

## pGL2 Luciferase Reporter Vectors

INSTRUCTIONS FOR USE OF PRODUCTS E1611, E1621, E1631 AND E1641.

## pGL2 Luciferase Reporter Vectors

All technical literature is available on the Internet at: [www.promega.com/tbs/](http://www.promega.com/tbs/)  
Please visit the web site to verify that you are using the most current version of this Technical Manual. Please contact Promega Technical Services if you have questions on use of this system. E-mail: [techserv@promega.com](mailto:techserv@promega.com)

1. Description .....	2
2. Product Components and Storage Conditions .....	2
3. pGL2 Vector Maps and Sequence Reference Points .....	3
A. pGL2-Basic Vector .....	3
B. pGL2-Promoter Vector .....	4
C. pGL2-Enhancer Vector .....	5
D. pGL2-Control Vector .....	6
4. Vector Manipulation .....	7
A. Cloning Strategies .....	7
B. Mapping Genetic Elements Located within DNA Fragments .....	8
C. Sequencing .....	9
5. Transfection of Eukaryotic Cells .....	9
6. Assays of Luciferase Activity .....	10
7. Appendix .....	12
A. Structure and Function .....	12
B. pGL3 and pGL4 Vectors .....	12
C. pGL2-Basic Vector Restriction Sites .....	13
D. pGL2-Promoter Vector Restriction Sites .....	15
E. pGL2-Enhancer Vector Restriction Sites .....	18
F. pGL2-Control Vector Restriction Sites .....	20
G. References .....	22
H. Related Products .....	23

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

## 1. Description

The pGL2 Luciferase Reporter Vectors<sup>(a)</sup> are designed for the quantitative analysis of factors that potentially regulate mammalian gene expression. These factors may be *cis*-acting, such as promoters and enhancers, or *trans*-acting, such as various DNA-binding factors. The pGL2 Vectors carry the coding region for wildtype firefly (*Photinus pyralis*) luciferase, used to monitor transcriptional activity in transfected eukaryotic cells. The assay of this genetic reporter is rapid, sensitive and quantitative. In addition, the pGL2 Vectors contain numerous features that aid the characterization and mutagenesis of the putative regulatory sequences.

## 2. Product Components and Storage Conditions

Product	Size	Cat.#
pGL2-Control Vector	20µg	E1611
pGL2-Enhancer Vector	20µg	E1621
pGL2-Promoter Vector	20µg	E1631
pGL2-Basic Vector	20µg	E1641

Information on related products, including the Luciferase Assay System, is provided in Sections 5, 6 and 7.H.

**Storage Conditions:** Store the vector DNA at -20°C.

## 3. pGL2 Vector Maps and Sequence Reference Points

This listings of restriction sites for the pGL2 Luciferase Reporter Vectors are provided in Sections 7.C-F.

### 3.A. pGL2 Basic Vector

The pGL2-Basic Vector lacks eukaryotic promoter and enhancer sequences, allowing maximum flexibility in cloning putative regulatory sequences. Expression of luciferase activity in cells transfected with this plasmid depends on insertion and proper orientation of a functional promoter upstream from *luc*. Potential enhancer elements can also be inserted upstream of the promoter or in the BamHI or Sall sites downstream of the luciferase gene.

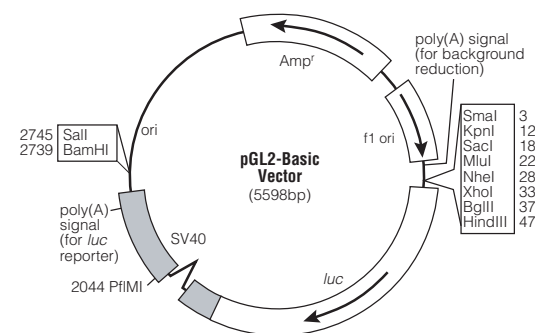


Figure 1. pGL2-Basic Vector map.

#### pGL2-Basic Vector Sequence Reference Points:

SV40 Promoter	(none)
SV40 Enhancer	(none)
Multiple cloning region	1-53
GLprimer2 binding site	77-99
Luciferase gene ( <i>luc</i> )	76-1728
SV40 late poly(A) signal	2518-2739
RVprimer4 binding site	2796-2815
β-lactamase ( <i>Amp<sup>r</sup></i> ) gene	3815-4675
ColE1-derived plasmid replication origin	3053
f1 origin	4807-5262
GLprimer1 binding site	5565-5587

### 3.B. pGL2-Promoter Vector

The pGL2-Promoter Vector contains the SV40 promoter upstream of the luciferase gene. DNA fragments containing putative enhancer elements can be inserted in either orientation and upstream or downstream of the promoter-*luc* transcriptional unit.

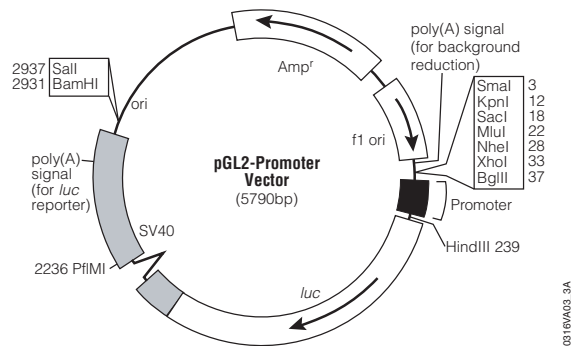


Figure 2. pGL2-Promoter Vector map.

#### pGL2-Promoter Vector Sequence Reference Points:

Multiple cloning region	1-49
SV40 Promoter	42-244
Luciferase gene ( <i>luc</i> )	268-1920
GLprimer2 binding site	269-291
RVprimer4 binding site	2988-3007
ColE1-derived plasmid replication origin	3245
β-lactamase ( <i>Amp<sup>r</sup></i> ) gene	4007-4867
f1 origin	4999-5454
GLprimer1 binding site	5757-5779

### 3.C. pGL2-Enhancer Vector

The pGL2-Enhancer Vector contains an SV40 enhancer, derived from the SV40 early promoter, located downstream of the luciferase gene after the poly(A) signal. This aids in the verification of functional promoter elements because the presence of the enhancer in many cases will result in transcription of *luc* at higher levels.

