

Promoter Analysis Tools

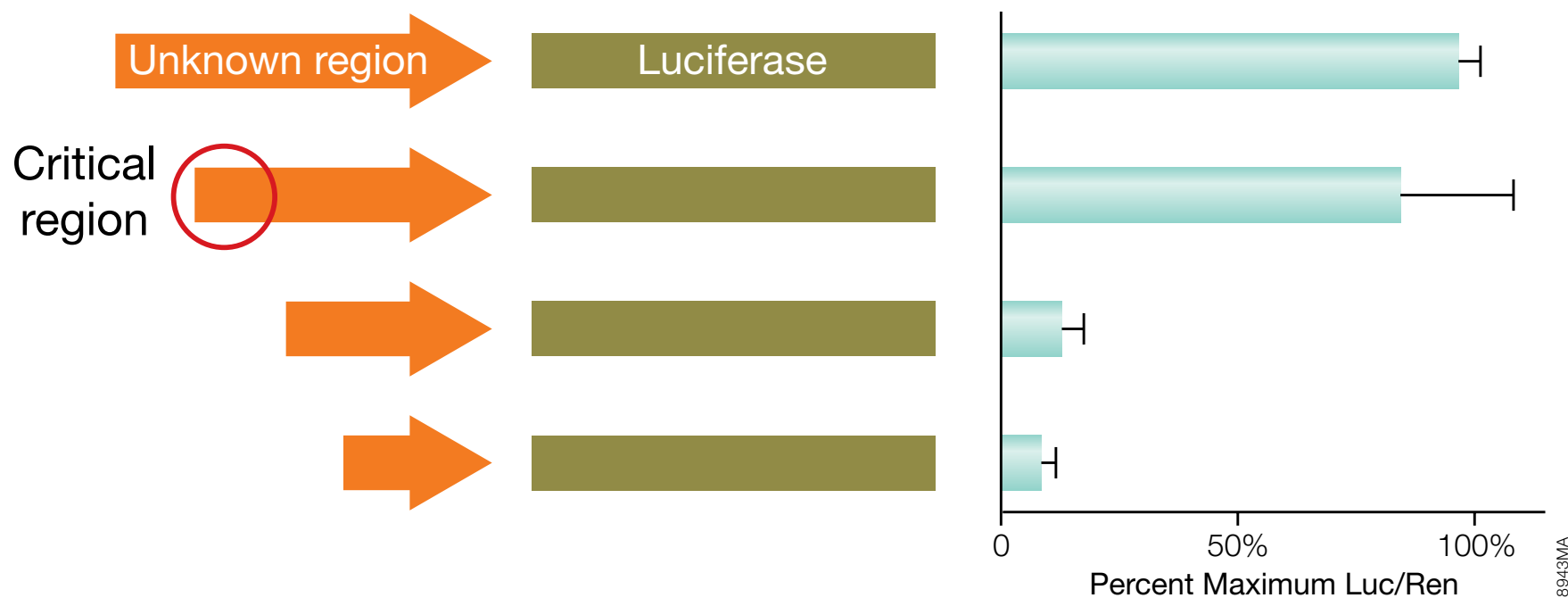
pGL4 Luciferase Vectors, Dual-Luciferase[®]
Assays and more...

