

Certificate of Analysis

pGL4.32[*luc2P/NF-κB-RE/Hygro*] Vector:

Part No. Size
E849A 20µg

Part# 9PIE849
Revised 7/16



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 pGL4.32[*luc2P/NF-κB-RE/Hygro*] Vector^(a-c) contains five copies of an NF-κB response element (NF-κB-RE) that drives transcription of the luciferase reporter gene *luc2P* (*Photinus pyralis*). *luc2P* is a synthetically derived luciferase sequence with humanized codon optimization that is designed for high expression and reduced anomalous transcription. The *luc2P* gene contains hPEST, a protein destabilization sequence. The protein encoded by *luc2P* responds more quickly than the protein encoded by the *luc2* gene upon induction. The vector backbone contains an ampicillin resistance gene to allow selection in *E. coli* and a mammalian selectable marker for hygromycin resistance.

Concentration: 1µg/µl.

GenBank® Accession Number: EU581860.

Storage Buffer: The pGL4.32[*luc2P/NF-κB-RE/Hygro*] 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 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 Enzyme Buffer at 37°C for 16–24 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

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

Sequence: The pGL4.32[*luc2P/NF-κB-RE/Hygro*] Vector has been completely sequenced and has 100% identity with the published sequence, available at: www.promega.com/vectors

Signed by:

R. Wheeler, Quality Assurance

^(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 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. 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 WARRANTIES OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING FOR MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE WITH REGARDS TO THE PRODUCT. The terms of this label license shall be governed under the laws of the State of Wisconsin, USA. This label license relates to Promega patents and/or patent applications on improvements to the luciferase gene.

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

^(c)U.S. Pat. No. 7,728,118.



AF9PIE849 0716E849



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

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.

© 2008–2016 Promega Corporation. All Rights Reserved.

Dual-Glo and Dual-Luciferase are registered trademarks of and Bright-Glo and ONE-Glo are trademarks of Promega Corporation.

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.

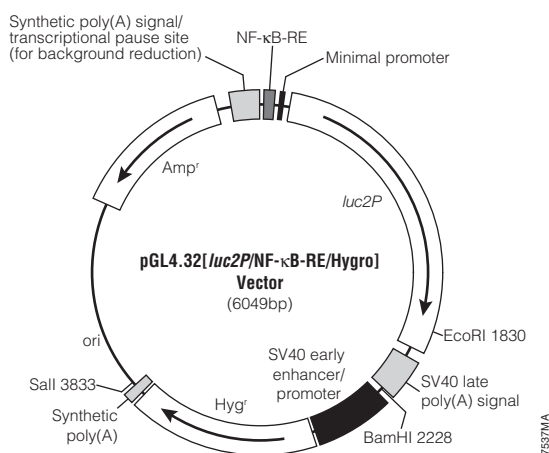
Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIE849
Printed in USA, Revised 7/16.



pGL4.32[*luc2P*/NF-κB-RE/Hygro] Vector Features List and Map:

NF-κB response element	33–84
Minimal promoter	117–147
<i>luc2P</i> reporter gene	180–1955
SV40 late poly(A) signal	1995–2216
SV40 early enhancer/promoter	2264–2682
Synthetic hygromycin (Hyg ^r) coding region	2707–3744
Synthetic poly(A) signal	3768–3816
Reporter Vector primer 4 (RVprimer4) binding region	3883–3902
<i>ColE1</i> -derived plasmid replication origin	4140
Synthetic β-lactamase (Amp ^r) coding region	4931–5791
Synthetic poly(A) signal/transcriptional pause site	5896–6049
Reporter Vector primer 3 (RVprimer3) binding region	5998–6017



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

Sample Protocol to Determine Induction of Luciferase by TNFα in HEK 293 Cells Transfected with the pGL4.32[*luc2P*/NF-κB-RE/Hygro] Vector

Materials to be Supplied by User

- Dulbecco's PBS (DPBS)
- 0.05% (w/v) trypsin in DPBS
- DMEM
- DMEM supplemented with 10% fetal bovine serum (DMEM/FBS)
- Tumor necrosis factor-α (Sigma T0157), 10μg/ml solution in PBS containing 1mg/ml BSA
- Bright-Glo™ Luciferase Assay System (Cat.# E2610)
- HEK 293 cells

Day 1: Plate Cells

1. Grow HEK 293 cells in DMEM/FBS to approximately 75% confluency.
2. Harvest cells via trypsinization: Remove the DMEM/FBS, wash the cells with DPBS and add the trypsin/DPBS (1X volume). After 2 minutes, add a 4X volume of DMEM/FBS, collect the cell suspension and pellet the cells by centrifugation. Aspirate the supernatant and resuspend in DMEM/FBS. We have routinely used a concentration of 15,000 viable cells/100μl DMEM/FBS.
3. Dispense 100μl of the cell suspension into the wells of a 96-well plate. Plate enough wells to perform each test condition in triplicate.
4. Cover the plate and place it in a tissue culture incubator at 37°C overnight (or for 24 hours).

Day 2: Transfect Cells

1. Transfect the cells using a high-efficiency transfection reagent. Each well of 96-well plate to be transfected requires 0.1μg pGL4.32[*luc2P*/NF-κB-RE/Hygro] plasmid DNA. Transfection conditions may require optimization.
2. Cover the plate and place it in a tissue culture incubator at 37°C overnight or as needed for cell recovery depending on the transfection method used. We have used 24 hours recovery time for lipid-mediated transfections.

Day 3: Induce Transfected Cells

1. Prepare 1X induction and 1X control solutions. Calculate the volume of 1X induction and 1X control solution by multiplying the number of wells needed for each solution by 100μl and prepare 110% of this amount.
 - 1X induction solution: Dilute 10μg/ml TNFα solution to 20ng/ml in DMEM/FBS. Final TNFα concentration will be 20ng/ml.
 - 1X control solution: DMEM/FBS.
2. Remove media from wells that will be treated with either 1X induction solution or 1X control solution.
3. Add 100μl of 1X induction solution to the cells to be induced and 100μl of 1X control solution to the control noninduced cells.
4. Return the plate to the tissue culture incubator and induce for 5 hours.
5. Analyze luciferase activity using an appropriate luciferase detection assay. We have observed comparable results for fold induction of the vector using a variety of luciferase reagents, including: Bright-Glo™ Luciferase Assay System (Cat.# E2610, Technical Manual #TM052); ONE-Glo™ Luciferase Assay System (Cat.# E6110, Technical Manual #TM292); Dual-Luciferase® Reporter Assay System (Cat.# E1910, Technical Manual #TM040); and Dual-Glo® Luciferase Assay System (Cat.# E2920, Technical Manual #TM058).
6. Calculate the fold induction as follows:

$$\text{Fold Induction} = \frac{\text{Average relative light units of induced cells}}{\text{Average relative light units of control cells}}$$