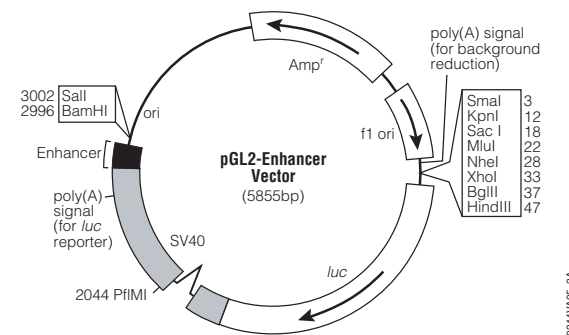


Figure 3. pGL2-Enhancer Vector map.

#### pGL2-Enhancer Vector Sequence Reference Points:

Multiple cloning region	1-52
Luciferase gene ( <i>luc</i> )	76-1728
GLprimer2 binding site	77-99
SV40 late poly(A) signal	2518-2739
SV40 Enhancer	2748-2984
RVprimer4 binding site	3053-3072
ColE1-derived plasmid replication origin	3310
β-lactamase ( <i>Amp<sup>r</sup></i> ) gene	4072-4932
f1 origin	5064-5519
GLprimer1 binding site	5822-5844



The SmaI/XmaI site upstream of the restriction sites mentioned above can be used to insert other putative regulatory elements without compromising the ability to make nested deletions. It can also be used to aid in mapping nested deletions.

Because the pGL2 Vectors are supplied as modified DNA, *E. coli* hosts may be either restriction+ or restriction-. The use of a *recA* host such as JM109 is preferred because this prevents undesirable recombination between the insert and the host chromosomal DNA. A strain that has an F' episome is required for ssDNA production. Grow JM109 on minimal plates (M-9) supplemented with thiamine-HCl prior to preparation of competent cells and transformation. This selects for the presence of the F' episome.

Protocols for restriction digestion, alkaline phosphatase treatment, linker ligation and transformation of competent cells can be found in reference 2.

#### 4.B. Mapping Genetic Elements Located within DNA Fragments

The locations of functional elements within a DNA fragment are often determined by making a set of unidirectional deletions following the method of Henikoff (3) and assaying for biological activity. This method takes advantage of the unique properties of Exonuclease III (ExoIII), which will digest 5'-overhangs, but not 3'-overhangs or  $\alpha$ -phosphorothioate nucleotide filled-in overhangs. Nested deletions of the insert can be made directly in the promoter regions of the pGL2-Promoter and pGL2-Control Luciferase Reporter Vectors using this method, eliminating the need for subcloning steps. The upstream MluI, NheI and XhoI sites, rare restriction sites in eukaryotic DNA, provide the potential 5'-starting points for ExoIII digestion at the upstream end of the insert; KpnI and SacI generate the 3'-overhangs resistant to ExoIII. After treatment with ExoIII, S1 nuclease is added to remove the resulting ssDNA, and T4 DNA ligase is added to reclose the vectors. Deletion clones can be screened by gel electrophoresis of miniprep DNA, and the precise deletion endpoints within the promoter region can be determined by DNA sequencing using primers designed for the pGL2 Vectors.

#### 4.C. Sequencing

For some applications, it may be desirable to sequence the DNA inserted into the pGL2 Vectors. Two examples of such applications are to determine the exact position of generated deletions (see Section 4.B) and to confirm production of a site-specific mutation. Two primers are available for this purpose: GLprimer1 for sequencing clockwise across the upstream cloning sites, and GLprimer2 for sequencing counterclockwise across the cloning sites.

GLprimer1 5'-TGTATCITTATGGTACTIGTAACTG-3'

GLprimer2 5'-CTTTATGTTTTGGCGCTTCCA-3'

GLprimer1 is especially useful for identifying positions of nested deletions. Note that both primers can be used for dsDNA sequencing, but GLprimer2 also may be used for ssDNA sequencing.

#### 5. Transfection of Eukaryotic Cells

Transfection of DNA into eukaryotic cells may be mediated by cationic lipid compounds (4,5), calcium phosphate (6,7), DEAE-dextran (6,8) or electroporation (7). Transfection systems based on cationic lipids (TransFast™ Transfection Reagent, Transfectam® Reagent and Tfx™ Reagents), calcium phosphate and DEAE-dextran (Profection® Mammalian Transfection Systems) are available from Promega. For more information on these transfection reagents, please request the *TransFast™ Transfection Reagent Technical Bulletin #TB260*, the *Transfectam® Reagent Technical Bulletin #TB116*, the *Tfx™-Reagents Technical Bulletin #TB216* or the *Profection® Mammalian Transfection Systems Technical Manual #TM012*. All of these documents are available on our web site at: [www.promega.com/tbs/](http://www.promega.com/tbs/)

**Note:** The specific transcriptional characteristics of the pGL2 Vectors are likely to vary for different cell types. This may be particularly true for COS cells that contain the SV40 large T antigen. The SV40 large T antigen promotes replication of the SV40 origin, found in the promoter region of the pGL2-Promoter and pGL2-Control Vectors. The combination of large T antigen and SV40 origin will result in a higher copy number of these vectors in COS cells, which in turn may result in increased expression of the reporter gene compared to expression levels from vectors lacking the SV40 origin.

## 6. Assays of Luciferase Activity

Experimental strategies using firefly luciferase may involve the analysis of a few samples per day or as many as several thousand samples per hour, and equipment used to measure luminescence may vary from inexpensive, single-sample luminometers to high-end CCD luminometers. To support this wide range of applications, Promega has developed three luciferase assays with different but complementary characteristics: Luciferase Assay System (Cat.# E1500), Bright-Glo™ Luciferase Assay System (Cat.# E2610), Steady-Glo® Luciferase Assay System (Cat.# E2510), and ONE-Glo™ Luciferase Assay System (Cat.# E6110). Reagent choice depends on the relative importance of experimental format, assay sensitivity and luminescence duration.

