

pCBG68-Basic 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 AY258593) and on the Internet at: www.promega.com/vectors/

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

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	218	EarI	4	172, 2136, 3940, 4578
AccI	2	1571, 2003	EclHKI	1	3145
Acc65I	1	1	Eco47III	1	2128
AcyI	3	215, 1622, 3682	Eco52I	3	1747, 1751, 4643
AflIII	2	15, 2252	EcoICRI	1	9
AgeI	1	1657	FseI	1	1753
Alw44I	2	2566, 3812	FspI	2	3367, 4540
AlwNI	1	2668	HincII	3	981, 1894, 2004
ApaI	2	1228, 1261	HindII	3	981, 1894, 2004
AvaI	2	26, 32	HindIII	1	53
AvaII	2	3283, 3505	HpaI	1	1894
BalI	1	1036	Hsp92I	3	215, 1622, 3682
BamHI	1	1996	KpnI	1	5
BanII	5	11, 33, 1228, 1261, 4223	MluI	1	15
BbsI	3	149, 1714, 2081	MspAII	5	154, 1088, 2594, 2839, 3780
BclI	1	932	NaeI	3	1751, 2122, 4191
BglI	2	3265, 4533	NcoI	1	86
BglII	1	36	NgoMIV	3	1749, 2120, 4189
BlpI	1	1061	NheI	1	21
Bpu1102I	1	1061	NotI	1	4643
BsaI	3	1201, 1272, 3206	NspI	1	2256
BsaAI	2	1281, 4294	PaeR7I	1	32
BsaBI	1	1995	PinAI	1	1657
BsaHI	3	215, 1622, 3682	PshAI	1	2067
BsaMI	3	60, 1815, 1908	PspAI	1	26
BsmI	3	60, 1815, 1908	PvuI	3	1386, 3515, 4561
Bsp120I	2	1224, 1257	SacI	1	11
BspHI	3	1725, 2972, 3980	SalI	1	2002
BspMI	1	4773	ScaI	2	3625, 4708
BsrBRI	1	1995	SinI	2	3283, 3505
BssSI	4	208, 752, 2425, 3809	SmaI	1	28
BstXI	2	292, 1044	SrfI	1	28
BstZI	3	1747, 1751, 4643	SspI	3	3949, 4502, 4617
ClaI	3	1989, 4701, 4805	StyI	1	86
Csp45I	1	254	TfiI	4	403, 1636, 2227, 4698
DraI	4	1955, 3011, 3030, 3722	VspI	1	3317
DraII	2	1257, 1258	XbaI	1	1734
DraIII	1	4297	XhoI	1	32
DrdI	3	1480, 2360, 4341	XmaI	1	26
DsaI	2	86, 558	XmnI	1	3744
EagI	3	1747, 1751, 4643			

Restriction Enzymes That Do Not Cut the pCBG68-Basic Vector.

AccB7I	BbuI	CspI	I-PpoI	PflMI	PvuII	SpeI	XcmI
AccIII	BsrGI	Eco72I	KasI	PmeI	RsrII	SphI	
AflII	BssHII	Eco81I	NarI	PmlI	SacII	SpII	
AscI	Bst1107I	EcoNI	NdeI	Ppu10I	SfiI	Sse8387I	
AvrII	Bst98I	EcoRI	NruI	PpuMI	SgfI	StuI	
BbeI	BstEII	EcoRV	NsiI	Psp5II	SgrAI	Swal	
BbrPI	Bsu36I	EheI	Pacl	PstI	SnaBI	Tth111I	

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