*Harnessing the power
of bioluminescence to
understand cellular physiology*

Tap screen to start

Traditional Promoter Dissection

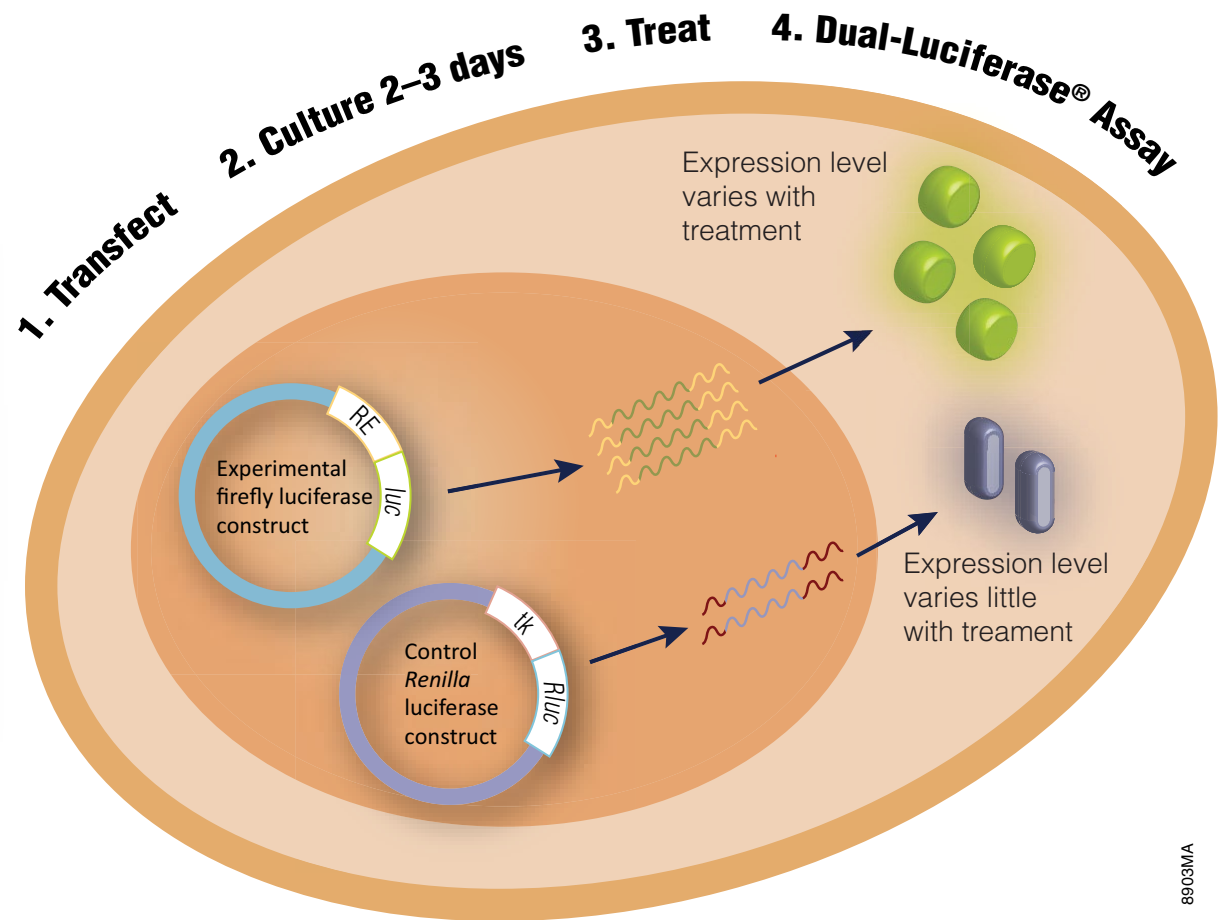
Deletion mutagenesis can be used to find functional elements responsible for the regulation of gene transcription. The cloned promoter is inserted upstream of the luciferase gene and the resulting construct transfected into cells to determine the function of the sequence.



Commonly called “promoter bashing”

