

TECHNICAL BULLETIN

# pGEM<sup>®</sup>-*luc* Vector

Instructions for Use of Product  
**E1541**



# pGEM<sup>®</sup>-*luc* Vector

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

1.	Description.....	1
2.	Product Components .....	2
3.	pGEM <sup>®</sup> - <i>luc</i> Vector Multiple Cloning Region and Circle Map .....	2
4.	pGEM <sup>®</sup> - <i>luc</i> Vector Restriction Sites.....	4
5.	Related Products.....	6
6.	References.....	6
7.	Summary of Changes .....	6

## 1. Description

The pGEM<sup>®</sup>-*luc* Vector is a cassette vector designed to be a source of the *luc* gene encoding firefly luciferase, which is found in the pGL2 Vectors. The plasmid is not intended for the expression of luciferase in eukaryotic cells.

The pGEM<sup>®</sup>-*luc* Vector was constructed by positioning the luciferase gene (*luc*) (1–3) in the center of the multiple cloning region of the pGEM<sup>®</sup>-11Zf(–) Vector, providing a number of unique restriction sites at both ends of the gene (Figure 1). Sites that are surrounded by parentheses are not unique, as additional sites for each also exists in the luciferase gene. To make use of these nonunique sites, a partial restriction enzyme digest should be performed. Note also that using HindIII or NsiI to clone the luciferase gene will include upstream ATG codons, which may reduce the efficiency of expression in eukaryotes.

Luciferase is a 61kDa monomeric protein that does not require post-translational modifications for enzymatic activity. Thus, it can function as a genetic reporter immediately upon translation (2,3). Luciferase synthesized by in vitro translation can be labeled with <sup>35</sup>S, as the protein contains 4 cysteine and 14 methionine residues. To ensure full enzymatic activity of luciferase, no more than 5 codons can be deleted from either the 5' - or 3' -end of the coding region.

For some experiments, antisense RNA to luciferase mRNA may be useful as a nucleic acid probe. Such antisense RNA can be generated from the T7 RNA polymerase promoter in pGEM<sup>®</sup>-*luc* Vector.

The sequences of Promega vectors are available online at: [www.promega.com/vectors/](http://www.promega.com/vectors/) and are also available from the GenBank<sup>®</sup> database (the GenBank<sup>®</sup>/EMBL Accession Number for the pGEM<sup>®</sup>-*luc* Vector is X65316).

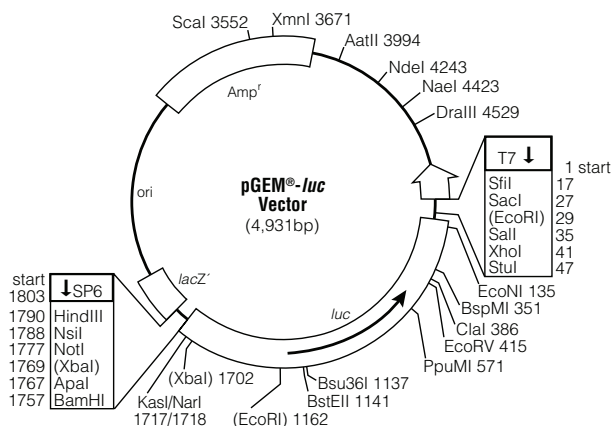
## 2. Product Components

PRODUCT	SIZE	CAT.#
pGEM <sup>®</sup> - <i>luc</i> DNA	20µg	E1541

The pGEM<sup>®</sup>-*luc* Vector is supplied with a glycerol stock of bacterial strain JM109. The JM109 cells do not contain the vector and are not competent cells.

**Storage Conditions:** Store the pGEM<sup>®</sup>-*luc* Vector at  $-30^{\circ}\text{C}$  to  $-10^{\circ}\text{C}$  and the glycerol stock of JM109 cells below  $-65^{\circ}\text{C}$ .

