

## Certificate of Analysis

### pGL4.36[*Luc2P*/MMTV/Hygro] Vector

Part No.  
E136A

Size  
20µg

Part# 9PIE136

Revised 1/18



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/](http://www.promega.com/protocols/)

**Description:** The pGL4.36[*Luc2P*/MMTV/Hygro] Vector<sup>(a-c)</sup> contains MMTV LTR (Murine Mammary Tumor Virus Long Terminal Repeat) that drives transcription of the luciferase reporter gene *Luc2P* in response to activation of several nuclear receptors such as glucocorticoid receptor and androgen receptor. *Luc2P* is a synthetically derived luciferase sequence with humanized codon optimization. The *Luc2P* gene also contains hPEST, a protein destabilization sequence. The protein encoded by *Luc2P* responds more quickly to induction than the protein encoded by the *Luc2* gene. The vector backbone contains an ampicillin resistance gene to allow selection in *E. coli* and the mammalian selectable marker for hygromycin resistance.

**Concentration:** 1µg/µl.

**GenBank® Accession Number:** FJ773214.

**Storage Buffer:** The pGL4.36[*Luc2P*/MMTV/Hygro] Vector is supplied in 10mM Tris-HCl, 1mM EDTA (pH 8.0).

**Storage Conditions:** See the Product Information Label for storage temperature recommendations and expiration date.

**Usage Note:** Mix well prior to use.



AF9PIE136 0118E136



## 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	<a href="http://www.promega.com">www.promega.com</a>

## Quality Control Assays

### Contaminant Assays

**Contaminating Nucleic Acid Assay:** RNA, single-stranded DNA and chromosomal DNA are not evident in a specified sample of this vector as determined by agarose gel electrophoresis.

**Nuclease Assay:** Following incubation of 1µg of this 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:**  $A_{260}/A_{280} \geq 1.80$ ,  $A_{260}/A_{250} \geq 1.05$ .

### Functional Assays

**Identity Assay:** The vector has been sequenced completely and has 100% identity with the published sequence available at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Restriction Digestion:** The functional purity of this vector DNA is verified by complete digestion with several restriction enzymes at 37°C for 1 hour. Samples are examined by agarose gel electrophoresis, and cut and uncut vector DNA are compared with marker DNA.

Signed by:

R. Wheeler, Quality Assurance

### 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.

©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 REGARD 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.

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

©U.S. Pat. No. 7,728,118.

© 2009, 2015, 2016, 2018 Promega Corporation. All Rights Reserved.

GenBank is a registered trademark of U.S. Department of Health and Human Services.

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

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.

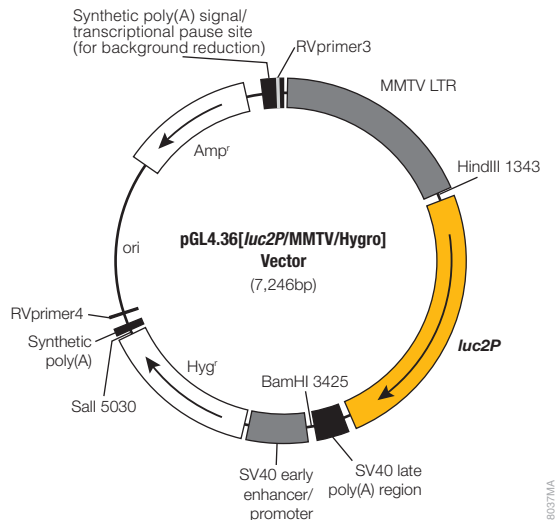
Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIE136  
Printed in USA Revised 1/18



## Features List and Map for the pGL4.36[*luc2P*/MMTV/Hygro] Vector

MMTV LTR	14–1328
<i>luc2P</i> reporter gene	1377–3152
SV40 late poly(A) signal	3192–3413
SV40 early poly(A) enhancer/promoter	3461–3879
synthetic hygromycin (Hyg <sup>r</sup> ) coding region	3904–4941
synthetic poly(A) region	4965–5013
reporter vector primer 4 (RVprimer4) binding region	5080–5099
ColE1-derived plasmid replication origin	5337
synthetic β-lactamase (Amp <sup>r</sup> ) coding region	6128–6988
synthetic poly(A) signal/transcription pause site	7093–7246
reporter vector primer 3 (RVprimer3) binding region	7195–7214



**Figure 1. pGL4.36[*luc2P*/MMTV/Hygro] Vector map.**

Sequence information, vector maps and restriction enzyme tables for the pGL4 Vectors are available online at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

Additional information is available in Technical Manual #TM259, available online at: [www.promega.com/protocols/](http://www.promega.com/protocols/)

## Sample Protocol to Determine Induction of Luciferase by Dexamethasone in HeLa Cells Transfected with the pGL4.36[*luc2P*/MMTV/Hygro] Vector

### Materials to Be Supplied by the User

- 1X PBS
- 0.05% (w/v) trypsin
- DMEM with 10% fetal bovine serum (growth medium)
- DMEM without phenol red
- DMEM without phenol red supplemented with 5% charcoal/dextran-treated fetal bovine serum (assay medium)
- dexamethasone (Sigma Cat.# D4902), 10mM solution in ethanol
- Bright-Glo™ Luciferase Assay System (Cat.# E2610)
- transfection reagent
- HeLa cells

### Day 1: Plate Cells

Seed HeLa cells at 10,000 cells/ well in a solid white 96-well tissue culture-treated plate using phenol red-free DMEM containing 5% charcoal/dextran-treated FBS (80µl/well).

### Day 2: Transfect Cells

1. Transfect the cells using a high-efficiency transfection reagent. Each well of cells in a 96-well plate requires 0.1µg of pGL4.36[*luc2P*/MMTV/Hygro] Vector DNA. Transfection conditions may require optimization. We have routinely added approximately 10µl/well of a transfection master mix.
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 use 24 hours of recovery time for lipid-mediated transfections.

### Day 3: Induce Transfected Cells

1. Prepare 10X induction and 10X control solution. Calculate the volume of 10X induction and 10X control solution by multiplying the number of wells needed for each solution by 10µl, and prepare 110% of this amount.
  - 10X induction solution: Dilute 10mM dexamethasone solution in DMEM without phenol red to 100µM (1:100 dilution). The final dexamethasone concentration will be 10µM.
  - Control solution: Add the same volume of ethanol only used in 10X induction solution to DMEM without phenol red.
2. Add 10µl of 10X induction solution to wells to be induced or control solution to control noninduced wells.
3. Return the plate to the tissue culture incubator, and induce for overnight to 24 hours.

### Day 4: Harvest and Analyze Luciferase Activity

1. 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).

2. Calculate the fold induction as follows:

$$\text{fold induction} = \frac{\text{average relative light units of induced cells}}{\text{average relative light units of control cells}}$$