Certificate of Analysis

pGL4.28[/uc2CP/minP/Hygro] Vector:

Part No. Size E846A 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 pGL4.28[/uc2CP/minP/Hygro] Vector(a-e) encodes the luciferase reporter gene /uc2CP and is designed for high expression and reduced anomalous transcription. The vector contains a multiple cloning region for insertion of a response element of interest upstream of a minimal promoter and the *luc2CP* gene. *luc2CP* is a synthetically derived luciferase sequence with humanized codon optimization. The *luc2CP* gene also contains hCL1 and hPEST, both of which are protein destabilization sequences. The protein encoded by Juc2CP responds more quickly than the protein encoded by the luc2 gene upon induction. The vector backbone contains an ampicillin resistance gene to allow for selection in E. coli and a mammalian selectable marker for hygromycin resistance.

Concentration: 1µg/µl.

GenBank® Accession Number: DQ904460.

Storage Buffer: The pGL4.28[/uc2CP/minP/Hygro] Vector is supplied in 10mM Tris-HCI (pH 7.4), 1mM EDTA.

Storage Conditions: See the Product Information Label for storage 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 Label

Usage Note: 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 the vector in restriction digest buffer B at 37°C for 16 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 pGL4.28[/uc2CP/minP/Hygro] Vector has been completely sequenced and has 100% identity with the published sequence, available at: www.promega.com/vectors

(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 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 exequences, 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 for usually 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 RE Researchers may use this product for research use only, no commercial use is allowed. "Commercial use" means any and all

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

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

(d)Patent Pending.

Signed by:

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

Ren Wheeler

R. Wheeler. Quality Assurance

Part# 9PIE846 Revised 10/16



AF9PIF846 1016F846



Promega Corporation	on
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, 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 the result of the failure of Promega products to perform in accordance with the stated specifications.

© 2006-2009, 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.

Part# 9PIE846 Printed in USA. Revised 10/16.



pGL4.28[luc2CP/minP/Hygro] Vector Features List and Maps

Minimal promoter	78–108
<i>luc2P</i> reporter gene	141-1970
SV40 late poly(A) region	2007–2228
SV40 early enhancer/promoter	2276-2694
Synthetic hygromycin (Hygr) coding region	2719-3756
Synthetic poly(A) signal	3780-3828
Reporter vector primer 4 (RVprimer4) binding region	3895-3914
Co/E1-derived plasmid replication origin	4152
Synthetic β -lactamase (Ampr) coding region	4943-5803
Synthetic poly(A) signal/transcriptional pause site	5908-6061
Reporter vector primer 3 (RVprimer3) binding region	6010-6029

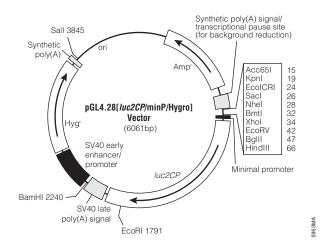


Figure 1. The pGL4.28[*luc2CP*/minP/Hygro] Vector map.

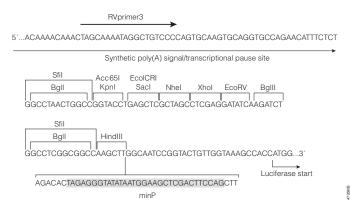


Figure 2. Multiple cloning region of the pGL4.28[/uc2CP/minP/Hygro] Vector.

Sequence information, vector maps 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, available online at: www.promega.com/protocols

Summary of Changes, 8/15 Revision

The following changes were made to the 8/15 version of this document: Legal disclaimers were updated to remove expired information.