## 3. pGEM<sup>®</sup>-*luc* Vector Multiple Cloning Region and Circle Map

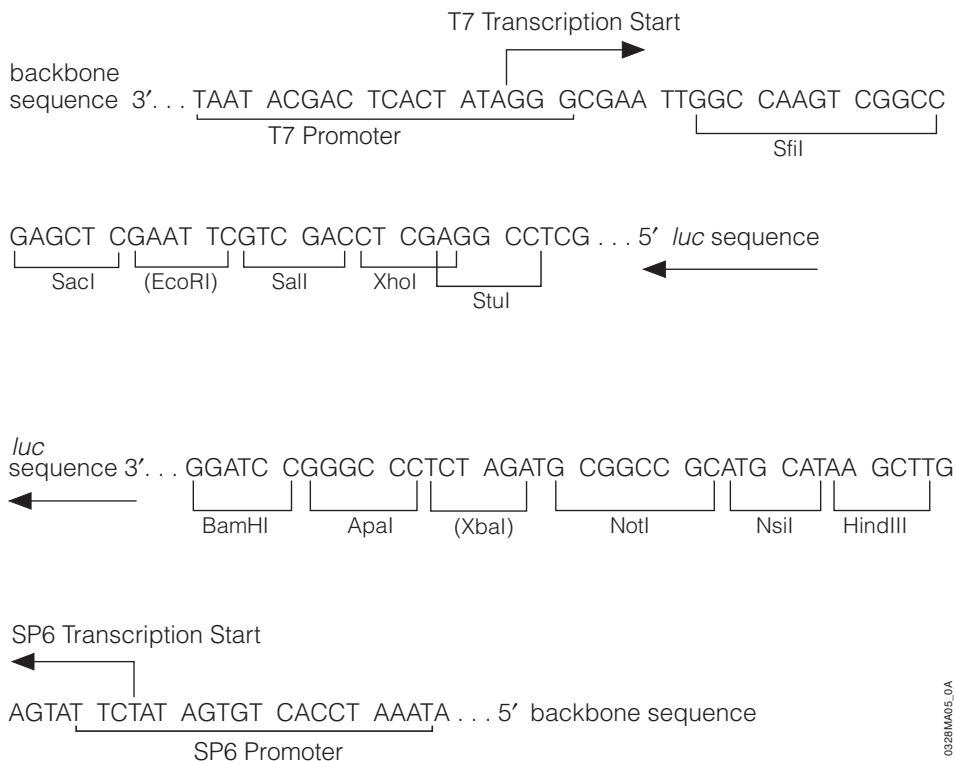


**Figure 1. pGEM<sup>®</sup>-*luc* Vector circle map. Sites shown in parentheses are not unique.**

### pGEM<sup>®</sup>-*luc* Vector sequence reference points:

T7 RNA polymerase transcription initiation site	1
multiple cloning region	10–50; 1757–1795
luciferase cDNA sequences	57–1756
luciferase coding region	105–1754
SP6 RNA polymerase promoter (–17 to +3)	1801–1820
SP6 RNA polymerase transcription initiation site	1803
<i>lac</i> operon sequences	1828–2057; 4754–4912
<i>lacZ</i> start codon	1842
<i>lac</i> operator	1853–1880
β-lactamase coding region	3002–3859
T7 RNA polymerase promoter (–17 to +3)	4915–3

**Note:** *lacZ* start codon is disrupted and therefore inactive.



**Figure 2. pGEM<sup>®</sup>-*luc* Vector promoter and adjacent unique restriction enzyme sites.** The sequence shown corresponds to RNA synthesized by T7 RNA polymerase and is complementary to RNA synthesized by SP6 RNA polymerase. The bold arrow indicates the orientation of the *luc* gene open reading frame.