**Table 1. Characteristics of Promega Luciferase Assay Reagents.**

	Bright-Glo™ Reagent	Steady-Glo® Reagent	Luciferase Assay Reagent	ONE-Glo™ Reagent
<b>Format</b>	NH or H	NH or H	NH	NH or H
<b>Process</b>	continuous	batch	bench scale	batch or continuous
<b>Number of Steps</b>	1	1	4	1
<b>Sensitivity</b>	highest	lower	higher	high
<b>Signal Half-Life</b>	~30 minutes	~5 hours	~12 minutes	~50 minutes
<b>Precision</b>	High	High	High	Highest
<b>Cell Lysis Time</b>	~2 minutes maximum	~5 minutes maximum	NA	~3 minutes

NH = nonhomogeneous (first create a lysate); H = homogeneous; NA = not applicable

The Luciferase Assay System has long been the standard reagent for routine laboratory analysis. Before using this reagent, cells from which the luciferase is to be measured must be washed and lysed. This reagent was optimized for high sensitivity in nonhomogeneous, single-sample measurements. The Luciferase Assay System requires a luminometer fitted with injectors to efficiently measure luminescence in 96-well plates.

The Bright-Glo™, Steady-Glo® and ONE-Glo™ Reagents were developed to perform assay reactions within multiwell plates and in the presence of complete cell culture medium: no cell preparation steps such as washing or lysing are required before the luminescence reaction is initiated. Both of these are single-step reagents, requiring only addition of the reagent before measuring luminescence. This makes them ideal reagents for efficient and precise quantitation in 96-, 384- and 1536-well plates.

The Bright-Glo™ and Steady-Glo® Reagents are complementary in their characteristics based on the inverse relationship between luminescence duration and assay sensitivity (9). Generally as the half-life of the luminescence increases, assay sensitivity decreases. The Steady-Glo® Reagent provides long

luminescence duration (changing only about 10% per hour); however, to achieve this long luminescence duration, the assay sensitivity must be reduced. This reagent was designed for experiments in which many microplates are processed as a batch.

In contrast, the Bright-Glo™ Reagent provides high assay sensitivity with shorter luminescence duration (<10% decrease per 5 minutes). This reagent is designed for general research applications and for experiments using robotics for continuous sample processing. Furthermore, as a result of increased sample capacity, the Bright-Glo™ Reagent provides greater assay sensitivity than the Luciferase Assay Reagent in most applications (9).

The ONE-Glo™ Reagent provides the ultimate performance for luciferase assays. It features a high-sensitivity assay with extended duration. The ONE-Glo™ Reagent also demonstrates more robust performance and provides reagent handling enhancements.

The Luciferase Assay System, Bright-Glo™ Reagent and Steady-Glo® Reagent and ONE-Glo™ Reagent provide the highest standards in assay quantitation, sensitivity and convenience. Since these reagents are based on the same underlying design principles, different reagents can be used as experimental needs change. For more information request the *Luciferase Assay System Technical Bulletin #TB281*, the *Steady-Glo® Luciferase Assay System Technical Manual #TM051*, the *Bright-Glo™ Luciferase Assay System Technical Manual #TM052*, or the *ONE-Glo™ Luciferase Assay System Technical Manual #TM292*.

When studying promoter functionalities, it is often desirable to include a second reporter (e.g., *Renilla* luciferase) as an internal control for normalization. Plasmids derived from pGL2, pGL3 or pGL4 vectors can be co-transfected with *Renilla* luciferase vectors, such as phRL-TK, and assayed using the Dual-Luciferase® Reporter Assay System (Cat.# E1910) or the Dual-Glo® Luciferase Assay System (Cat.# E2920).

**Table 2. Characteristics of Promega Dual-Luciferase Assays.**

	Dual-Luciferase® Assay	Dual-Glo® Assay
<b>Format</b>	NH	H
<b>Process</b>	bench scale	batch
<b>Number of Steps</b>	5	2
<b>Sensitivity</b>	higher	lower
<b>Signal Half-Life – firefly</b>	~9 minutes	~2 hours
<b>Signal Half-Life – Renilla</b>	~2 minutes	~2 hours
<b>Precision</b>	high	high
<b>Cell Lysis Time</b>	~10 minutes	~15 minutes

NH = nonhomogeneous (lysis step not part of assay);

H = homogeneous (lysis step included in assay)

## 7. Appendix

### 7.A. Structure and Function

Promega supplies four different pGL2 Vectors: pGL2-Basic, pGL2-Promoter, pGL2-Enhancer and pGL2-Control. Each vector carries the luciferase gene (*luc*) followed by the SV40 small t antigen intron and early polyadenylation [poly(A)] signals. A second copy of the poly(A) site, located upstream of the luciferase coding region, limits background transcription from spurious promoters in the plasmids. The vectors also contain a high copy number prokaryotic origin of replication for maintenance in *E. coli*, an ampicillin-resistance gene for selection and a filamentous phage origin of replication (f1 ori) for single-stranded DNA (ssDNA) production. Restriction sites for insertion of DNA fragments are located upstream and downstream of the luciferase gene. Two of the upstream sites (XhoI and BglIII) yield cohesive ends compatible with the downstream sites (Sall and BamHI, respectively), allowing the interchange of the DNA insert for rapid analysis of positional effects. Except for the inclusion of promoters and enhancers, the four pGL2 Vectors are structurally identical. Specific features of each plasmid are summarized in Section 3. Maps of the pGL2-Basic, pGL2-Promoter, pGL2-Enhancer and pGL2-Control Vectors are shown in Figures 1-5. The list of restriction sites for these vectors is provided in Section 7.

