Product Contents

pGL4.20[/uc2/Puro] Vector:

Part No. Size E675A 20μg



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 pGL4.20[/uc2/Puro] Vector(a-d) encodes the luciferase reporter gene /uc2 (Photinus pyralis) and is designed for high expression and reduced anomalous transcription. This vector also contains a mammalian selectable marker for puromycin resistance in which the number of transcription factor binding sites has been reduced and mammalian codon usage optimized. This vector is engineered with fewer consensus regulatory sequences than the pGL3 Vectors and a synthetic reporter gene, which has been codon optimized for mammalian expression.

The pGL4.20[luc2/Puro] Vector is a basic vector with no promoter. However, it contains a multiple cloning region to allow for the cloning of a promoter of choice.

Concentration: 1µg/µl.

GenBank® Accession Number: DQ188840

Storage Buffer: The pGL4.20[luc2/Puro] Vector is supplied in 10mM Tris-HCI (pH 7.4), 1mM EDTA.

Storage Conditions: Store the pGL4.20[/uc2/Puro] Vector at -20°C. 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:

- 1. The method used to purify this vector yields DNA that may be suitable for transfection of mammalian cells without further manipulation
- 2. 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.
- 3. 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.

Summary of Changes:

The following change was made to the 6/15 revision of this document: Expired product disclaimers were removed.

Part# 9PIE675 Revised 10/16





Promega Corporation	n
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 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 WAR-RANTIES 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 lim-Each Promega product is shipped with documentation able, 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

© 2005–2009, 2013, 2015, 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

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:

fler Wheeler

Part# 9PIF675

Printed in USA Revised 10/16.



pGL4.20[luc2/Puro] Vector Features List and Maps

Multiple cloning region	1-70
luc2 reporter gene	100-1752
SV40 late poly(A) signal	1787-2008
SV40 early enhancer/promoter	2056-2474
Synthetic puromycin-N-acetyltransferase (Puror) coding region	2499-3098
Synthetic poly(A) signal	3123-3171
Reporter Vector primer 4 binding region	3238-3257
Col El-derived plasmid replication origin	3495
Synthetic β-lactamase (Amp ^r) coding region	4286-5146
Synthetic poly(A) signal/transcriptional pause site	5251-5404
Reporter Vector primer 3 binding region	5353-5372

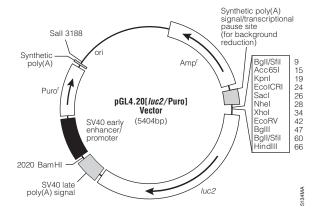


Figure 1. pGL4.20[/uc2/Puro] Vector circle map.

Sfil Acc65I EcolCRI BgIl Kpnl Sacl Nhel Xhol EcoRV BgIll GGCCTAACTGGCCGGTACCTGAGCTCGCAGGTACAGATCT Sfil BgIl HindIII GGCCTCGGCGGCCAAGCTTGGCAATCCGGTACTGTTGGTAAAGCCACCATGG...3' Luciferase start

Figure 2. Multiple cloning region of the pGL4.20[luc2/Puro] Vector.

Sequence information and restriction enzyme tables for the pGL4 Vectors are available online at: www.promega.com/vectors

Further information on use of the pGL4 Vectors is available in Technical Manual #TM259 available online at: www.promega.com/protocols

Summary of Changes:

The following change was made to the 6/15 revision of this document: Expired product disclaimers were removed.

(a)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 luminescence activity of this product and its derivatives; or (2) contact 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

(b)U.S. Pat. No. 5,670,356

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

(d)Patent Pending

Part# 9PIE675
Printed in USA Revised 10/16