#### 4. pGEM<sup>®</sup>-*luc* 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. Vector sequences are available in the GenBank<sup>®</sup> database (GenBank<sup>®</sup>/EMBL Accession Number X65316) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Table 1. Restriction Enzymes That Cut the pGEM<sup>®</sup>-*luc* Vector Between 1 and 5 Times.**

<b>Enzyme</b>	<b># of Sites</b>	<b>Location</b>	<b>Enzyme</b>	<b># of Sites</b>	<b>Location</b>
<b>AatII</b>	1	3994	<b>BstEII</b>	1	1141
<b>AccI</b>	1	36	<b>BstZI</b>	2	18, 1777
<b>AccIII</b>	2	538, 1054	<b>Bsu36I</b>	1	1137
AcyI	5	325, 1718, 1744, 3609, 3991	Cfr10I	4	321, 1480, 3152, 4421
AflIII	2	1256, 2179	<b>ClaI</b>	1	386
<b>Alw44 I</b>	3	2493, 3739, 4236	<b>Csp45I</b>	2	794, 1582
AlwNI	1	2595	<b>DraI</b>	3	2938, 2957, 3649
<b>ApaI</b>	1	1767	DraII	3	571, 1764, 4048
<b>AvaI</b>	2	41, 693	DraIII	1	4529
<b>AvaII</b>	3	571, 3210, 3432	DrdI	3	2287, 4156, 4573
<b>BalI</b>	1	11	EagI	2	18, 1777
<b>BamHI</b>	1	1757	<b>EclHKI</b>	1	3072
<b>BanI</b>	4	1717, 1923, 3020, 4485	Eco52I	2	18, 1777
<b>BanII</b>	4	27, 733, 1767, 4455	Eco81I	1	1137
BbeI	1	1721	<b>EcoICRI</b>	1	25
BbsI	3	345, 461, 1739	EcoNI	1	135
<b>BbuI</b>	2	1094, 1786	<b>EcoRI</b>	2	29, 1162
<b>BglI</b>	3	17, 3192, 4764	<b>EcoRV</b>	1	415
BsaI	1	3133	EheI	1	1719
BsaAI	2	1599, 4526	FspI	2	3294, 4771
BsaHI	5	325, 1718, 1744, 3609, 3991	<b>HaeII</b>	5	1721, 2057, 2427, 4371, 4379
Bsp120I	1	1763	<b>HincII</b>	2	37, 449
BspHI	3	2899, 3907, 4012	HindII	2	37, 449
BspMI	1	351	<b>HindIII</b>	1	1790
BsrGI	1	1259	<b>Hsp92I</b>	5	325, 1718, 1744, 3609, 3991
BssSI	3	235, 3736, 4043	KasI	1	1717

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

**Table 1. Restriction Enzymes That Cut the pGEM<sup>®</sup>-*luc* Vector Between 1 and 5 Times. (continued)**

<b>Enzyme</b>	<b># of Sites</b>	<b>Location</b>	<b>Enzyme</b>	<b># of Sites</b>	<b>Location</b>
<b>NaeI</b>	1	4423	<b>SalI</b>	1	35
<b>NarI</b>	1	1718	<b>ScaI</b>	1	3552
<b>NdeI</b>	1	4243	<b>SfiI</b>	1	17
<b>NotI</b>	1	1777	SgrAI	1	321
<b>NsiI</b>	1	1788	<b>SinI</b>	3	571, 3210, 3432
NspI	4	1094, 1786, 2183, 4100	<b>SphI</b>	2	1094, 1786
PacI	1	431	SplI	1	1595
PaeR7I	1	41	<b>SspI</b>	2	3876, 4734
Ppu10I	1	1784	<b>StuI</b>	1	47
PpuMI	1	571	<b>VspI</b>	3	1950, 2009, 3244
Psp5II	1	571	<b>XbaI</b>	2	1702, 1769
<b>PvuI</b>	2	3442, 4792	XcmI	1	1019
<b>PvuII</b>	2	2003, 4821	<b>XhoI</b>	1	41
<b>SacI</b>	1	27	<b>XmnI</b>	1	3671