### 7.B. pGL3 and pGL4 Vectors

Promega also provides two improved series of luciferase reporter vectors, the pGL3 Luciferase Reporter Vectors and pGL4 Luciferase Reporter Vectors. The pGL3 Vectors provide significant advances over the pGL2 Vectors (1). The pGL3 Vectors contain a modified firefly luciferase cDNA, designated *luc+*, and a redesigned vector backbone. These changes were made to increase luciferase expression, improve in vivo vector stability, and provide greater flexibility in performing genetic manipulations. The modified reporter vectors have resulted in luciferase expression levels dramatically higher than those obtained with the pGL2 Reporter Vectors, while maintaining relatively low background luciferase expression. For further information, see the *pGL3 Luciferase Reporter Vectors Technical Manual #TM033* available at: [www.promega.com/tbs/](http://www.promega.com/tbs/)

The pGL4 Luciferase Reporter Vectors are the latest generation of reporter gene vectors optimized for expression in mammalian cells. Numerous configurations of pGL4 Vectors are available, including those with the synthetic firefly *luc2* (*Photinus pyralis*) and *Renilla hRLuc* (*Renilla reniformis*) genes, which have been codon optimized for more efficient expression in mammalian cells. Furthermore, both the reporter genes and the vector backbone, including the *bla* ( $\beta$ -lactamase or Amp<sup>r</sup>) and mammalian selectable marker genes for hygromycin (Hygro or Hyg<sup>r</sup>), neomycin (Neo or Neo<sup>r</sup>) and puromycin (Puro or Puro<sup>r</sup>), have been engineered to reduce the number of consensus transcription factor binding sites, reducing background and the risk of anomalous transcription. For more information, see the *pGL4 Luciferase Reporter Vectors Technical Manual #TM259* available at: [www.promega.com/tbs/](http://www.promega.com/tbs/)

### 7.C. pGL2-Basic Vector Restriction Sites

The following restriction enzyme tables were constructed using DNASTAR<sup>®</sup> sequence analysis software. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3' end of the cut DNA (the base to the left of the cut site). For more information on the cut sites of these enzymes, or if you identify a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. The pGL2-Basic Vector sequence is also available in the GenBank<sup>®</sup> database (GenBank<sup>®</sup>/EMBL Accession Number **X65323**) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Table 3. Restriction Enzymes That Cut the pGL2-Basic Vector 1-5 Times.**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
<b>AccB7I</b>	1	2044	BssSI	2	3168, 4552
<b>AccI</b>	1	2746	<b>BstEII</b>	1	683
<b>AccIII</b>	2	771, 1287	<b>Bsu36I</b>	1	689
<b>Acc65I</b>	1	8	Cfr10I	5	345, 1504, 2863, 3968, 4932
AcyI	4	83, 109, 1502, 4425	<b>ClaI</b>	1	1441
AflIII	3	22, 569, 2995	<b>Csp45I</b>	2	245, 1033
<b>Alw26I</b>	4	1099, 1331, 3949, 4725	DraII	1	1255
<b>Alw44I</b>	2	3309, 4555	DraIII	1	5040
AlwNI	1	3411	DrdI	2	3103, 5084
AspHI	5	18, 1541, 3313, 4474, 4559	EaeI	1	4276
<b>AvaI</b>	3	1, 33, 1132	<b>EclHKI</b>	1	3888
<b>AvaII</b>	3	1255, 4026, 4248	<b>Eco47III</b>	1	2871
<b>BamHI</b>	1	2739	Eco81I	1	689
<b>BanII</b>	3	18, 1100, 4966	<b>EcoICRI</b>	1	16
BbeI	1	112	EcoNI	1	1693
BbsI	5	86, 1364, 1480, 2192, 2824	<b>EcoRI</b>	1	663
<b>BbuI</b>	1	739	<b>EcoRV</b>	1	1414
<b>BglI</b>	2	4008, 5276	EheI	1	110
<b>BglIII</b>	1	37	FspI	2	4110, 5283
BsaI	1	3949	<b>HaeII</b>	5	112, 2873, 3243, 4882, 4890
BsaAI	3	230, 5037, 5370	<b>HincII</b>	4	1380, 2608, 2747, 5451
BsaBI	1	2507	HindII	4	1380, 2608, 2747, 5451
BsaHI	4	83, 109, 1502, 4425	<b>HindIII</b>	1	47
BspHI	2	3715, 4723	<b>HpaI</b>	2	2608, 5451
BspMI	1	1474	<b>Hsp92I</b>	4	83, 109, 1502, 4425
BsrGI	1	566	KasI	1	108

**Note:** The enzymes listed in boldface type are available from Promega.

### 7.C. pGL2-Basic Vector Restriction Sites (continued)

**Table 3. Restriction Enzymes That Cut the pGL2-Basic Vector 1-5 Times (continued).**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
KpnI	1	12	SacI	1	18
MluI	1	22	SalI	1	2745
MspAII	5	139, 1429, 3337, 3582, 4523	ScaI	1	4368
NaeI	2	2865, 4934	SgrAI	1	1504
NarI	1	109	SinI	3	1255, 4026, 4248
NgoMIV	2	2863, 4932	SmaI	1	3
NheI	1	28	SphI	1	739
NspI	2	739, 2999	SplI	1	230
PacI	1	1400	SspI	5	2323, 2476, 4692, 5245, 5360
Paer7I	1	33	StyI	1	2192
PfiMI	1	2044	VspI	2	2411, 4060
PpuMI	1	1255	XbaI	1	123
PshAI	1	2810	XcmI	1	811
Psp5II	1	1255	XhoI	1	33
PspAI	1	1	XmaI	1	1
PvuI	2	4258, 5304	XmnI	1	4487

