



Promega

Technical Bulletin

pGEM[®]-3Z Vector

INSTRUCTIONS FOR USE OF PRODUCT P2151.



www.promega.com

PRINTED IN USA.
Revised 5/09

Part# TB033

pGEM[®]-3Z Vector

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

1. Description	1
2. Product Components and Storage Conditions	1
3. pGEM [®] -3Z Vector Multiple Cloning Region and Circle Map	2
4. pGEM [®] -3Z Vector Restriction Sites	3
5. Related Products	6
6. Reference	7

1. Description

The pGEM[®]-3Z Vector is intended for use as a standard cloning vector, as well as for the highly efficient synthesis of RNA in vitro. The vector carries the *lacZ* α -peptide and multiple cloning region arrangement from pUC18 (1). In addition, the vector contains both the SP6 and T7 RNA polymerase promoters flanking the multiple cloning region. This arrangement gives rise to a functional α -peptide that is capable of complementing the product of the *lacZ* Δ M15 gene to produce functional β -galactosidase. Cells with the genotype, *lacZ* Δ M15, and also containing the pGEM[®]-3Z Vector will be blue when plated on indicator media containing IPTG and X-Gal. However, when the *lacZ* α -peptide sequence is disrupted by cloning into the pGEM[®]-3Z multiple cloning region, complementation does not occur and no β -galactosidase activity is produced. Therefore, bacterial colonies harboring recombinant pGEM[®]-3Z Vector constructs remain white.

The sequences of Promega vectors are available online at:
www.promega.com/vectors/ and are also available from the GenBank[®]
 database.

2. Product Components and Storage Conditions

Product	Size	Cat.#
pGEM [®] -3Z Vector	20 μ g	P2151

Storage Conditions: Store the pGEM[®]-3Z Vector at -20°C.

3. pGEM[®]-3Z Vector Multiple Cloning Region and Circle Map

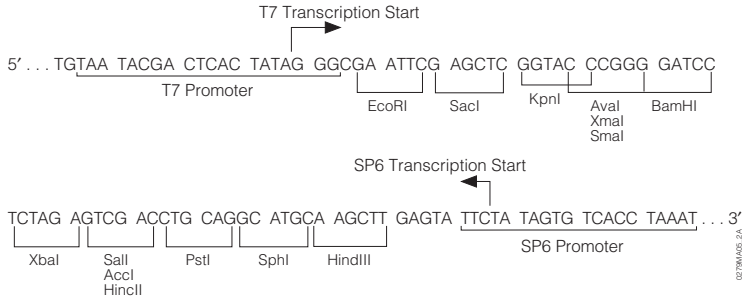


Figure 1. pGEM[®]-3Z Vector promoter and multiple cloning region sequence. The sequence shown corresponds to RNA synthesized by T7 RNA polymerase and is complementary to RNA synthesized by SP6 RNA polymerase.



Note: The pGEM[®]-3Z and pGEM[®]-4Z Vectors are identical except for the orientation of the SP6 and T7 promoters.

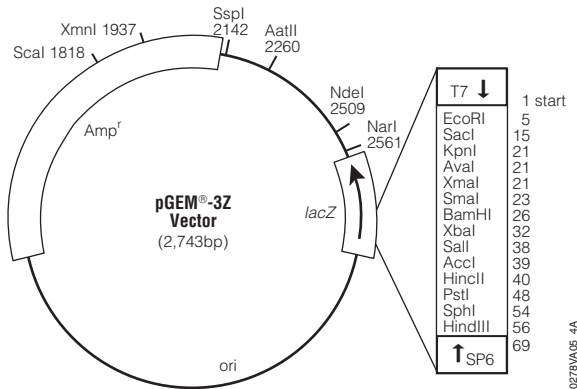


Figure 2. pGEM[®]-3Z Vector circle map and sequence reference points.

pGEM[®]-3Z Vector sequence reference points.

T7 RNA polymerase transcription initiation site	1
multiple cloning region	5-61
SP6 RNA polymerase promoter (-17 to +3)	67-86
SP6 RNA polymerase transcription initiation site	69
<i>lac</i> operon sequences	94-323; 2564-2724
binding site of pUC/M13 Reverse Sequencing Primer	104-125
<i>lacZ</i> start codon	108
<i>lacZ</i> operator	128-144
β -lactamase (<i>Amp^r</i>) coding region	1265-2125
binding site of pUC/M13 Forward Sequencing Primer	2677-2700
T7 RNA polymerase promoter (-17 to +3)	2727-3

Specialized applications of the pGEM[®]-3Z Vector.

- Blue/white screening for recombinants.
- Transcription in vitro from dual, opposed promoters. (For protocol information, please see the *Riboprobe[®] in vitro Transcription Systems Technical Manual*, #TM016.)

4. pGEM[®]-3Z 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 vector sequence is available in the GenBank[®] database (GenBank[®]/EMBL Accession Number X65304) and on the Internet at: www.promega.com/vectors/

