



Promega

Technical Bulletin

pGEM[®]-3Zf(+) Vector

INSTRUCTIONS FOR USE OF PRODUCT P2271.



www.promega.com

PRINTED IN USA.
Revised 11/07



AF9TB086 1107TB086

Part# TB086

pGEM[®]-3Zf(+) 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

I. Description.....	1
II. Product Components and Storage Conditions	2
III. pGEM [®] -3Zf(+) Vector Multiple Cloning Region and Circle Map.....	2
IV. pGEM [®] -3Zf(+) Vector Restriction Sites	4
V. Related Products	6
VI. Reference	7

I. Description

The pGEM[®]-3Zf(+) Vector is a derivative of the pGEM[®]-3Z Vector and contains the origin of replication of the filamentous phage f1. The plasmid serves as a standard cloning vector, as a template for in vitro transcription and as a template for the production of circular ssDNA.

The pGEM[®]-3Zf(+) Vector contains SP6 and T7 RNA polymerase promoters flanking a multiple cloning region within the α -peptide coding region of β -galactosidase (1). Insertional inactivation of the α -peptide allows recombinant clones to be directly identified by color screening on indicator plates. The multiple cloning region is unique and includes restriction sites for EcoRI, SacI, KpnI, Aval, SmaI, BamHI, XbaI, Sall, Accl, HincII, PstI, SphI and HindIII.

For induction of ssDNA, bacterial cells with the F' episome (e.g., JM109, XL-1 Blue, DH5 α [™]) containing pGEM[®]-3Zf(+) Vector recombinants are infected with an appropriate helper phage. The plasmid then enters the f1 replication mode, and the resulting ssDNA is exported from the cell as an encapsidated virus-like particle. The sequence of the ssDNA rescued upon infection with helper phage is complementary to the sequence shown in Figure 1. The exported ssDNA can be used for mutagenesis in vitro or can be sequenced using the T7 Promoter Primer or pUC/M13 Forward Primer.

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

II. Product Components and Storage Conditions

Product	Size	Cat.#
pGEM [®] -3Zf(+) Vector	20µg	P2271

Storage Conditions: Store the pGEM[®]-3Zf(+) Vector at -20°C.

III. pGEM[®]-3Zf(+) Vector Multiple Cloning Region and Circle Map

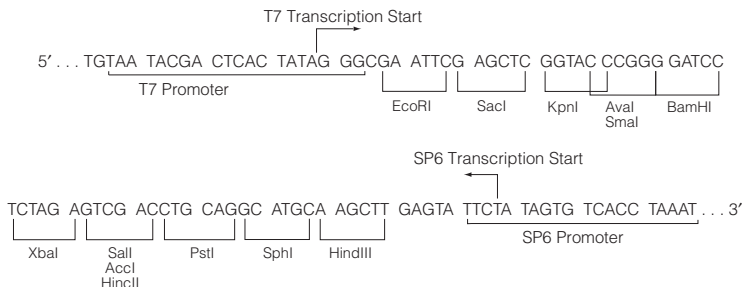


Figure 1. pGEM[®]-3Zf(+) Vector promoter and multiple cloning region sequence. The sequence shown corresponds to RNA synthesized by T7 RNA polymerase and is complementary to RNA synthesized by SP6 RNA polymerase. The strand shown is complementary to the ssDNA strand produced by this vector.

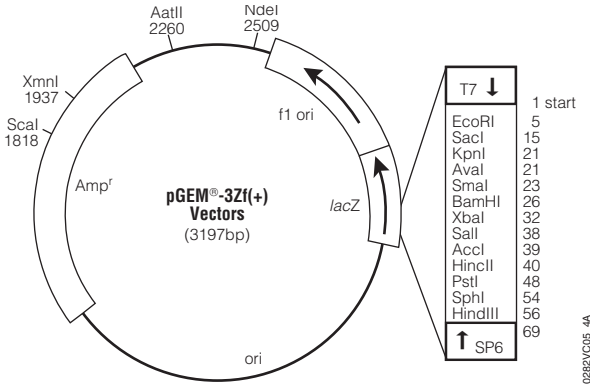


Figure 2. pGEM®-3Zf(+) Vector circle map and sequence reference points. The pGEM®-3Zf(+) and pGEM®-3Zf(-) Vectors are identical except for the orientation of the f1 origin. Use the T7 Promoter Primer or pUC/M13 Forward Primer to sequence ssDNA produced by the pGEM®-3Zf(+) Vector.

pGEM®-3Zf(+) Vector sequence reference points:

T7 RNA polymerase transcription initiation site	1
SP6 RNA polymerase transcription initiation site	69
T7 RNA polymerase promoter (-17 to +3)	3181-3
SP6 RNA polymerase promoter (-17 to +3)	67-86
multiple cloning region	5-61
<i>lacZ</i> start codon	108
<i>lac</i> operon sequences	3018-3178; 94-323
<i>lac</i> operator	128-144
β -lactamase (Amp ^r) coding region	1265-2125
phage f1 region	2562-3017
binding site of pUC/M13 Forward Sequencing Primer	3138-3154
binding site of pUC/M13 Reverse Sequencing Primer	104-120

Specialized applications of the pGEM®-3Zf(+) Vector:

