



**Promega**

# Technical Bulletin

---

## **pSP72 Vector**

INSTRUCTIONS FOR USE OF PRODUCT P2191.



[www.promega.com](http://www.promega.com)

PRINTED IN USA.  
Revised 9/06



AF9TB040 0906TB040

Part# TB040

# pSP72 Vector

All technical literature is available on the Internet at: [www.promega.com/tbs/](http://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](mailto:techserv@promega.com)

I. Description.....	1
II. Product Components and Storage Conditions .....	1
III. pSP72 Vector Multiple Cloning Region and Circle Map .....	2
IV. pSP72 Vector Restriction Sites .....	4
V. Related Products .....	6
VI. Reference .....	6

## I. Description

The pSP72 Vector (1) can be used as a standard cloning vector and can also be used for transcription of RNA in vitro. The pSP72 Vector contains the SP6 and T7 RNA polymerase promoters flanking a unique multiple cloning region, which includes restriction sites for XhoI, PvuII, HindIII, SphI, PstI, SalI, AccI, XbaI, BamHI, SmaI, KpnI, SacI, EcoRI, ClaI, EcoRV and BglIII.

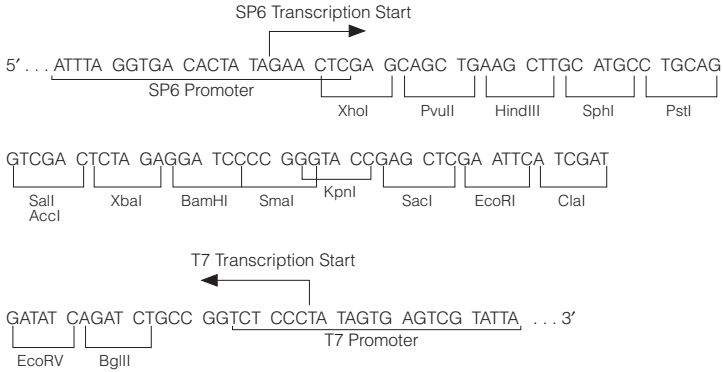
The sequences of Promega vectors are available online at: [www.promega.com/vectors/](http://www.promega.com/vectors/) and are also available from the GenBank® database.

## II. Product Components and Storage Conditions

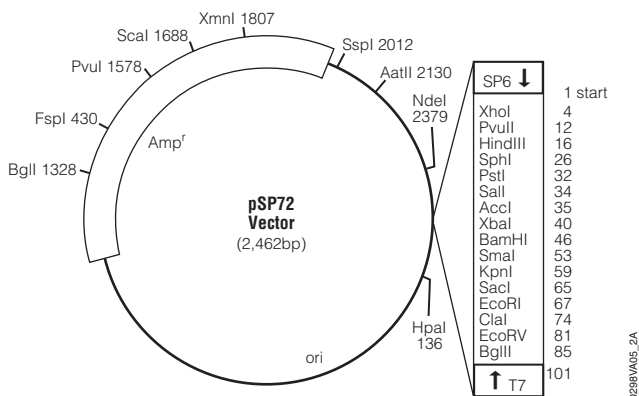
Product	Size	Cat.#
pSP72 Vector	20µg	P2191

**Storage Conditions:** Store the pSP72 Vector at -20°C.

### III. pSP72 Vector Multiple Cloning Region and Circle Map



**Figure 1. pSP72 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.



0298/VA05\_2A

**Figure 2. pSP72 Vector circle map and sequence reference points.**

**pSP72 Vector sequence reference points.**

SP6 RNA polymerase transcription initiation site	1
T7 RNA polymerase transcription initiation site	101
SP6 RNA polymerase promoter (-17 to +3)	2446-3
T7 RNA polymerase promoter (-17 to +3)	99-118
multiple cloning region	4-90
β-lactamase coding region	1135-1995

**!** 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 pSP72 Vector.

**Specialized application of the pSP72 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/](http://www.promega.com/tbs/)

#### IV. pSP72 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 X65332) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

