

## Product Contents

### pGL4.30[*luc2P*/NFAT-RE/Hygro] Vector:

<b>Part No.</b>	<b>Size</b>
E848A	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/tbs](http://www.promega.com/tbs)

**Description:** The pGL4.30[*luc2P*/NFAT-RE/Hygro] Vector<sup>(a,b,c,d,e)</sup> contains an NFAT response element (NFAT-RE) that drives the 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.

See the *pGL4 Luciferase Reporter Vectors Technical Manual #TM259* for more information.

**Concentration:** 1µg/µl.

**GenBank® Accession Number:** DQ904462.

**Storage Buffer:** The pGL4.30[*luc2P*/NFAT-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 digest buffer B at 37°C for 16 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

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

**Sequence:** The pGL4.30[*luc2P*/NFAT-RE/Hygro] Vector has been completely sequenced and has 100% identity with the published sequence, available at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

Part# 9PIE848  
Revised 10/09



**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>

### 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.

<sup>(a)</sup>READ THIS FIRST BEFORE OPENING PRODUCT.

The sale of this product and its use by the purchaser are subject to the terms of a limited use label license, the full text of which is shipped with this product and also available at: [www.promega.com/LULL](http://www.promega.com/LULL). That text must be read by the purchaser prior to opening this product to determine whether the purchaser agrees that all use of the product shall be in accordance with the license terms. If the purchaser is not willing to accept the terms of the limited use label license, Promega is willing to accept the return of the unopened product and provide the purchaser with a full refund. However, if the product is opened for any reason, then the purchaser agrees to be bound by the terms of the limited use label license.

<sup>(b)</sup>Patents Pending.

<sup>(c)</sup>Australian Pat. No. 2001-285278 and other patents pending.

<sup>(d)</sup>The method of recombinant expression of *Coleoptera* luciferase is covered by U.S. Pat. Nos. 5,583,024, 5,674,713 and 5,700,673. A license (from Promega for research reagent products and from The Regents of the University of California for all other fields) is needed for any commercial sale of nucleic acid contained within or derived from this product.

<sup>(e)</sup>The NFAT response element and its use are licensed under one or more of the following patents: U.S. Pat. Nos. 5,837,840, 6,197,925, 5,989,810, 6,096,515, 6,388,052, 6,150,099, 6,171,781, 6,352,830, 6,312,899 and 6,875,571.

© 2006, 2008, 2009 Promega Corporation. All Rights Reserved.

Bright-Glo is a trademark of Promega Corporation.

GenBank is a registered trademark of the U.S. Department of Health and Human Services. *TransIT* is a registered trademark of Mirus Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

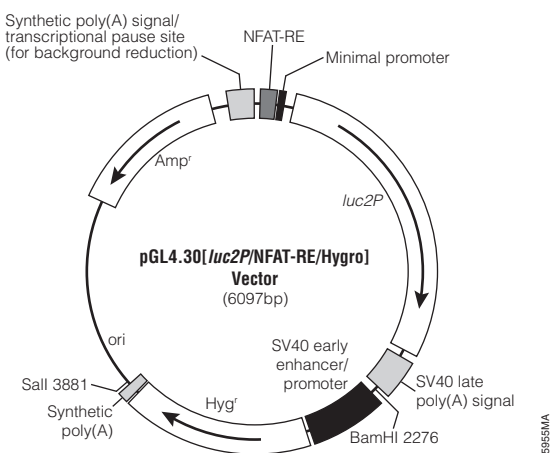
All specifications are subject to change without prior notice.

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# 9PIE848  
Printed in USA, Revised 10/09.

### pGL4.30[*luc2P*/NFAT-RE/Hygro] Vector Features List and Map:

NFAT response element	33–122
Minimal promoter	136–166
<i>luc2P</i> reporter gene	228–2003
SV40 late poly(A) signal	2043–2264
SV40 early enhancer/promoter	2312–2730
Synthetic hygromycin (Hyg <sup>r</sup> ) coding region	2755–3792
Synthetic poly(A) signal	3816–3864
Reporter Vector primer 4 (RVprimer4) binding region	3931–3950
<i>ColE1</i> -derived plasmid replication origin	4188
Synthetic β-lactamase (Amp <sup>r</sup> ) coding region	4979–5839
Synthetic poly(A) signal/transcriptional pause site	5944–6097
Reporter Vector primer 3 (RVprimer3) binding region	6046–6065



**Figure 1. Vector map.**

### Sample Protocol to Determine Luciferase Induction by Ionomycin/PMA in HEK 293 Cells Transfected with pGL4.30[*luc2P*/NFAT-RE/Hygro] Vector

#### Materials to be Supplied by User

- Dulbecco's PBS (DPBS)
- 0.05% (w/v) trypsin in DPBS
- DMEM supplemented with 10% fetal bovine serum (DMEM/FBS)
- *TransIT*<sup>®</sup>-LT1 (Mirus Part# MIR2304)
- DMSO
- 1mM ionomycin in DMSO
- 1mg/ml phorbol 12-myristate 13-acetate (PMA) in DMSO
- Bright-Glo<sup>™</sup> 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 the cells via trypsinization. Remove the DMEM/FBS, wash the cells with DPBS and add the trypsin/DPBS (1X volume). After 2 minutes, add 4X volume of DMEM/FBS, collect the cell suspension and pellet the cells by centrifugation. Aspirate supernatant and resuspend in DMEM/FBS at a concentration of 20,000 viable cells/90μl DMEM/FBS.
3. Dispense 90μl of the cell suspension to 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. Prepare the DNA transfection master mix. Each well of 96-well plate to be transfected requires 10μl DMEM, 0.3μl *TransIT*<sup>®</sup>-LT1 and 0.1μg pGL4.30[*luc2P*/NFAT-RE/Hygro] plasmid DNA. To prepare the master mix, calculate the total number of wells that will be transfected and prepare 110% of this amount. It is recommended that at least 10 wells of master mix be prepared.
  - For each well, mix 10μl DMEM and 0.3μl *TransIT*<sup>®</sup>-LT1 in a microcentrifuge tube, briefly vortex at maximum setting and incubate at room temperature for 15 minutes.
  - For each well to be transfected, add 0.1μg of pGL4.30[*luc2P*/NFAT-RE/Hygro] Vector to the DMEM/*TransIT*<sup>®</sup>-LT1, vortex briefly and incubate at room temperature for 15 minutes.
2. Add 10μl of master mix to each well that is to be transfected.
3. Cover the plate and place it in a tissue culture incubator at 37°C overnight (or for 24 hours).

#### Day 3: Induce Transfected Cells and Measure Luciferase Activity

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 1mM stock ionomycin solution to 1μM (1:1,000) and 1mg/ml stock solution of PMA to 10ng/ml (1:100,000) in DMEM/FBS.
  - 1X control solution: Dilute DMSO 1:1,000 in DMEM/FBS.
2. Remove medium from wells that will be treated with either 1X induction or 1X control solutions.
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 17 hours.

#### Day 4: Analyze Luciferase Activity

1. Analyze luciferase activity using the Bright-Glo<sup>™</sup> Luciferase Assay System as described in Technical Manual #TM052.
2. Calculate the fold induction as follows:

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