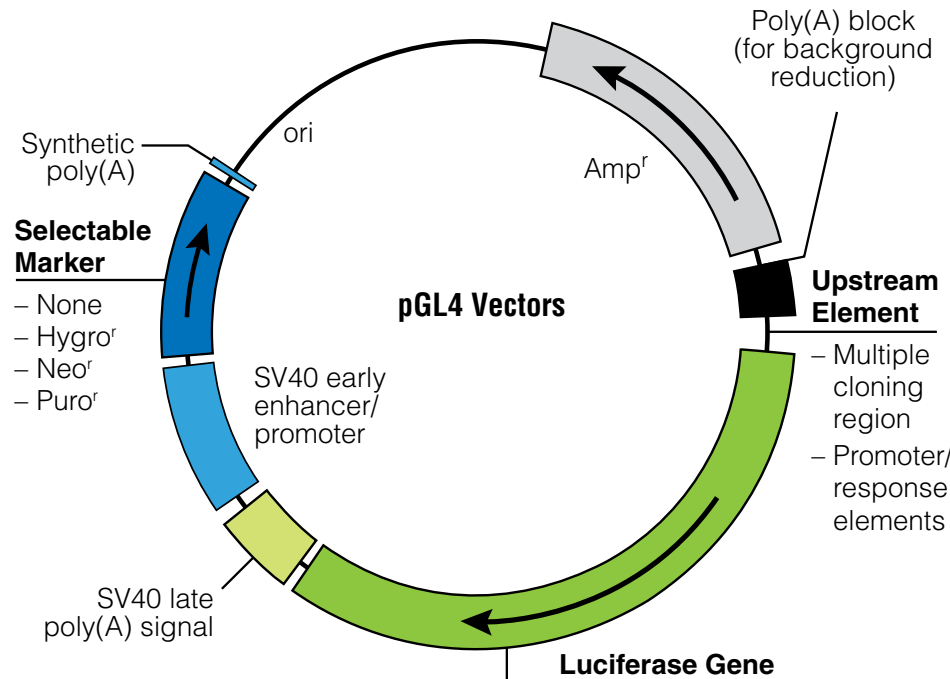
Assay Principle

- In a Dual-Luciferase[®] Assay, both the experimental construct with the promoter of interest and a control construct are introduced into the cell
- The *Renilla* control vector provides a consistent expression of *hRLuc* for normalization of the Firefly luciferase experimental vector
- The *hRLuc* activity will normalize for conditions of poor transfection efficiency and/or cytotoxicity
- Firefly and *Renilla* luciferase expression is easily assayed with the Dual-Luciferase[®] or Dual-Glo[®] Luciferase Assay Systems
- The roles of the firefly and *Renilla* luciferases can be reversed.

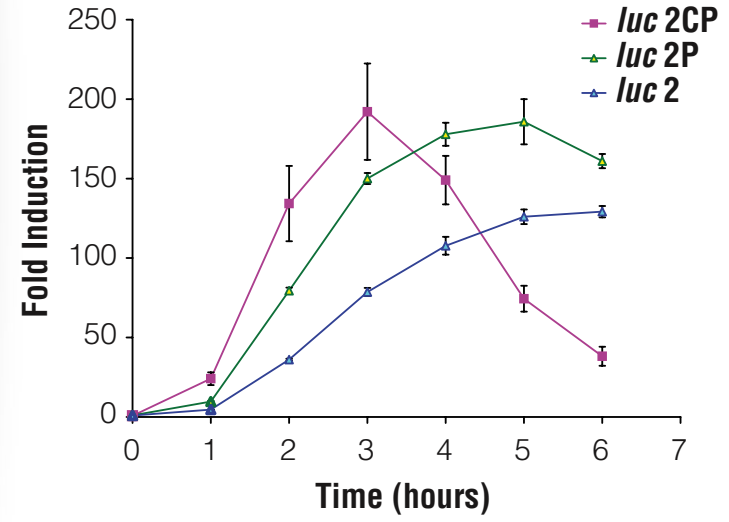


Experimental pGL4 Firefly Luciferase Vectors

Choose the functionality you need in your vector



pGL4 Vector backbone was redesigned to remove 75% of cryptic transcription factor binding sites in the pGL3 Vector backbone.



Luc2 Firefly Luciferase is the best

- Codon optimized for mammalian cells
- Removal of 90% of cryptic transcription factor binding sites found in *Luc+* of pGL3

[pGL4 Technical Manual](#) 

[online pGL4 Vector Selector](#)

RapidResponse™ *Luc2* genes have shortened half-lives and the shortened half-lives allow closer coupling to triggering event.