**Table 1. Restriction Enzymes That Cut the pSP72 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	2130	<b>DraI</b>	3	1074, 1093, 1785
<b>AccI</b>	1	35	<b>DraII</b>	1	2184
<b>Acc65I</b>	1	55	<b>DrdI</b>	2	423, 2292
<b>AcyI</b>	2	1745, 2127	<b>EaeI</b>	2	154, 1596
<b>AflIII</b>	1	315	<b>EarI</b>	2	199, 2003
<b>Alw26I</b>	5	100, 1269, 2045, 2198, 2240	<b>EclHKI</b>	2	1208, 2389
<b>Alw44I</b>	3	629, 1875, 2372	<b>EcoICRI</b>	1	63
<b>AlwNI</b>	1	731	<b>EcoRI</b>	1	67
<b>AspHI</b>	5	65, 633, 1794, 1879, 2376	<b>EcoRV</b>	1	81
<b>AvaI</b>	2	4, 51	<b>FokI</b>	4	1174, 1355, 1642, 2285
<b>AvaII</b>	2	1346, 1568	<b>FspI</b>	1	1430
<b>BamHI</b>	1	46	<b>HaeII</b>	2	193, 563
<b>BanI</b>	2	55, 1156	<b>HgaI</b>	4	426, 1004, 1734, 2292
<b>BanII</b>	1	65	<b>HincII</b>	2	36, 136
<b>BbuI</b>	1	26	<b>HindII</b>	2	36, 136
<b>BglI</b>	1	1328	<b>HindIII</b>	1	16
<b>BglIII</b>	1	85	<b>HpaI</b>	1	136
<b>BsaI</b>	2	100, 1269	<b>Hsp92I</b>	2	1745, 2127
<b>BsaOI</b>	4	231, 655, 1578, 1727	<b>KpnI</b>	1	59
<b>BsaHI</b>	2	1745, 2127	<b>MaeI</b>	4	41, 810, 1063, 1398
<b>BsaJI</b>	3	50, 51, 475	<b>MaeII</b>	4	1018, 1434, 1807, 2127
<b>Bsp1286I</b>	5	65, 633, 1794, 1879, 2376	<b>MspAII</b>	5	12, 657, 902, 1843, 2309
<b>BspHI</b>	3	1035, 2043, 2148	<b>NdeI</b>	1	2379
<b>BspMI</b>	2	21, 147	<b>NspI</b>	3	26, 319, 2236
<b>BssSI</b>	3	488, 1872, 2179	<b>PaeR7I</b>	1	4
<b>BstOI</b>	4	131, 343, 464, 477	<b>PleI</b>	5	31, 116, 209, 694, 1197
<b>Cfr10I</b>	2	91, 1288			
<b>ClaI</b>	1	74			

**Table 1. Restriction Enzymes That Cut the pSP72 Vector Between 1 and 5 Times (continued).**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
PspAI	1	51	SmaI	1	53
PstI	1	32	SphI	1	26
PvuI	1	1578	Sse8387I	1	32
PvuII	1	12	SspI	1	2012
RsaI	3	57, 1688, 2364	TfiI	2	150, 290
SacI	1	65	VspI	4	116, 145, 1380, 2413
SalI	1	34	XbaI	1	40
Sau96I	5	1250, 1329, 1346, 1568, 2184	XhoI	1	4
ScaI	1	1688	XmaI	1	51
SinI	2	1346, 1568	XmnI	1	1807

**Table 2. Restriction Enzymes That Do Not Cut the pSP72 Vector.**

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

**Table 3. Restriction Enzymes That Cut the pSP72 Vector 6 or More Times.**

AcI	<b>CfoI</b>	<b>HinFI</b>	MnII	<b>Sau3AI</b>
AluI	<b>DdeI</b>	<b>HpaII</b>	MseI	ScrFI
BbvI	<b>DpnI</b>	HphI	<b>MspI</b>	SfaNI
BsrI	DpnII	<b>Hsp92II</b>	<b>NciI</b>	<b>TaqI</b>
<b>BsrSI</b>	Fnu4HI	MaeIII	<b>NdeII</b>	<b>Tru9I</b>
Bst7II	<b>HaeIII</b>	MboI	NlaIII	<b>XhoII</b>
BstUI	<b>HhaI</b>	<b>MboII</b>	NlaIV	

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

## V. Related Products

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

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.

DNAS<sup>®</sup> is a registered trademark of DNAS<sup>®</sup>, 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.