

TECHNICAL BULLETIN

pGEM[®]-4Z Vector

Instructions for Use of Product
P2161



pGEM[®]-4Z Vector

All technical literature is available at: 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

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

1. Description

The pGEM[®]-4Z 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*DM15 gene to produce functional β -galactosidase. Cells with the genotype, *lacZ*DM15, and also containing the pGEM[®]-4Z Vector will be blue in color when plated on indicator media containing IPTG and X-Gal. However, when the *lacZ* α -peptide is disrupted by cloning into the pGEM[®]-4Z multiple cloning region, complementation does not occur, and no β -galactosidase activity is produced. Therefore, bacterial colonies harboring recombinant pGEM[®]-4Z Vector constructs remain white.

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



2. Product Components and Storage Conditions

PRODUCT	SIZE	CAT.#
pGEM [®] -4Z Vector	20µg	P2161

The pGEM[®]-4Z Vector is provided with a glycerol stock of bacterial strain JM109. The JM109 cells do not contain the vector and are not competent.

Storage Conditions: Store the pGEM[®]-4Z Vector at -20°C and the glycerol stock of JM109 cells at -70°C.

3. pGEM[®]-4Z Vector Multiple Cloning Region and Circle Map

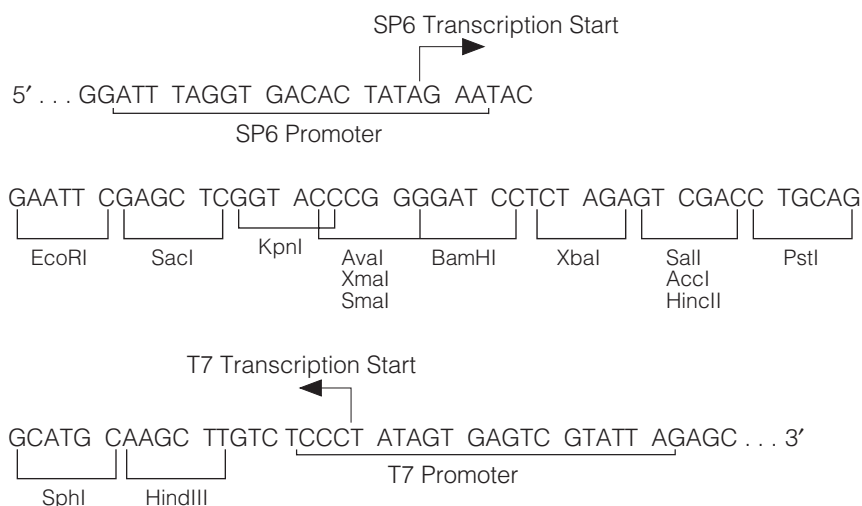
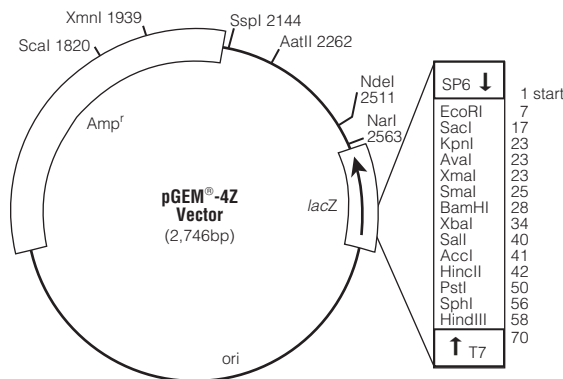


Figure 1. pGEM[®]-4Z Vector promoter and multiple cloning region sequence. The sequence shown corresponds to RNA synthesized by SP6 RNA polymerase and is complementary to RNA synthesized by T7 RNA polymerase.



0280VAGE_4A

Figure 2. pGEM®-4Z Vector circle map and sequence reference points. The pGEM®-3Z and pGEM®-4Z Vectors are identical except for the orientation of the SP6 and T7 promoters.

pGEM®-4Z Vector sequence reference points:

SP6 RNA polymerase transcription initiation site	1
multiple cloning region	7–63
T7 RNA polymerase promoter (–17 to +3)	68–87
T7 RNA polymerase transcription initiation site	70
<i>lac</i> operon sequences	96–325; 2566–2726
<i>lacZ</i> start codon	110
<i>lacZ</i> operator	130–146
β-lactamase (Amp ^r) coding region	1267–2127
SP6 RNA polymerase promoter (–17 to +3)	2730–3

