

Certificate of Analysis

pGL4.11[*luc2P*] Vector:

Part No. E666A
Size 20µg

Part# 9PIE666
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

Description: The pGL4.11[*luc2P*] Vector^(a,b,c) encodes the luciferase reporter gene *luc2P* (*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 pGL4.11[*luc2P*] Vector is a basic vector with no promoter. However, it contains a multiple cloning region to allow cloning of a promoter of choice. The *luc2P* reporter gene contains hPEST, a protein destabilization sequence. The protein encoded by *luc2P* responds more quickly and with greater magnitude to changes in transcriptional activity than the *luc2* gene, its more stable counterpart.

Concentration: 1µg/µl.

GenBank® Accession Number: AY738223.

Storage Buffer: The pGL4.11[*luc2P*] 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 expiration date on the product information label.

Usage Notes:

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



AF9PIE666 1016E666



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	www.promega.com

Quality Control Assays

Nuclease Assay: Following incubation of 1µg of pGL4.11[*luc2P*] Vector in standard restriction digest buffers at 37°C for 16–24 hours, no evidence of nuclease activity was detected by agarose gel electrophoresis.

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

Sequence: The pGL4.11[*luc2P*] Vector has been completely sequenced and has 100% identity with the published sequence, available at: www.promega.com/vectors

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.

© 2004–2016 Promega Corporation. All Rights Reserved.

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.

Signed by:

R. Wheeler, Quality Assurance

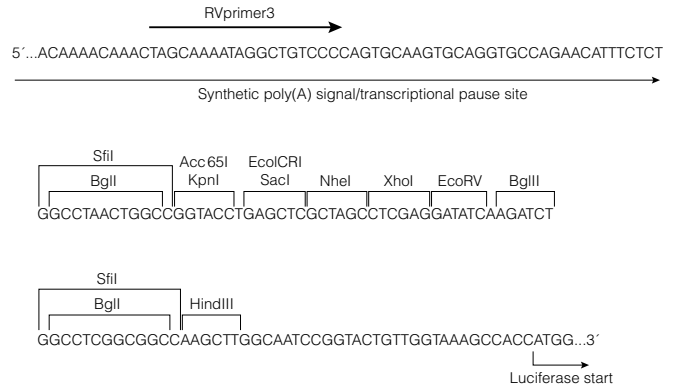
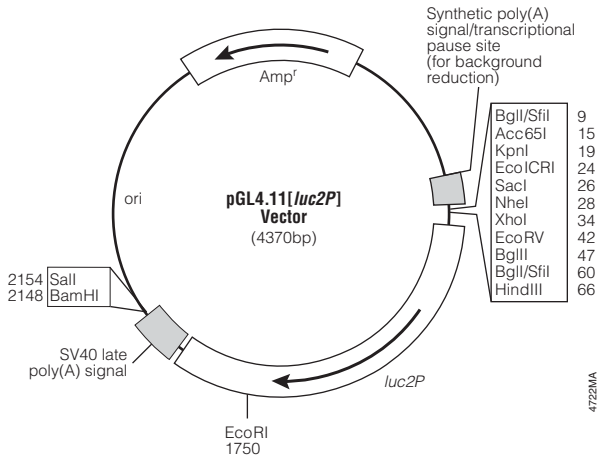
Part# 9PIE666
Printed in USA. Revised 10/16.



Promega

pGL4.11[*luc2P*] Vector Features List and Map

Multiple cloning region	1-70
<i>luc2</i> reporter gene (synthetic firefly luciferase; includes hPEST)	100-1875
SV40 late poly(A) region	1915-2136
Reporter Vector primer 4 (RVprimer4) binding region	2204-2223
<i>ColE1</i> -derived plasmid replication origin	2461
Synthetic β -lactamase (<i>Amp^r</i>) coding region	3252-4112
Synthetic poly(A) signal/transcriptional pause site	4217-4370
Reporter Vector primer 3 (RVprimer3) binding region	4319-4338



Multiple cloning region of the pGL4.11[*luc2P*] Vector.

©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 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.

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

©Patents Pending.

Part# 9PIE666

Printed in USA. Revised 10/16.