**Table 4. Restriction Enzymes That Do Not Cut the pGL2-Basic Vector.**

AatII	Bpu1102I	Eco52I	PmeI	SpeI
AflIII	Bsp120I	Eco72I	PmlI	SrfI
AgeI	BssHII	FseI	Ppu10I	Sse8387I
ApaI	Bst1107I	<b>I-PpoI</b>	PstI	StuI
AscI	<b>Bst98I</b>	NcoI	PvuII	SwaI
AvrII	<b>BstXI</b>	NdeI	RsrII	Tth111I
BalI	<b>BstZI</b>	NotI	SacII	
BbrPI	<b>CspI</b>	NruI	SfiI	
BclI	DsaI	NsiI	Sgfl	
BplI	EagI	PinAI	<b>SnaBI</b>	

**Note:** The enzymes listed in boldface type are available from Promega.

**Table 5. Restriction Enzymes That Cut the pGL2-Basic Vector Six or More Times.**

AcI	Bst71I	HaeIII	MboII	Sau96I
AluI	<b>BstOI</b>	HgaI	MnlI	ScrFI
BanI	BstUI	<b>HhaI</b>	MseI	SfaNI
BbvI	<b>CfoI</b>	<b>HinfI</b>	<b>MspI</b>	<b>TaqI</b>
BsaOI	<b>DdeI</b>	<b>HpaII</b>	<b>NciI</b>	TfiI
BsaJI	<b>DpnI</b>	HphI	<b>NdeII</b>	<b>Tru9I</b>
<b>BsaMI</b>	DpnII	<b>Hsp92II</b>	NlaIII	<b>XhoII</b>
BsmI	<b>DraI</b>	MaeI	NlaIV	
<b>Bsp1286I</b>	EarI	MaeII	PleI	
BsrI	Fnu4HI	MaeIII	<b>RsaI</b>	
<b>BsrSI</b>	<b>FokI</b>	<b>MboI</b>	<b>Sau3AI</b>	

### 7.D. pGL2-Promoter Vector Restriction Sites

The following restriction enzyme tables were constructed using DNASTAR® sequence analysis software. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3' end of the cut DNA (the base to the left of the cut site). For more information on the cut sites of these enzymes, or if you identify a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. The pGL2-Promoter Vector sequence is also available in the GenBank® database (GenBank®/EMBL Accession Number **X65326**) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Table 6. Restriction Enzymes That Cut the pGL2-Promoter Vector 1-5 Times.**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccB7I	1	2236	BbsI	5	278, 1556, 1672, 2384, 3016
AccI	1	2938	<b>BbuI</b>	1	931
AccIII	2	963, 1479	<b>BglI</b>	3	176, 4200, 5468
Acc65I	1	8	<b>BglII</b>	1	37
AcyI	4	275, 301, 1694, 4617	BsaI	1	4141
AflIII	3	22, 761, 3187	BsaAI	3	422, 5229, 5562
Alw26I	4	1291, 1523, 4141, 4917	BsaBI	2	42, 2699
<b>Alw44I</b>	2	3501, 4747	BsaHI	4	275, 301, 1694, 4617
AlwNI	1	3603	BspHI	2	3907, 4915
AspHI	5	18, 1733, 3505, 4666, 4751	BspMI	1	1666
<b>AvaI</b>	3	1, 33, 1324	BsrGI	1	758
<b>AvaII</b>	3	1447, 4218, 4440	BssSI	2	3360, 4744
AvrII	1	223	<b>BstEII</b>	1	875
<b>BamHI</b>	1	2931	<b>Bsu36I</b>	1	881
<b>BanII</b>	3	18, 1292, 5158	Cfr10I	5	537, 1696, 3055, 4160, 5124
BbeI	1	304	<b>ClaI</b>	1	1633

7.D. pGL2-Promoter Vector Restriction Sites (continued)

Table 6. Restriction Enzymes That Cut the pGL2-Promoter Vector 1-5 Times (continued).

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
<b>Csp45I</b>	2	437, 1225	<b>NcoI</b>	1	130
<b>DraII</b>	1	1447	<b>NgoMIV</b>	2	3055, 5124
<b>DraIII</b>	1	5232	<b>NheI</b>	1	28
<b>DrdI</b>	2	3295, 5276	<b>NspI</b>	2	931, 3191
<b>DsaI</b>	1	130	<b>Pacl</b>	1	1592
<b>EaeI</b>	1	4468	<b>PaeR7I</b>	1	33
<b>EclHKI</b>	1	4080	<b>PflMI</b>	1	2236
<b>Eco47III</b>	1	3063	<b>PpuMI</b>	1	1447
<b>Eco81I</b>	1	881	<b>PshAI</b>	1	3002
<b>EcoICRI</b>	1	16	<b>Psp5II</b>	1	1447
<b>EcoNI</b>	1	1885	<b>PspAI</b>	1	1
<b>EcoRI</b>	1	855	<b>PvuI</b>	2	4450, 5496
<b>EcoRV</b>	1	1606	<b>SacI</b>	1	18
<b>EheI</b>	1	302	<b>SalI</b>	1	2937
<b>FspI</b>	2	4302, 5475	<b>ScaI</b>	1	4560
<b>HaeII</b>	5	304, 3065, 3435, 5074, 5082	<b>SfiI</b>	1	176
<b>HincII</b>	4	1572, 2800, 2939, 5643	<b>SgrAI</b>	1	1696
<b>HindII</b>	4	1572, 2800, 2939, 5643	<b>SinI</b>	3	1447, 4218, 4440
<b>HindIII</b>	1	239	<b>SmaI</b>	1	3
<b>HpaI</b>	2	2800, 5643	<b>SphI</b>	1	931
<b>Hsp92I</b>	4	275, 301, 1694, 4617	<b>SpII</b>	1	422
<b>KasI</b>	1	300	<b>SspI</b>	5	2515, 2668, 4884, 5437, 5552
<b>KpnI</b>	1	12	<b>StuI</b>	1	222
<b>MluI</b>	1	22	<b>StyI</b>	3	130, 223, 2384
<b>MspAII</b>	5	331, 1621, 3529, 3774, 4715	<b>VspI</b>	2	2603, 4252
<b>NaeI</b>	2	3057, 5126	<b>XbaI</b>	1	315
<b>NarI</b>	1	301	<b>XcmI</b>	1	1003
			<b>XhoI</b>	1	33
			<b>XmaI</b>	1	1
			<b>XmnI</b>	1	4679

Table 7. Restriction Enzymes That Do Not Cut the pGL2-Promoter Vector.