Specialized applications of the pGEM®-4Z 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[®]-4Z 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. Vector sequences are available in the GenBank[®] database (GenBank[®]/EMBL Accession Number X65305) and on the Internet at: www.promega.com/vectors/

Table 1. Restriction Enzymes That Cut the pGEM[®]-4Z Vector Between 1 and 5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	2262	BstXI	1	2725
AccI	1	41	Cfr10I	1	1420
Acc65I	1	19	DraI	3	1206, 1225, 1917
AcyI	3	1877, 2259, 2563	DraII	1	2316
AflIII	1	447	DrdI	2	555, 2424
Alw26I	5	69, 1401, 2177, 2330, 2372	EaeI	3	286, 1728, 2715
Alw44I	3	761, 2007, 2504	EarI	3	331, 2135, 2623
AlwNI	1	863	EclHKI	1	1340
AspHI	5	17, 765, 1926, 2011, 2508	EcoICRI	1	15
AvaI	1	23	EcoRI	1	7
AvaII	2	1478, 1700	EheI	1	2564
BamHI	1	28	FokI	5	1306, 1487, 1774, 2417, 2661
BanI	4	19, 191, 1288, 2562	FspI	2	1562, 2585
BanII	1	17	HaeII	3	325, 695, 2566
BbeI	1	2566	HgaI	4	558, 1136, 1866, 2424
BbuI	1	56	HincII	1	42
BglI	2	1460, 2578	HindII	1	42
BsaI	1	1401	HindIII	1	58
BsaHI	3	1877, 2259, 2563	Hsp92I	3	1877, 2259, 2563
BsaJI	5	23, 24, 186, 607, 2681	KasI	1	2562
BsaOI	5	363, 787, 1710, 1859, 2606	KpnI	1	23
Bsp1286I	5	17, 765, 1926, 2011, 2508	MaeI	4	35, 942, 1195, 1530
BspHI	3	1167, 2175, 2280	MaeII	5	1150, 1566, 1939, 2259, 2701
BspMI	1	53	NarI	1	2563
BssSI	3	620, 2004, 2311			
BstOI	5	187, 475, 596, 609, 2682			

Table 1. Restriction Enzymes That Cut the pGEM[®]-4Z Vector Between 1 and 5 Times (continued).

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
NdeI	1	2511	SinI	2	1478, 1700
NspI	3	56, 451, 2368	SmaI	1	25
PleI	5	46, 85, 341, 826, 1329	SphI	1	56
PspAI	1	23	Sse8387I	1	50
PstI	1	50	SspI	1	2144
PvuI	2	1710, 2606	TaqI	4	11, 41, 547, 1991
PvuII	2	271, 2635	TfiI	2	282, 422
RsaI	3	21, 1820, 2496	VspI	3	218, 277, 1512
SacI	1	17	XbaI	1	34
SalI	1	40	XmaI	1	23
ScaI	1	1820	XmnI	1	1939

Table 2. Restriction Enzymes That Do Not Cut the pGEM[®]-4Z Vector.

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



4. pGEM[®]-4Z Vector Restriction Sites (continued)

Table 3. Restriction Enzymes That Cut the pGEM[®]-4Z Vector 6 or More Times.

AclI	CfoI	HinfI	MnlI	NlaIV
AluI	DdeI	HpaII	MseI	Sau3AI
BbvI	DpnI	HphI	MspI	Sau96I
BsrI	DpnII	Hsp92II	MspAII	ScrFI
BsrSI	Fnu4HI	MaeIII	NciI	SfaNI
Bst71I	HaeIII	MboI	NdeII	Tru9I
BstUI	HhaI	MboII	NlaIII	XhoII

5. Related Products

pGEM[®] Vectors

Product	Size	Cat.#
pGEM [®] -3Z Vector	20µg	P2151
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 [®] -13Zf(+) Vector	20µg	P2541

All pGEM[®] Vectors are provided with a glycerol stock of bacterial strain JM109. The JM109 cells do not contain the vector and are not competent.

Other Vectors

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

Riboprobe® in vitro Transcription Systems

Product	Cat.#
Riboprobe® System—SP6	P1420
Riboprobe® System—T7	P1440

RiboMAX™ Large-Scale RNA Production Systems

Product	Cat.#
RiboMAX™ Large Scale RNA Production System—SP6	P1280
RiboMAX™ Large Scale RNA Production System—T7	P1300
T7 RiboMAX™ Express Large Scale RNA Production System	P1320

© 1988, 1992–1995, 2000, 2004, 2007, 2009, 2017 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. Department 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.