4861MB

Experimental pGL4 Firefly Luciferase Vectors

| Vector | MCS | Luciferase Gene | Selectable Marker | Cat. # |
|---|-----|-----------------|-------------------|--------|
| pGL4.10 [luc2] Vector | Y | <i>luc2</i> | No | E6651 |
| pGL4.11 [luc2P] Vector | Y | <i>luc2P</i> | No | E6661 |
| pGL4.12 [luc2CP] Vector | Y | <i>luc2CP</i> | No | E6671 |
| pGL4.14 [luc2/Hygro] Vector | Y | <i>luc2</i> | Hygromycin | E6691 |
| pGL4.15 [luc2P/Hygro] Vector | Y | <i>luc2P</i> | Hygromycin | E6701 |
| pGL4.16 [luc2CP/Hygro] Vector | Y | <i>luc2CP</i> | Hygromycin | E6711 |
| pGL4.17 [luc2/Neo] Vector | Y | <i>luc2</i> | Neomycin | E6721 |
| pGL4.18 [luc2P/Neo] Vector | Y | <i>luc2P</i> | Neomycin | E6731 |
| pGL4.19 [luc2CP/Neo] Vector | Y | <i>luc2CP</i> | Neomycin | E6741 |
| pGL4.20 [luc2/Puro] Vector | Y | <i>luc2</i> | Puromycin | E6751 |
| pGL4.21 [luc2P/Puro] Vector | Y | <i>luc2P</i> | Puromycin | E6761 |
| pGL4.22 [luc2CP/Puro] Vector | Y | <i>luc2CP</i> | Puromycin | E6771 |
| Constitutively Expressed | | | | |
| pGL4.13 [luc2/SV40] Vector | N | <i>luc2</i> | No | E6681 |
| pGL4.50 [luc2/CMV/Hygro] Vector | N | <i>luc2</i> | Hygromycin | E1310 |
| pGL4.51 [luc2/CMV/Neo] Vector | N | <i>luc2</i> | Neomycin | E1320 |

[Renilla Luciferase Control Vectors
Ordering Information >](#)



