

## pCBG68-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 AY258594) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pCBG68-Control Vector 1-5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
<b>AatII</b>	1	410	<b>EclHKI</b>	1	3583
<b>AccI</b>	2	1763, 2441	<b>Eco47III</b>	1	2566
<b>Acc65I</b>	1	1	Eco52I	3	1939, 1943, 5081
AcyI	3	407, 1814, 4120	<b>EcoICRI</b>	1	9
AflIII	2	15, 2690	FseI	1	1945
<b>AgeI</b>	1	1849	FspI	2	3805, 4978
<b>Alw44I</b>	2	3004, 4250	<b>HincII</b>	3	1173, 2086, 2442
AlwNI	1	3106	HindII	3	1173, 2086, 2442
<b>ApaI</b>	2	1420, 1453	<b>HindIII</b>	1	245
<b>AvaI</b>	2	26, 32	<b>HpaI</b>	1	2086
<b>AvaII</b>	2	3721, 3943	<b>Hsp92I</b>	3	407, 1814, 4120
AvrII	1	229	<b>KpnI</b>	1	5
<b>BalI</b>	1	1228	<b>MluI</b>	1	15
<b>BamHI</b>	1	2434	<b>NaeI</b>	3	1943, 2560, 4629
<b>BanII</b>	5	11, 33, 1420, 1453, 4661	<b>NcoI</b>	1	278
BbsI	3	341, 1906, 2519	<b>NgoMIV</b>	3	1941, 2558, 4627
<b>BbuI</b>	2	2292, 2364	<b>NheI</b>	1	21
<b>BclI</b>	1	1124	<b>NotI</b>	1	5081
<b>BglI</b>	3	182, 3703, 4971	<b>NsiI</b>	2	2290, 2362
<b>BglII</b>	1	36	NspI	3	2292, 2364, 2694
BlpI	1	1253	PaeR7I	1	32
Bpu1102I	1	1253	PinAI	1	1849
BsaI	3	1393, 1464, 3644	Ppu10I	2	2286, 2358
BsaAI	2	1473, 4732	PshAI	1	2505
BsaBI	2	48, 2187	PspAI	1	26
BsaHI	3	407, 1814, 4120	<b>PvuI</b>	3	1578, 3953, 4999
<b>BsaMI</b>	3	252, 2007, 2100	<b>SacI</b>	1	11
BsmI	3	252, 2007, 2100	<b>SalI</b>	1	2440
Bsp120I	2	1416, 1449	<b>Scal</b>	2	4063, 5146
BspHI	3	1917, 3410, 4418	<b>SfiI</b>	1	182
BspMI	1	5211	<b>SinI</b>	2	3721, 3943
BsrBRI	2	48, 2187	<b>SmaI</b>	1	28
BssSI	4	400, 944, 2863, 4247	<b>SphI</b>	2	2292, 2364
<b>BstXI</b>	2	484, 1236	SrfI	1	28
<b>BstZI</b>	3	1939, 1943, 5081	<b>SspI</b>	3	4387, 4940, 5055
<b>ClaI</b>	3	2181, 5139, 5243	<b>StuI</b>	1	228
<b>Csp45I</b>	1	446	<b>StyI</b>	2	229, 278
<b>DraI</b>	4	2147, 3449, 3468, 4160	TfiI	4	595, 1828, 2665, 5136
DraII	2	1449, 1450	<b>VspI</b>	1	3755
DraIII	1	4735	<b>XbaI</b>	1	1926
DrdI	3	1672, 2798, 4779	<b>XhoI</b>	1	32
DsaI	2	278, 750	<b>XmaI</b>	1	26
EagI	3	1939, 1943, 5081	<b>XmnI</b>	1	4182
EarI	4	364, 2574, 4378, 5016			

### Restriction Enzymes That Do Not Cut the pCBG68-Control Vector.

<b>AccB7I</b>	BsrGI	<b>CspI</b>	EheI	Pacl	<b>PstI</b>	<b>SnaBI</b>	XcmI
<b>AccIII</b>	<b>BssHII</b>	Eco72I	<b>I-PpoI</b>	PflMI	<b>PvuII</b>	<b>SpeI</b>	
AflII	Bst1107I	Eco81I	KasI	PmeI	RsrII	SplI	
AscI	<b>Bst98I</b>	EcoNI	<b>NarI</b>	PmlI	<b>SacII</b>	Sse8387I	
BbeI	<b>BstEII</b>	EcoRI	<b>NdeI</b>	PpuMI	<b>SgfI</b>	Swal	
BbrPI	<b>Bsu36I</b>	EcoRV	<b>NruI</b>	Psp5II	SgrAI	<b>Tth111I</b>	

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