



Promega

Technical Bulletin

pSP73 Vector

INSTRUCTIONS FOR USE OF PRODUCT P2221.



www.promega.com

PRINTED IN USA.
Revised 9/06



AF9TE041 0906TE041

Part# TB041

pSP73 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	1
III. pSP73 Vector Multiple Cloning Region and Circle Map.....	2
IV. pSP73 Vector Restriction Sites.....	4
V. Related Products	6
VI. Reference	7

I. Description

The pSP73 Vector (1) offers a wide range of restriction sites, providing greater versatility in cloning and transcription of RNA *in vitro*. The pSP73 Vector contains the SP6 and T7 RNA polymerase promoters and a unique multiple cloning region, which includes restriction sites for BglII, EcoRV, ClaI, EcoRI, SacI, KpnI, SmaI, BamHI, XbaI, AccI, Sall, PstI, SphI, HindIII, PvuII and XhoI.

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.#
pSP73 Vector	20µg	P2221

Storage Conditions: Store the pSP73 Vector at -20°C.

III. pSP73 Vector Multiple Cloning Region and Circle Map

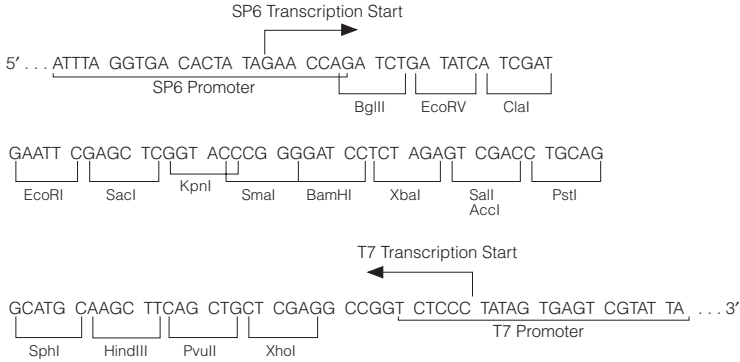


Figure 1. pSP73 Vector promoter and multiple cloning region sequence. The sequence shown corresponds to RNA synthesized by SP6 RNA polymerase and is complementary to RNA synthesized by T7 RNA polymerase.

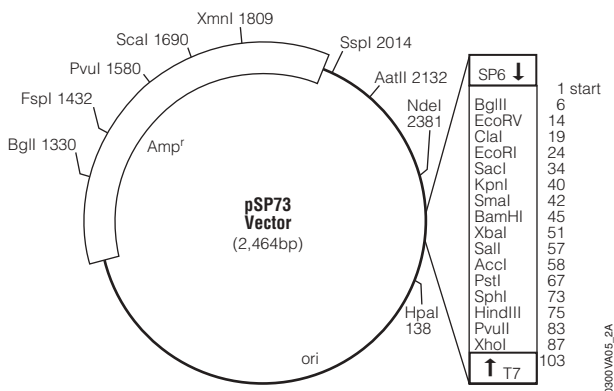


Figure 2. pSP73 Vector circle map and sequence reference points.

pSP73 Vector sequence reference points.

SP6 RNA polymerase transcription initiation site	1
T7 RNA polymerase transcription initiation site	103
SP6 RNA polymerase promoter (-17 to +3)	2448-3
T7 RNA polymerase promoter (-17 to +3)	107-120
multiple cloning region	6-92
β -lactamase coding region	1137-1997

! The pSP72 and pSP73 Vectors are identical except for the orientation of the multiple cloning region.

Note: Blue/white screening of recombinants is **not** possible with the pSP73 Vector.

Specialized application of the pSP73 Vector.

- Transcription in vitro from dual-opposed promoters (For protocol information, please request the Riboprobe® in vitro Transcription Systems Technical Manual, #TM016) available at: www.promega.com/tbs/

IV. pSP73 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 also available in the GenBank® database (GenBank®/EMBL Accession Number X65333) and on the Internet at: www.promega.com/vectors/

Table 1. Restriction Enzymes That Cut the pSP73 Vector Between 1 and 5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	2132	ClaI	1	19
AccI	1	58	DraI	3	1076, 1095, 1787
Acc65I	1	36	DraII	1	2186
AcyI	2	1747, 2129	DrdI	2	425, 2294
AflIII	1	317	EaeI	2	156, 1598
Alw26I	5	102, 1271, 2047, 2200, 2242	EarI	2	201, 2005
Alw44I	3	631, 1877, 2374	EclHKI	2	1210, 2391
AlwNI	1	733	EcoICRI	1	32
AspHI	5	34, 635, 1796, 1881, 2378	EcoRI	1	24
AvaI	2	40, 87	EcoRV	1	14
AvaII	2	1348, 1570	FokI	4	1176, 1357, 1644, 2287
BamHI	1	45	FspI	1	1432
BanI	2	36, 1158	HaeII	2	195, 565
BanII	1	34	HgaI	4	428, 1006, 1736, 2294
BbuI	1	73	HincII	2	59, 138
BglI	1	1330	HindII	2	59, 138
BglIII	1	6	HindIII	1	75
BsaI	2	102, 1271	HpaI	1	138
BsaOI	4	233, 657, 1580, 1729	Hsp92I	2	1747, 2129
BsaHI	2	1747, 2129	KpnI	1	40
BsaJI	3	40, 41, 477	MaeI	4	52, 812, 1065, 1400
Bsp1286I	5	34, 635, 1796, 1881, 2378	MaeII	4	1020, 1436, 1809, 2129
BspHI	3	1037, 2045, 2150	MspAII	5	83, 659, 904, 1845, 2311
BspMI	2	70, 149	NdeI	1	2381
BssSI	3	490, 1874, 2181	NspI	3	73, 321, 2238
BstOI	4	133, 345, 466, 479	Paer7I	1	87
Cfr10I	2	93, 1290			

Table 1. Restriction Enzymes That Cut the pSP73 Vector Between 1 and 5 Times (continued).

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
PleI	5	63, 118, 211, 696, 1199	SinI	2	1348, 1570
PspAI	1	40	SmaI	1	42
PstI	1	67	SphI	1	73
PvuI	1	1580	Sse8387I	1	67
PvuII	1	83	SspI	1	2014
RsaI	3	38, 1690, 2366	TfiI	2	152, 292
SacI	1	34	VspI	4	118, 147, 1382, 2415
SalI	1	57	XbaI	1	51
Sau96I	5	1252, 1331, 1348, 1570, 2186	XhoI	1	87
ScaI	1	1690	XmaI	1	40
			XmnI	1	1809

Table 2. Restriction Enzymes That Do Not Cut the pSP73 Vector.

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

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

IV. pSP73 Vector Restriction Sites (continued)

Table 3. Restriction Enzymes That Cut the pSP73 Vector 6 or More Times.

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

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

V. Related Products

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

Product	Size	Cat.#
pGEM®-3Z Vector	20µg	P2151
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

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

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

For Laboratory Use.

VI. Reference

1. Krieg, P.A. and Melton, D.A. (1987) In vitro RNA synthesis with SP6 RNA polymerase. *Methods Enzymol.* **155**, 397-415.

© 1988-2006 Promega Corporation. All Rights Reserved.

pGEM and RiboProbe are registered trademarks of Promega Corporation.

DNASar is a registered trademark of DNASar, 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.