**Table 2. Restriction Enzymes That Do Not Cut the pGEM<sup>®</sup>-*luc* Vector.**

<b>Acc B7I</b>	Bpu1102I	<b>Eco47III</b>	PflMI	<b>SmaI</b>
<b>Acc65I</b>	BsaBI	Eco72I	PinAI	<b>SnaBI</b>
AflII	<b>BsaMI</b>	FseI	PmeI	<b>SpeI</b>
<b>AgeI</b>	BsmI	<b>HpaI</b>	PmlI	SrfI
AscI	<b>BssHIII</b>	<b>I-PpoI</b>	PshAI	Sse8387I
AvrII	Bst1107I	<b>KpnI</b>	PspAI	<b>StyI</b>
BbrPI	<b>Bst98I</b>	<b>MluI</b>	<b>PstI</b>	Swal
<b>BclI</b>	<b>BstXI</b>	<b>NcoI</b>	RsrII	<b>Tth111I</b>
<b>BglII</b>	<b>CspI</b>	<b>NheI</b>	<b>SacII</b>	<b>XmaI</b>
BlpI	DsaI	<b>NruI</b>	<b>SgfI<sup>(a)</sup></b>	

**Table 3. Restriction Enzymes That Cut the pGEM<sup>®</sup>-*luc* Vector 6 or More Times.**

AcI	Bst71I	<b>FokI</b>	MboI	PleI
<b>AluI</b>	BstOI	<b>HaeIII</b>	<b>MboII</b>	<b>RsaI</b>
Alw26I	BstUI	HgaI	MnlI	<b>Sau3AI</b>
AspH	<b>CfoI</b>	<b>HhaI</b>	MseI	Sau96I
Bbv	DdeI	<b>Hinfl</b>	<b>MspI</b>	ScrFI
BsaOI	<b>DpnI</b>	<b>HpaII</b>	<b>MspA1I</b>	SfaNI
BsaJI	DpnII	HphI	<b>NciI</b>	<b>TaqI</b>
<b>Bsp1286I</b>	EaeI	<b>Hsp92II</b>	<b>NdeII</b>	TfiI
BsrI	EarI	MaeI	NlaIII	<b>Tru9I</b>
<b>BsrSI</b>	Fnu4HI	MaeIII	NlaIV	<b>XhoII</b>

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



## 5. Related Products

Product	Size	Cat.#
Luciferase Assay System	100 Assays	E1500
Luciferase Assay System with Reporter Lysis Buffer	100 assays	E4030
Luciferase Assay System, 10-Pack	1,000 assays	E1501
Luciferase Assay System Freezer Pack	1,000 assays	E4530
Luciferase 1000 Assay System	1,000 assays	E4550
Luciferase Assay Reagent	1,000 assays	E1483
Steady-Glo® Luciferase Assay System	10ml	E2510
	100ml	E2520
	10 × 100ml	E2550
Dual-Glo® Luciferase Assay System	10ml	E2920
	100ml	E2940
	10 × 100ml	E2980
Bright-Glo™ Luciferase Assay System	10ml	E2610
	100ml	E2620
	10 × 100ml	E2650

## 6. References

1. Ow, D. *et al.* (1986) Transient and stable expression of the firefly luciferase gene in plant cells and transgenic plants. *Science* **234**, 856–9.
2. de Wet, J.R. *et al.* (1987) Firefly luciferase gene: structure and expression in mammalian cells. *Mol. Cell Biol.* **7**, 725–37.
3. Wood, K.V. (1990) Firefly Luciferase: A new tool for molecular biologists. *Promega Notes* **28**, 1–3.

## 7. Summary of Changes

The following change was made to the 6/17 revision of this document:

The f1 origin of replication was removed from the reported vector sequence.

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

© 1988–2017 Promega Corporation. All Rights Reserved.

pGEM is a registered trademark of Promega Corporation.

DNASTAR is a registered trademark of DNASTAR, Inc. GenBank is a registered trademark of the U.S. Department of Health and Human Services.

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.