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

Note: All Promega technical literature is available on the Internet at:
www.promega.com

IV. pGEM[®]-3Zf(+) 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 X65306) and on the Internet at: www.promega.com/vectors/

Table 1. Restriction Enzymes That Cut the pGEM[®]-3Zf(+) Vector Between 1 and 5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	2260	Cfr10I	2	1418, 2887
AccI	1	39	DraI	3	1204, 1223, 1915
Acc65I	1	17	DraII	1	2314
AcyI	2	1875, 2257	DraIII	1	2786
AflIII	1	445	DrdI	3	553, 2422, 2741
Alw26I	4	1399, 2175, 2328, 2370	EaeI	3	284, 1726, 3167
Alw44I	3	759, 2005, 2502	EarI	3	329, 2133, 3075
AlwNI	1	861	EclHKI	1	1338
AspHI	5	15, 763, 1924, 2009, 2506	EcoICRI	1	13
AvaI	1	21	EcoRI	1	5
AvaII	2	1476, 1698	FokI	5	1304, 1485, 1772, 2415, 3113
BamHI	1	26	FspI	2	1560, 3037
BanI	4	17, 189, 1286, 2823	HaeII	4	323, 693, 2937, 2945
BanII	2	15, 2861	HgaI	5	556, 1134, 1864, 2422, 3003
BbuI	1	54	HincII	1	40
BglI	2	1458, 3030	HindII	1	40
BsaI	1	1399	HindIII	1	56
BsaOI	5	361, 785, 1708, 1857, 3058	Hsp92I	2	1875, 2257
BsaAI	1	2786	KpnI	1	21
BsaHI	2	1875, 2257	MaeI	5	33, 940, 1193, 1528, 2937
BsaJI	5	21, 22, 184, 605, 3133	NaeI	1	2889
BspHI	3	1165, 2173, 2278	NdeI	1	2509
BspMI	1	51	NgoMIV	1	2887
BssSI	3	618, 2002, 2309	NspI	3	54, 449, 2366
BstOI	5	185, 473, 594, 607, 3134	PspAI	1	21
			PstI	1	48
			PvuI	2	1708, 3058

Table 1. Restriction Enzymes That Cut the pGEM[®]-3Zf(+) Vector Between 1 and 5 Times (continued).

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
PvuII	2	269, 3087	SspI	2	2142, 2578
RsaI	3	19, 1818, 2494	TaqI	5	9, 39, 545, 1989, 2819
SacI	1	15	TfiI	2	280, 420
Sall	1	38	VspI	3	216, 275, 1510
ScaI	1	1818	XbaI	1	32
SinI	2	1476, 1698	XmaI	1	21
SmaI	1	23	XmnI	1	1937
SphI	1	54			
Sse8387I	1	48			

Table 2. Restriction Enzymes That Do Not Cut the pGEM[®]-3Zf(+) Vector.

AccB7I	BsaMI	EagI	NotI	SgfI
AccIII	BsmI	Eco47III	NruI	SgrAI
AflIII	Bsp120I	Eco52I	NsiI	SnaBI
AgeI	BsrBRI	Eco72I	PaeR7I	SpeI
Apal	BsrGI	Eco81I	PacI	SpII
AscI	BssHII	EcoNI	PfiMI	SrfI
AvrII	Bst1107I	EcoRV	PinAI	StuI
BalI	Bst98I	EheI	PmeI	StyI
BbeI	BstEII	FseI	PmlI	Swal
BbrPI	BstXI	HpaI	Ppu10I	Tth111I
BbsI	BstZI	I-PpoI	PpuMI	XcmI
BclI	Bsu36I	KasI	PshAI	XhoI
BglII	Clal	MluI	Psp5II	
BlpI	CspI	NarI	RsrII	
Bpu1102I	Csp45I	NcoI	SacII	
BsaBI	DsaI	NheI	SfiI	

Table 3. Restriction Enzymes That Cut the pGEM[®]-3Zf(+) Vector 6 or More Times.

AcI	DdeI	Hsp92II	NciI	Tru9I
AluI	DpnI	MaeII	NdeII	XhoII
BbvI	DpnII	MaeIII	NlaIII	
Bsp1286I	Fnu4HI	MboI	NlaIV	
BsrI	HaeIII	MboII	PleI	
BsrSI	HhaI	MnlI	Sau3AI	
Bst7II	HinfI	MseI	Sau96I	
BstUI	HpaII	MspI	ScrFI	
CfoI	HphI	MspAII	SfaNI	

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

V. Related Products

Product	Size	Cat.#
pGEM [®] -3Z Vector	20µg	P2151
pGEM [®] -4Z Vector	20µg	P2161
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 – T3	P1430
Riboprobe [®] System – T7	P1440

For Laboratory Use.

VI. Reference

1. Yanish-Perron, C. *et al.* (1985) Improved M13 phage cloning vectors and host strains: Nucleotide sequences of the M13mp18 and pUC19 vectors. *Gene* **33**, 103-19.

© 1990-2007 Promega Corporation. All Rights Reserved.

pGEM and Riboprobe are registered trademarks of Promega Corporation.

DH5 α is a trademark of Life Technologies, Inc. DNASTAR is a registered trademark of DNASTAR, Inc. GenBank is a registered trademark of the 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.