<b>AatII</b>	Bpu1102I	Eco52I	PmeI	<b>SpeI</b>
<b>AflIII</b>	Bsp120I	Eco72I	PmlI	SrfI
<b>AgeI</b>	<b>BssHIII</b>	FseI	Ppu10I	Sse8387I
<b>ApaI</b>	Bst1107I	<b>I-PpoI</b>	<b>PstI</b>	Swal
<b>AscI</b>	<b>Bst98I</b>	<b>NdeI</b>	<b>PvuII</b>	<b>Tth111I</b>
<b>BalI</b>	<b>BstXI</b>	<b>NotI</b>	RsrII	
BbrPI	<b>BstZI</b>	<b>NruI</b>	<b>SacII</b>	
<b>BclI</b>	<b>CspI</b>	<b>NsiI</b>	<b>SgfI</b>	
<b>BlpI</b>	EagI	PinAI	<b>SnaBI</b>	

Table 8. Restriction Enzymes That Cut the pGL2-Promoter Vector More Than Six Times.

<b>Acil</b>	<b>Bst71I</b>	<b>HaeIII</b>	<b>MboII</b>	Sau96I
<b>AluI</b>	<b>BstOI</b>	HgaI	MnlI	ScrFI
<b>BanI</b>	BstUI	<b>HhaI</b>	MseI	SfaNI
BbvI	<b>CfoI</b>	<b>Hinfl</b>	<b>MspI</b>	<b>TaqI</b>
BsaOI	<b>DdeI</b>	<b>HpaII</b>	<b>NciI</b>	TfiI
BsaJI	<b>DpnI</b>	HphI	<b>NdeII</b>	<b>Tru9I</b>
<b>BsaMI</b>	DpnII	<b>Hsp92II</b>	NlaIII	<b>XhoII</b>
BsmI	<b>DraI</b>	MaeI	NlaIV	
<b>Bsp1286I</b>	EarI	MaeII	PleI	
BsrI	Fnu4HI	MaeIII	<b>RsaI</b>	
<b>BsrSI</b>	<b>FokI</b>	<b>MboI</b>	<b>Sau3AI</b>	

Note: The enzymes listed in boldface type are available from Promega.

### 7.E. pGL2-Enhancer Vector Restriction Sites

The following restriction enzyme tables were constructed using DNASTAR® sequence analysis software. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3' end of the cut DNA (the base to the left of the cut site). For more information on the cut sites of these enzymes, or if you identify a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. The pGL2-Enhancer Vector sequence is also available in the GenBank® database (GenBank®/EMBL Accession Number X65325) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Table 9. Restriction Enzymes That Cut the pGL2-Enhancer Vector 1-5 Times.**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccB7I	1	2044	Cfr10I	5	345, 1504, 3120, 4225, 5189
AccI	1	3003	ClaI	1	1441
AccIII	2	771, 1287	Csp45I	2	245, 1033
Acc65I	1	8	DraII	1	1255
AcyI	4	83,109,1502, 4682	DraIII	1	5297
AflIII	3	22, 569, 3252	DrdI	2	3360, 5341
Alw26I	4	1099, 1331, 4206, 4982	DsaI	1	2748
Alw44I	2	3566, 4812	EaeI	1	4533
AlwNI	1	3668	EclHKL	1	4145
AspHI	5	18, 1541, 3570, 4731, 4816	Eco47III	1	3128
AvaI	3	1, 33, 1132	Eco8II	1	689
AvaII	3	1255, 4283, 4505	EcoICRI	1	16
BamHI	1	2996	EcoNI	1	1693
BanII	3	18, 1100, 5223	EcoRI	1	663
BbeI	1	112	EcoRV	1	1414
BbsI	5	86, 1364, 1480, 2192, 3081	EheI	1	110
BbuI	3	739, 2843, 2915	FspI	2	4367, 5540
BglII	2	4265, 5533	HaeII	5	112, 3130, 3500, 5139, 5147
BglIII	1	37	HincII	4	1380, 2608, 3004, 5708
BsaI	1	4206	HindII	4	1380, 2608, 3004, 5708
BsaAI	3	230, 5294, 5627	HindIII	1	47
BsaBI	1	2507	HpaI	2	2608, 5708
BsaHI	4	83, 109,1502, 4682	Hsp92I	4	83, 109, 1502, 4682
BspHI	2	3972, 4980	KasI	1	108
BspMI	1	1474	KpnI	1	12
BsrGI	1	566	MluI	1	22
BssSI	2	3425, 4809	MspAII	5	139, 1429, 3594, 3839, 4780
BstEII	1	683			
Bsu36I	1	689			

**Table 9. Restriction Enzymes That Cut the pGL2-Enhancer Vector 1-5 Times (continued).**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
NaeI	2	3122, 5191	SacI	1	18
NarI	1	109	SalI	1	3002
NcoI	1	2748	Scal	1	4625
NgoMIV	2	3120, 5189	SgrAI	1	1504
NheI	1	28	SinI	3	1255, 4283, 4505
NsiI	2	2841, 2913	SmaI	1	3
NspI	4	739, 2843, 2915, 3256	SphI	3	739, 2843, 2915
PacI	1	1400	SplI	1	230
PaeR7I	1	33	SspI	5	2323, 2476, 4949, 5502, 5617
PflMI	1	2044	StyI	2	2192, 2748
Ppu10I	2	2837, 2909	VspI	2	2411, 4317
PpuMI	1	1255	XbaI	1	123
PshAI	1	3067	XcmI	1	811
Psp5II	1	1255	XhoI	1	33
PspAI	1	1	XmaI	1	1
PvuI	2	4515, 5561	XmnI	1	4744