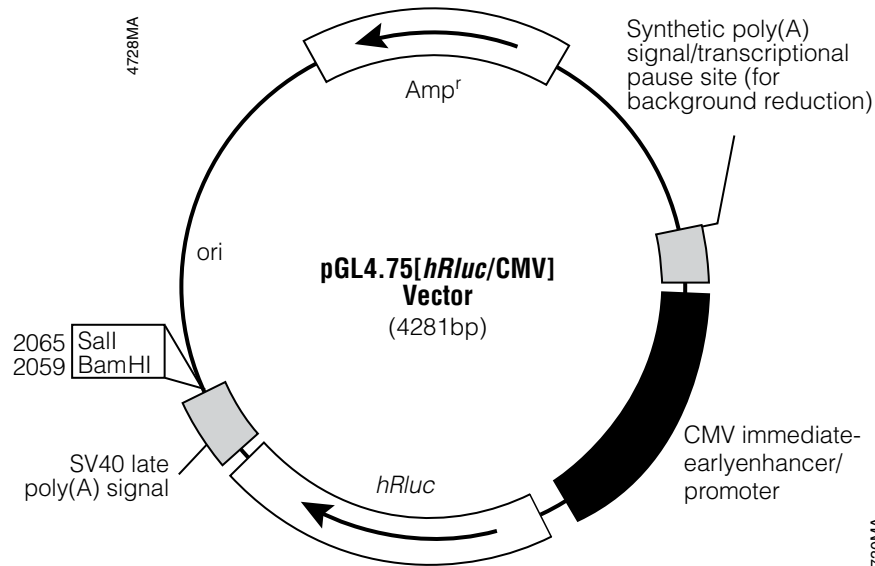
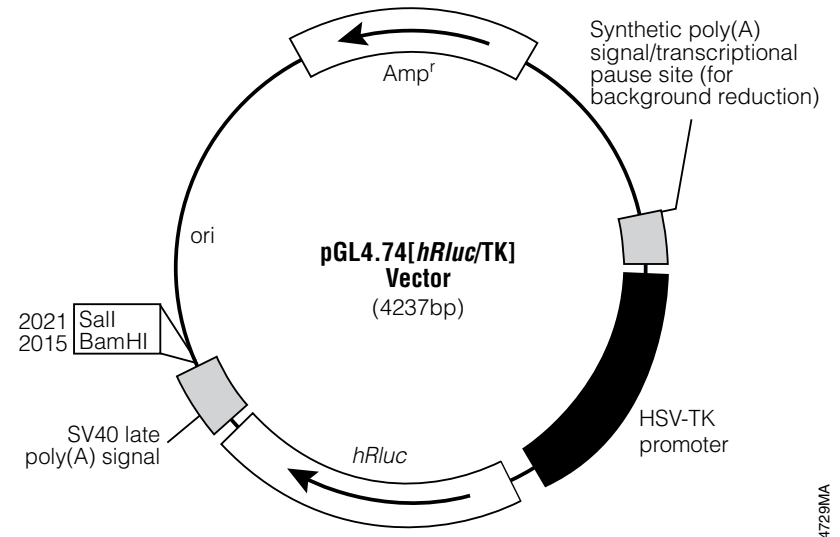
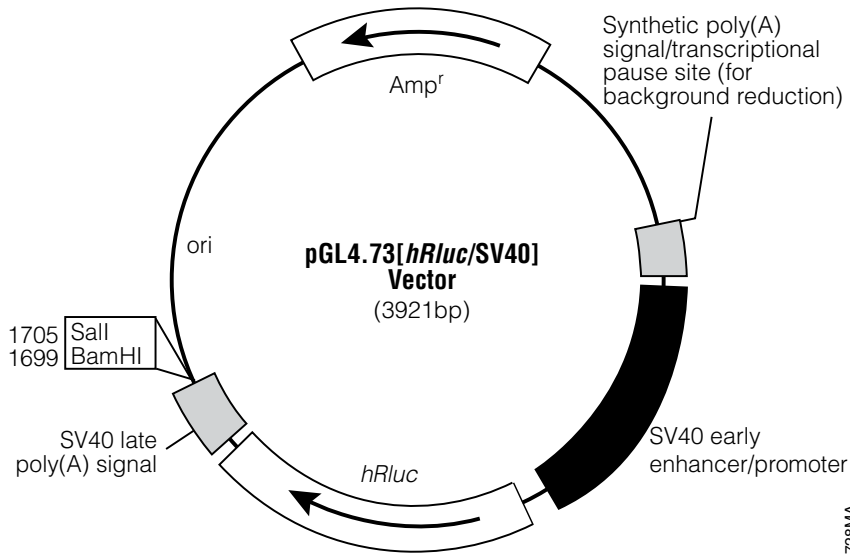
TechServ@promega.com



www.promega.com

Renilla Luciferase Control Vectors

The best companions for normalizing firefly luciferase reporters



hRLuc Improvements

- Codon optimized for mammalian cells
- Removal of 95% of cryptic transcription factor binding sites found in *Rluc* of pRL

[pGL4 Technical Manual](#) 

[online pGL4 Vector Selector](#)

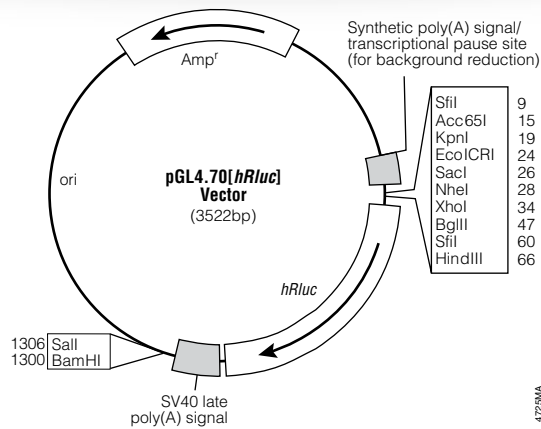
Renilla Luciferase Control Vectors

| Vector | MCS | Luciferase Gene | Selectable Marker | Cat. # |
|---|-----|-----------------|-------------------|--------|
| pGL4.73 [hRluc/SV40] Vector | N | <i>hRluc</i> | No | E6911 |
| pGL4.74 [hRluc/TK] Vector | N | <i>hRluc</i> | No | E6921 |
| pGL4.75 [hRluc/CMV] Vector | N | <i>hRluc</i> | No | E6931 |

