

pCBR-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 AY258592) and on the Internet at: www.promega.com/vectors/

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

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	506	FspI	2	3791, 4964
AccB7I	1	330	HaeII	4	2554, 2924, 4563, 4571
AccI	2	986, 2427	HgaI	5	910, 2787, 3365, 4095, 4496
Acc65I	1	1	HincII	4	591, 1827, 2072, 2428
AcyI	2	503, 4106	HindII	4	591, 1827, 2072, 2428
AflIII	3	15, 1782, 2676	HindIII	1	245
Alw26I	3	1051, 3630, 4406	HpaI	1	2072
Alw44I	2	2990, 4236	Hsp92I	2	503, 4106
AlwNI	1	3092	KpnI	1	5
ApaI	1	1420	MluI	1	15
AspHI	5	11, 369, 2994, 4155, 4240	NciI	5	27, 28, 3056, 3752, 4103
AvaI	3	26, 32, 1327	NcoI	1	278
AvaII	2	3707, 3929	NheI	1	21
AvrII	1	229	NotI	1	5067
BamHI	1	2420	NruI	2	1397, 1664
BanII	5	11, 33, 1165 1420, 4647	NsiI	3	1660, 2276, 2348
BbsI	3	645, 1181, 2505	NspI	3	2278, 2350, 2680
BbuI	2	2278, 2350	PaeR7I	1	32
BclI	2	1074, 1592	PfiMI	1	330
BglI	3	182, 3689, 4957	Ppu10I	3	1656, 2272, 2344
BglIII	1	36	PshAI	1	2491
BsaI	2	1051, 3630	PspAI	1	26
BsaAI	1	4718	PvuI	2	3939, 4985
BsaBI	2	48, 2173	PvuII	2	960, 1624
BsaHI	2	503, 4106	SacI	1	11
BstXI	1	484	SalI	1	2426
BstZI	3	1925, 1929, 5067	ScaI	3	1801, 4049, 5132
ClaI	3	2167, 5125, 5229	SfiI	1	182
DraI	5	813, 2133, 3435, 3454, 4146	SinI	2	3707, 3929
DraII	1	1450	SmaI	1	28
DraIII	2	1198, 4721	SphI	2	2278, 2350
DrdI	2	2784, 4765	SrfI	1	28
DsaI	1	278	SspI	5	669, 747, 4373, 4926, 5041
EaeI	5	1247, 1925, 1929, 3957, 5067	StuI	1	228
EagI	3	1925, 1929, 5067	StyI	4	229, 278, 398, 1493
EarI	3	2560, 4364, 5002	VspI	1	3741
EclHKI	1	3569	XbaI	1	1912
Eco47III	1	2552	XhoI	1	32
Eco52I	3	1925, 1929, 5067	XmaI	1	26
EcoICRI	1	9	XmnI	1	4168
FseI	1	1931			

Restriction Enzymes That Do Not Cut the pCBR-Control Vector.

AccIII	BbrPI	Bst1107I	Eco72I	I-PpoI	PmeI	SacII	Sse8647I
AflII	BlpI	Bst98I	Eco81I	KasI	PmlI	SgfI	Swal
AgeI	Bpu1102I	BstEII	EcoNI	NarI	PpuMI	SgrAI	Tth111I
AscI	Bpu1268I	Bsu36I	EcoRI	NdeI	Psp5II	SnaBI	XcmI
BalI	BsrGI	CspI	EcoRV	Pacl	PstI	SpeI	
BbeI	BssHII	Csp45I	EheI	PinAI	RsrII	SplI	

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