**Table 10. Restriction Enzymes That Do Not Cut the pGL2-Enhancer Vector.**

AatII	BlpI	EagI	PmeI	SpeI
AflIII	Bpu1102I	Eco52I	PmlI	SrfI
AgeI	Bsp120I	Eco72I	PstI	Sse8387I
ApaI	BssHIII	FseI	PvuII	StuI
AscI	Bst1107I	I-PpoI	RsrII	Swal
AvrII	Bst98I	NdeI	SacII	Tth111I
BalI	BstXI	NotI	SfiI	
BbrPI	BstZI	NruI	SgfI	
BclI	CspI	PinAI	SnaBI	

**Table 11. Restriction Enzymes That Cut the pGL2-Enhancer Vector More Than Six Times.**

AcI	BsrSI	Fnu4HI	MaeII	PleI
AluI	Bst71I	FokI	MaeIII	RsaI
BanI	BstOI	HaeIII	MboI	Sau3AI
BbvI	BstUI	HgaI	MboII	Sau96I
BsaOI	CfoI	HhaI	MnlI	ScrFI
BsaJI	DdeI	Hinfl	MseI	SfaNI
BsaMI	DpnI	HpaII	MspI	TaqI
BsmI	DpnII	HphI	NciI	TfiI
Bsp1286I	DraI	Hsp92II	NdeII	Tru9I
BsrI	EarI	MaeI	NlaIII	XhoII

**Note:** The enzymes listed in boldface type are available from Promega.

### 7.F. pGL2-Control Vector Restriction Sites

The following restriction enzyme tables were constructed using DNASTAR® sequence analysis software. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3' end of the cut DNA (the base to the left of the cut site). For more information on the cut sites of these enzymes, or if you identify a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. The pGL2-Control Vector sequence is also available in the GenBank® database (GenBank®/EMBL Accession Number X65324) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Table 12. Restriction Enzymes That Cut the pGL2-Control Vector 1-5 Times.**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccB7I	1	2236	BstEII	1	875
AccI	1	3195	Bsu36I	1	881
AccIII	2	963, 1479	Cfr10I	5	537, 1696, 3312, 4417, 5381
Acc65I	1	8	ClaI	1	1633
AcyI	4	275, 301, 1694, 4874	Csp45I	2	437, 1225
AflIII	3	22, 761, 3444	DraII	1	1447
Alw26I	4	1291, 1523, 4398, 5174	DraIII	1	5489
Alw44I	2	3758, 5004	DrdI	2	3552, 5533
AlwNI	1	3860	DsaI	2	130, 2940
AspHI	5	18, 1733, 3762, 4923, 5008	EaeI	1	4725
AvaI	3	1, 33, 1324	EclHKL	1	4337
AvaII	3	1447, 4475, 4697	Eco47III	1	3320
AvrII	1	223	Eco8II	1	881
BamHI	1	3188	EcoICRI	1	16
BanII	3	18, 1292, 5415	EcoNI	1	1885
BbeI	1	304	EcoRI	1	855
BbsI	5	278, 1556, 1672, 2384, 3273	EcoRV	1	1606
BbuI	3	931, 3035, 3107	EheI	1	302
BglI	3	176, 4457, 5725	HaeII	5	304, 3322, 3692, 5331, 5339
BglIII	1	37	HincII	4	1572, 2800, 3196, 5900
BsaI	1	4398	HindII	4	1572, 2800, 3196, 5900
BsaAI	3	422, 5486, 5819	HindIII	1	239
BsaBI	2	42, 2699	HpaI	2	2800, 5900
BsaHI	4	275, 301, 1694, 4874	Hsp92I	4	275, 301, 1694, 4874
BspHI	2	4164, 5172	KasI	1	300
BspMI	1	1666	KpnI	1	12
BsrGI	1	758	MluI	1	22
BssSI	2	3617, 5001			

**Table 12. Restriction Enzymes That Cut the pGL2-Control Vector 1-5 Times (continued).**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
MspAII	5	331, 1621, 3786, 4031, 4972	SallI	1	3194
NaeI	2	3314, 5383	ScaI	1	4817
NarI	1	301	SfilI	1	176
NcoI	2	130, 2940	SgrAI	1	1696
NgoMIV	2	3312, 5381	SinI	3	1447, 4475, 4697
NheI	1	28	SmaI	1	3
NsiI	2	3033, 3105	SphI	3	931, 3035, 3107
NspI	4	931, 3035, 3107, 3448	SpII	1	422
PacI	1	1592	SspI	5	2515, 2668, 5141, 5694, 5809
PaeR7I	1	33	StuI	1	222
PfIMI	1	2236	StyI	4	130, 223, 2384, 2940
Ppu10I	2	3029, 3101	VspI	2	2603, 4509
PpuMI	1	1447	XbaI	1	315
PshAI	1	3259	XcmI	1	1003
Psp5II	1	1447	XhoI	1	33
PspAI	1	1	XmaI	1	1
PvuI	2	4707, 5753	XmnI	1	4936
SacI	1	18			

**Table 13. Restriction Enzymes That Do Not Cut the pGL2-Control Vector.**

AatII	BlpI	CspI	NruI	SgfI
AflIII	Bpu1102I	EagI	PinAI	SnaBI
AgeI	Bsp120I	Eco52I	PmeI	SpeI
ApaI	BssHII	Eco72I	PmlI	SrfI
AscI	Bst1107I	FseI	PstI	Sse8387I
BalI	Bst98I	I-PpoI	PvuII	SwaI
BbrPI	BstXI	NdeI	RsrII	Tth111I
BclI	BstZI	NotI	SacII	

