multiple cloning region to allow cloning of a promoter of choice.

Concentration: 1µg/µl.

GenBank® Accession Number: DQ188837.

Storage Buffer: The pGL4.17[/uc2/Neo] Vector is supplied in 10mM Tris-HCI (pH 7.4), 1mM EDTA.

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

Usage Notes:

- 1. For easy transfer from one pGL4 Vector to another, the multiple cloning region is consistent throughout the pGL4 Vector series. For easy transfer between pGL3 Vectors and pGL4 Vectors, many of the restriction enzyme sites present in the pGL3 Vectors are also present in the pGL4 Vectors.
- 2. Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.

Quality Control Assays

Nuclease Assay: Following incubation of 1µg of pGL4.17[/uc2/Neo] Vector in standard restriction digest buffers at 37°C for 16-24 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

Physical Purity: $A_{260}/A_{280} \ge 1.80$, $A_{260}/A_{250} \ge 1.05$ at pH 7.4.

Sequence: The pGL4.17[/uc2/Neo] Vector has been completely sequenced and has 100% identity with the published sequence, available at: www.promega.com/vectors/

(a) BY LISE OF THIS PRODUCT. RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED LISE 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 seguences, and (2) insert and remove nucleic acid seguences in splicing research. No other use or transfer of this product or derivatives is authorized. 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. 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)Patent Pending.

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

(d)U.S. Pat. No. 7,728,118.

Ren Whele Signed by:

Part# 9PIE672 Revised 10/16



AF9PIF672 1016F672



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 harmful if misused. Due care should be exercised with all Promega products to prevent direct human

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed 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 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 one 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, tord (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tord (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tord (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tord (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tord (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tord (including but not limited to loss of use, revenue or profit) whether based upon warranty, contract, tord (including the result of the failure of Promega products to perform in accordance with the stated specifications.

© 2005–2009, 2013, 2015, 2016 Promega Corporation. All Rights Reserved.

GenBank 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.

Part# 9PIE672 Printed in USA. Revised 10/16.



pGL4.17[/uc2/Neo] Vector Features List and Maps Multiple cloning region 1 - 70luc2 reporter gene 100-1752 SV40 late poly(A) signal 1787-2008 SV40 early enhancer/promoter 2056-2474 Synthetic neomycin phosphotransferase (Neor) coding region 2499-3293 3318-3366 Synthetic poly(A) signal Reporter Vector primer 4 (RVprimer4) binding region 3433-3452 ColEI-derived plasmid replication origin 3690 Synthetic B-lactamase (Ampr) coding region 4481-5341 Synthetic poly(A) signal/transcriptional pause site 5446-5599 Reporter Vector primer 3 binding region 5548-5567

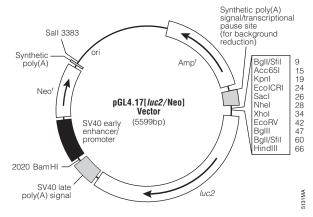


Figure 1. pGL4.17 Vector circle map.

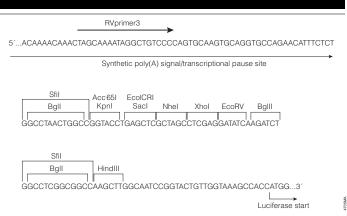


Figure 2. The multiple cloning region of the pGL4 Vectors.

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: www.promega.com/protocols/