Table 1. Restriction Enzymes That Cut the pGEM®-3Z Vector Between 1 and 5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	2260	EheI	1	2562
AccI	1	39	FokI	5	1304, 1485, 1772, 2415, 2659
Acc65I	1	17	FspI	2	1560, 2583
AcyI	3	1875, 2257, 2561	HaeII	3	323, 693, 2564
AflIII	1	445	HgaI	4	556, 1134, 1864, 2422
Alw26I	4	1399, 2175, 2328, 2370	HincII	1	40
Alw44I	3	759, 2005, 2502	HindII	1	40
AlwNI	1	861	HindIII	1	56
AspHI	5	15, 763, 1924, 2009, 2506	Hsp92I	3	1875, 2257, 2561
AvaI	1	21	KasI	1	2560
AvaII	2	1476, 1698	KpnI	1	21
BamHI	1	26	MaeI	4	33, 940, 1193, 1528
BanI	4	17, 189, 1286, 2560	MaeII	5	1148, 1564, 1937, 2257, 2699
BanII	1	15	NarI	1	2561
BbeI	1	2564	NdeI	1	2509
BbuI	1	54	NspI	3	54, 449, 2366
BglI	2	1458, 2576	PleI	5	44, 339, 824, 1327, 2727
BsaI	1	1399	PspAI	1	21
BsaOI	5	361, 785, 1708, 1857, 2604	PstI	1	48
BsaHI	3	1875, 2257, 2561	PvuI	2	1708, 2604
BsaJI	5	21, 22, 184, 605, 2679	PvuII	2	269, 2633
Bsp1286I	5	15, 763, 1924, 2009, 2506	RsaI	3	19, 1818, 2494
BspHI	3	1165, 2173, 2278	SacI	1	15
BspMI	1	51	Sall	1	38
BssSI	3	618, 2002, 2309	ScaI	1	1818
BstOI	5	185, 473, 594, 607, 2680	SinI	2	1476, 1698
Cfr10I	1	1418	SmaI	1	23
DraI	3	1204, 1223, 1915	SphI	1	54
DraII	1	2314	Sse8387I	1	48
DrdI	2	553, 2422	SspI	1	2142
EaeI	3	284, 1726, 2713	TaqI	4	9, 39, 545, 1989
EarI	3	329, 2133, 2621	TfiI	2	280, 420
EclHKI	1	1338	VspI	3	216, 275, 1510
EcoICRI	1	13	XbaI	1	32
EcoRI	1	5	XmaI	1	21
			XmnI	1	1937

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

Table 2. Restriction Enzymes That Do Not Cut the pGEM®-3Z Vector.

AccB7I	BsaBI	DraIII	NgoMIV	RsrII
AccIII	BsaMI	DsaI	NheI	SacII
AflIII	BsmI	EagI	NotI	SfiI
AgeI	Bsp120I	Eco47III	NruI	SgfI
ApaI	BsrGI	Eco52I	NsiI	SgrAI
AscI	BssHIII	Eco72I	Pacl	SnaBI
AvrII	Bst1107I	Eco81I	PaeR7I	SpeI
Ball	Bst98I	EcoNI	PflMI	SpII
BbrPI	BstEII	EcoRV	PinAI	SrfI
BbsI	BstXI	FseI	PmeI	StuI
BclI	BstZI	HpaI	PmlI	StyI
BglII	Bsu36I	I-PpoI	Ppu10I	Swal
BlpI	Clal	MluI	PpuMI	Tth111I
Bpu1102I	CspI	NaeI	PshAI	XcmI
BsaAI	Csp45I	NcoI	Psp5II	XhoI

Table 3. Restriction Enzymes Cut the pGEM®-3Z Vector 6 or More Times.

AccI	CfoI	HinfI	MnII	NlaIV
AluI	DdeI	HpaII	MseI	Sau3AI
BbvI	DpnI	HphI	MspI	Sau96I
BsrI	DpnII	Hsp92II	MspA1I	ScrFI
BsrSI	Fnu4HI	MaellI	NciI	SfaNI
Bst71I	HaeIII	MboI	NdeII	Tru9I
BstUI	HhaI	MboII	NlaIII	XhoII

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

5. Related Products

Product	Size	Cat.#
pGEM [®] -4Z Vector	20µg	P2161
pGEM [®] -3Zf(+) Vector	20µg	P2271
pGEM [®] -3Zf(-) Vector	20µg	P2261
pGEM [®] -5Zf(+) Vector	20µg	P2241
pGEM [®] -5Zf(-) Vector	20µg	P2351
pGEM [®] -7Zf(+) Vector	20µg	P2251
pGEM [®] -7Zf(-) Vector	20µg	P2371
pGEM [®] -9Zf(-) Vector	20µg	P2391
pGEM [®] -11Zf(+) Vector	20µg	P2411
pGEM [®] -11Zf(-) Vector	20µg	P2421
pGEM [®] -13Zf(+) Vector	20µg	P2541

Product	Size	Cat.#
pSP64 Poly(A) Vector	20µg	P1241
pSP72 Vector	20µg	P2191
pSP73 Vector	20µg	P2221

Sequencing Primers

Product	Size	Cat.#
SP6 Promoter Primer	2µg	Q5011
T7 Promoter Primer	2µg	Q5021
pUC/M13 Primer, Reverse (17mer)	2µg	Q5401
pUC/M13 Primer, Forward (17mer)	2µg	Q5391
pUC/M13 Primer, Forward (24mer)	2µg	Q5601
pUC/M13 Primer, Reverse (22mer)	2µg	Q5421

Riboprobe[®] in vitro Transcription Systems

Product	Cat.#
Riboprobe [®] System – SP6	P1420
Riboprobe [®] System – T7	P1440

For Laboratory Use.

RiboMAX[™] Large Scale RNA Production Systems

Product	Cat.#
RiboMAX [™] Large Scale RNA Production System – SP6	P1280
RiboMAX [™] Large Scale RNA Production System – T7	P1300

For Laboratory Use.

6. Reference

1. Yanisch-Perron, C., Vieira, J. and Messing, J. (1985) Improved M13 phage cloning vectors and host strains: Nucleotide sequences of the M13mp18 and pUC19 vectors. *Gene* **33**, 103–19.

© 1988, 1992–1994, 2000, 2004, 2006, 2007, 2009 Promega Corporation. All Rights Reserved.

pGEM and Riboprobe are registered trademarks of Promega Corporation. RiboMAX is a trademark of Promega Corporation. DNASTAR is a registered trademark of DNASTAR, Inc. GenBank is a registered trademark of 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.