Promoterless vectors with RapidResponse™ Gene options

| | | | | |
|--|---|----------------|----|-------|
| pGL4.70 [hRluc] Vector | Y | <i>hRluc</i> | No | E6881 |
| pGL4.71 [hRlucP] Vector | Y | <i>hRlucP</i> | No | E6891 |
| pGL4.72 [hRlucCP] Vector | Y | <i>hRlucCP</i> | No | E6901 |

Products may be covered by pending or issued patents or may have certain limitations. Please visit www.promega.com for more information. The method of recombinant Coleoptera luciferases is covered by U.S. patent Nos. 5,583,024; 5,674,713; and 5,700,673.



Experimental pGL4 Firefly Luciferase Vectors
[Ordering Information >](#)



TechServ@promega.com



www.promega.com

Related Products

| Product | Size | Cat. # |
|---|------------|--------|
| Dual Reporter Assays (larger sizes available) | | |
| Dual-Luciferase® Reporter Assay System 5 Step assay requiring lysate production. Use in multi-well plates requires dual-injectors | 100 assays | E1910 |
| Dual-Glo® Luciferase Assay System 2 step assay that lyses cells directly. Use in multiwell plates does not require injectors. | 10ml | E2920 |
| Rapid, Transfection-Grade Plasmid Preps | | |
| PureYield™ Plasmid Miniprep System | 100 preps | A1223 |
| PureYield™ Plasmid Midiprep System | 25 preps | A2492 |
| PureYield™ Plasmid Maxiprep System | 10 preps | A2392 |
| Transfection Reagent | | |
| FuGENE® HD Transfection Reagent* | 1ml | E2311 |
| | 5x1ml | E2312 |

Dual-Glo, Dual-Luciferase and GloMax are registered trademarks; GloResponse, RapidResponse and PureYield are trademarks of Promega Corporation. HighWire Press is a registered trademark of the Board of Trustees of the Leland Stanford Junior University

* FuGENE HD is sold only for research use at non-profit entities. See terms of use at www.promega.com/lul



Introduction to Reporter Gene Assays Animation

Online pGL4 Vector Selector

GloMax® Multi+ Detection System

Need a luminometer?

Go to www.promega.com/glomax to learn more and request a demo



References

Kopish, K. (2007) Deciphering the pGL4 vector code. *Promega Notes* **96**, 6-7.



Paguio, A., et al. (2005) pGL4 Vectors: A new generation of luciferase reporter vectors. *Promega Notes* **89**, 7-10.



Almond, B.D., et al. (2004) Introducing the Rapid Response™ Reporter Vectors. *Promega Notes* **87**, 18-22.



Zhuang, Y., et al. (2001) New synthetic Renilla gene and assay system increase expression, reliability and sensitivity. *Promega Notes* **79**, 6-11.



Schagat, T., Paguio, A. and Kopish, K. (2007) Normalizing genetic reporter assays: Approaches and considerations for increasing consistency and statistical significance. *Cell Notes* **17**, 9-12.



Allard, S.T.M. (2008) Bioluminescent reporter genes. *eNotes/fe0030*.



Protocols & Applications Guide: Bioluminescent Reporters

Citations

pGL4.10-pGL4.12 HighWire Press®

pGL4.14-pGL4.22 HighWire Press®

pGL4.70-pGL4.72 HighWire Press®

pGL4.73-pGL4.75 HighWire Press®



TechServ@promega.com



www.promega.com