**Note:** The enzymes listed in boldface type are available from Promega.

## 7.F. pGL2-Control Vector Restriction Sites (continued)

**Table 14. Restriction Enzymes That Cut the pGL2-Control Vector More Than Six Times.**

Acil	Bst71I	<b>HaeIII</b>	<b>MboII</b>	Sau96I
<b>AluI</b>	<b>BstOI</b>	HgaI	MnII	ScrFI
<b>BanI</b>	BstUI	<b>HhaI</b>	MseI	SfaNI
BbvI	<b>CfoI</b>	<b>HinfI</b>	<b>MspI</b>	<b>TaqI</b>
BsaOI	<b>DdeI</b>	<b>HpaII</b>	<b>NciI</b>	TfiI
BsaJI	<b>DpnI</b>	HphI	<b>NdeII</b>	<b>Tru9I</b>
<b>BsaMI</b>	DpnII	<b>Hsp92II</b>	NlaIII	<b>XhoII</b>
BsmI	<b>DraI</b>	MaeI	NlaIV	
<b>Bsp1286I</b>	EarI	MaeII	PleI	
BsrI	Fnu4HI	MaeIII	<b>RsaI</b>	
<b>BsrSI</b>	<b>FokI</b>	<b>MboI</b>	<b>Sau3AI</b>	

**Note:** The enzymes listed in boldface type are available from Promega.

## 7.G. References

- Groskreutz, D. *et al.* (1995) Increased expression and convenience with the new pGL3 Luciferase Reporter Vectors. *Promega Notes* **50**, 2-8.
- Sambrook, J. *et al.* (1989) *Molecular Cloning, A Laboratory Manual*, Cold Spring Harbor Press, Cold Spring Harbor, NY.
- Henikoff, S. (1987) Unidirectional digestion with exonuclease III in DNA sequence analysis. *Methods Enzymol.* **155**, 156-65.
- Schenborn, E. *et al.* (1995) Tfx™ Reagent: A new transfection reagent for eukaryotic cells. *Promega Notes* **52**, 2-7.
- Schenborn, E. and Goiffon, V. (1991) Optimization of Transfectam®-mediated transfection using a luciferase reporter system. *Promega Notes* **33**, 8-11.
- Cullen, B. (1987) Use of eukaryotic expression technology in functional analysis of cloned genes. *Methods Enzymol.* **152**, 684-704.
- Ausubel, F.M. *et al.* (1988) *Current Protocols in Molecular Biology*, John Wiley and Sons, NY.
- Rosenthal, N. (1987) Identification of regulatory elements in cloned genes with functional assays. *Methods Enzymol.* **152**, 704-20.
- Hawkins, E., Butler, B. and Wood, K.V. (2000) Bright-Glo™ and Steady-Glo™ Luciferase Assay Systems: Reagents for academic and industrial applications. *Promega Notes* **75**, 3-6.

## 7.H. Related Products

Product	Size	Cat.#
GLprimer2 (counter clockwise)	2µg	E1661
RVprimer3 (clockwise)	2µg	E4481
RVprimer4 (counter clockwise)	2µg	E4491
PureYield™ Plasmid Midiprep System	25 preps	A2492
	100 preps	A2495

## Luciferase Assay Systems

Product	Size	Cat.#
Luciferase Assay System	100 assays	E1500
Bright-Glo™ Luciferase Assay System	10ml	E2610
Steady-Glo® Luciferase Assay System	10ml	E2510
Dual-Luciferase® Reporter Assay System	100 assays	E1910
Dual-Glo® Luciferase Assay System	10ml	E2920
ONE-Glo™ Luciferase Assay System*	10ml	E6110

Available in additional sizes. \*For Laboratory Use.

## pGL4 Luciferase Vectors

Product	Size	Cat.#
pGL4.10[ <i>luc2</i> ] Vector	20µg	E6651
pGL4.11[ <i>luc2P</i> ] Vector	20µg	E6661
pGL4.12[ <i>luc2CP</i> ] Vector	20µg	E6671
pGL4.13[ <i>luc2/SV40</i> ] Vector	20µg	E6681
pGL4.14[ <i>luc2/Hygro</i> ] Vector	20µg	E6691
pGL4.17[ <i>luc2/Neo</i> ] Vector	20µg	E6721
pGL4.20[ <i>luc2/Puro</i> ] Vector	20µg	E6751
pGL4.23[ <i>luc2/minP</i> ] Vector	20µg	E8411
pGL4.26[ <i>luc2/minP/Hygro</i> ] Vector	20µg	E8441
pGL4.29[ <i>luc2P/CRE/Hygro</i> ] Vector	20µg	E8471
pGL4.30[ <i>luc2P/NFAT-RE/Hygro</i> ] Vector	20µg	E8481
pGL4.31[ <i>luc2P/GAL4UAS/Hygro</i> ] Vector	20µg	C9351
pGL4.32[ <i>luc2P/NF-κB-RE/Hygro</i> ] Vector	20µg	E8491

The complete listing of pGL4 Luciferase Vectors can be found at:

[www.promega.com/pgl4/](http://www.promega.com/pgl4/)

## 7.H. Related Products (continued)

### Luminometers

Product	Cat.#
GloMax®-Multi Base Instrument*	E7031
GloMax®-Multi Luminescence Module	E7041
GloMax®-Multi Fluorescence Module	E7051
GloMax®-Multi Absorbance Module	E7061
GloMax® 20/20 Luminometer	E5311
GloMax® 20/20 Luminometer with Single Auto-Injector	E5321
GloMax® 20/20 Luminometer with Dual Auto-Injector	E5331
GloMax® 96 Microplate Luminometer	E6501
GloMax® 96 Microplate Luminometer with Single Reagent Injector	E6511
GloMax® 96 Microplate Luminometer with Dual Reagent Injectors	E6521

\*Cat.# E7031 cannot be sold separately and must be purchased with at least one detection module (Cat.# E7041, E7051 or E7061).

<sup>(4)</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.

© 1992-2009 Promega Corporation. All Rights Reserved.

Dual-Glo, Dual-Luciferase, GloMax, ProFection, Steady-Glo and Transfectam are registered trademarks of Promega Corporation. Bright-Glo, ONE-Glo, PureYield and Tfx are trademarks of Promega Corporation.

DNASTAR is a registered trademark of DNASTAR, Inc. GenBank is a registered trademark of the U.S. Dept. 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 prices and 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.