

## pCBG99-Control 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 AY258596) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pCBG99-Control Vector 1-5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
<b>AatII</b>	1	506	Eco52I	3	1922, 1926, 5064
<b>AccB7I</b>	1	330	<b>EcoICRI</b>	1	9
<b>AccI</b>	2	986, 2424	FseI	1	1928
<b>Acc65I</b>	1	1	FspI	2	3788, 4961
AcyI	2	503, 4103	<b>HaeII</b>	5	1313, 2551, 2921, 4560, 4568
AflIII	4	15, 945, 1782, 2673	HgaI	5	910, 2784, 3362, 4092, 4493
<b>Alw26I</b>	3	1051, 3627, 4403	<b>HincII</b>	4	591, 1827, 2069, 2425
<b>Alw44I</b>	2	2987, 4233	HindIII	4	591, 1827, 2069, 2425
AlwNI	1	3089	<b>HindIII</b>	1	245
<b>ApaI</b>	1	1420	<b>HpaI</b>	1	2069
AspHI	5	11, 369, 2991, 4152, 4237	<b>Hsp92I</b>	2	503, 4103
<b>AvaI</b>	3	26, 32, 1327	<b>KpnI</b>	1	5
<b>AvaII</b>	2	3704, 3926	<b>MluI</b>	2	15, 945
AvrII	1	229	<b>NciI</b>	5	27, 28, 3053, 3749, 4100
<b>BamHI</b>	1	2417	<b>NcoI</b>	1	278
<b>BanII</b>	5	11, 33, 1165, 1420, 4644	<b>NheI</b>	1	21
BbsI	3	645, 1181, 2502	<b>NotI</b>	1	5064
<b>BbuI</b>	2	2275, 2347	<b>NruI</b>	2	1397, 1664
<b>BclI</b>	2	1074, 1592	<b>NsiI</b>	3	1660, 2273, 2345
<b>BglI</b>	3	182, 3686, 4954	NspI	3	2275, 2347, 2677
<b>BglIII</b>	1	36	PaeR7I	1	32
BsaI	2	1051, 3627	PflMI	1	330
BsaAI	1	4715	Ppu10I	3	1656, 2269, 2341
BsaBI	2	48, 2170	PshAI	1	2488
BsaHI	2	503, 4103	PspAI	1	26
Bsp120I	1	1416	<b>PvuI</b>	2	3936, 4982
BspHI	2	3393, 4401	<b>PvuII</b>	2	960, 1624
BspMI	1	5194	<b>SacI</b>	1	11
BsrBRI	2	48, 2170	<b>Sall</b>	1	2423
BssSI	3	1168, 2846, 4230	<b>Scal</b>	3	1801, 4046, 5129
<b>BstXI</b>	1	484	<b>SfiI</b>	1	182
<b>BstZI</b>	3	1922, 1926, 5064	<b>SinI</b>	2	3704, 3926
<b>ClaI</b>	3	2164, 5122, 5226	<b>SmaI</b>	1	28
<b>DraI</b>	5	813, 2130, 3432, 3451, 4143	<b>SphI</b>	2	2275, 2347
DraII	1	1450	SrfI	1	28
DraIII	2	1198, 4718	<b>SspI</b>	5	669, 747, 4370, 4923, 5038
DrdI	2	2781, 4762	<b>StuI</b>	1	228
DsaI	1	278	<b>StyI</b>	4	229, 278, 398, 1493
EaeI	5	1247, 1922, 1926, 3954, 5064	<b>VspI</b>	1	3738
EagI	3	1922, 1926, 5064	<b>XbaI</b>	1	1909
EarI	3	2557, 4361, 4999	<b>XhoI</b>	1	32
<b>EclHKI</b>	1	3566	<b>XmaI</b>	1	26
<b>Eco47III</b>	2	1311, 2549	<b>XmnI</b>	1	4165

### Restriction Enzymes That Do Not Cut the pCBG99-Control Vector.

<b>AccIII</b>	BbrPI	<b>Bst98I</b>	Eco81I	KasI	PmlI	<b>SgfI</b>	Swal
AflIII	BlpI	<b>BstEII</b>	EcoNI	<b>NarI</b>	PpuMI	SgrAI	<b>Tth111I</b>
<b>AgeI</b>	Bpu1102I	<b>Bsu36I</b>	<b>EcoRI</b>	<b>NdeI</b>	Psp5II	<b>SnaBI</b>	XcmI
AscI	BsrGI	<b>CspI</b>	<b>EcoRV</b>	Pacl	<b>PstI</b>	<b>SpeI</b>	
<b>BaI</b>	<b>BssHII</b>	<b>Csp45I</b>	EheI	PinAI	RsrII	SpII	
BbeI	Bst1107I	Eco72I	<b>I-PpoI</b>	PmeI	<b>SacII</b>	Sse8387I	

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