

## pSP-*luc*+NF Fusion Vector Restriction Enzyme Tables

The following restriction enzyme tables were constructed using DNASTAR® sequence analysis software. This information has not been verified 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). Vector sequences are also available in the GenBank® database (GenBank®/EMBL Accession Number U47123) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pSP-*luc*+NF Fusion Vector 1-5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
<b>AatII</b>	1	3771	EarI	5	132, 307, 850, 1840, 3644
<b>AccIII</b>	2	743, 1259	<b>EclHKI</b>	2	2849, 4030
<b>Acc65I</b>	1	11	EcoNI	3	605, 1005, 1665
AcyI	5	55, 81, 1474, 3386, 3768	<b>EcoRI</b>	1	1708
AflIII	2	541, 1956	<b>EcoRV</b>	1	1716
<b>Alw44I</b>	3	2270, 3516, 4013	EheI	1	82
AlwNI	1	2372	FspI	1	3071
AspHI	5	1513, 2274, 3435, 3520, 4017	<b>HaeII</b>	4	84, 997, 1834, 2204
<b>AvaI</b>	2	1104, 1726	<b>HincII</b>	2	1352, 1777
<b>AvaII</b>	3	1227, 2987, 3209	HindII	2	1352, 1777
AvrII	1	30	<b>HindIII</b>	1	36
<b>BanI</b>	4	11, 80, 908, 2797	<b>HpaI</b>	1	1777
<b>BanII</b>	1	1072	<b>Hsp92I</b>	5	55, 81, 1474, 3386, 3768
BbeI	1	84	KasI	1	80
BbsI	2	1336, 1452	<b>KpnI</b>	1	15
<b>BbuI</b>	1	711	<b>NarI</b>	1	81
<b>BclI</b>	1	628	<b>NcoI</b>	1	43
<b>BglI</b>	1	2969	<b>NdeI</b>	1	4020
<b>BglII</b>	1	24	<b>NheI</b>	1	17
BsaI	2	1741, 2910	NspI	3	711, 1960, 3877
BsaOI	5	1132, 1872, 2296, 3219, 3368	PaeR7I	1	1726
BsaHI	5	55, 81, 1474, 3386, 3768	PpuMI	1	1227
BspHI	4	631, 2676, 3684, 3789	Psp5II	1	1227
BspMI	3	1437, 1446, 1788	<b>PvuI</b>	1	3219
BsrGI	1	538	<b>ScaI</b>	2	213, 3329
BssSI	3	2129, 3513, 3820	SgrAI	1	1476
<b>BstEII</b>	1	47	<b>SinI</b>	3	1227, 2987, 3209
Cfr10I	4	317, 1476, 1732, 2929	<b>SphI</b>	1	711
<b>Clal</b>	1	1721	<b>SspI</b>	1	3653
<b>Csp45I</b>	1	217	<b>StyI</b>	2	30, 43
<b>DraI</b>	3	2715, 2734, 3426	<b>VspI</b>	4	1757, 1786, 3021, 4054
DraII	2	1227, 3825	<b>XbaI</b>	1	1702
DrdI	3	1449, 2064, 3933	XcmI	1	783
DsaI	2	43, 418	<b>XhoI</b>	1	1726
EaeI	2	1795, 3237	<b>XmnI</b>	1	3448

### Restriction Enzymes That Do Not Cut the pSP-*luc*+NF Fusion Vector .

<b>AccB7I</b>	BbrPI	<b>BssHII</b>	EagI	<b>MluI</b>	PinAI	RsrII	<b>SpeI</b>
<b>AccI</b>	BlpI	Bst1107I	<b>Eco47III</b>	<b>NaeI</b>	PmeI	<b>SacI</b>	SpII
AflII	Bpu1102I	<b>Bst98I</b>	Eco52I	<b>NgoMIV</b>	PmlI	<b>SacII</b>	SrfI
<b>AgeI</b>	BsaAI	<b>BstXI</b>	Eco72I	<b>NotI</b>	Ppu10I	<b>SalI</b>	Sse8387I
<b>ApaI</b>	BsaBI	<b>BstZI</b>	Eco81I	<b>NruI</b>	PshAI	<b>SfiI</b>	<b>StuI</b>
AscI	<b>BsaMI</b>	<b>Bsu36I</b>	<b>EcoICRI</b>	<b>NsiI</b>	PspAI	<b>SgfI</b>	Swal
<b>BalI</b>	BsmI	<b>CspI</b>	FseI	Pacl	<b>PstI</b>	<b>SmaI</b>	<b>Tth111I</b>
<b>BamHI</b>	Bsp120I	DraIII	<b>I-PpoI</b>	PflMI	<b>PvuII</b>	<b>SnaBI</b>	<b>XmaI</b>

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