

pCBR-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 AY258591) and on the Internet at: www.promega.com/vectors/

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

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	314	Eco52I	3	1733, 1737, 4629
AccB7I	1	138	EcoICRI	1	9
AccI	2	794, 1989	FseI	1	1739
Acc65I	1	1	FspI	2	3353, 4526
AcyI	2	311, 3668	HaeII	4	2116, 2486, 4125, 4133
AflIII	3	5, 1590, 2238	HgaI	5	718, 2349, 2927, 3657, 4058
Alw26I	3	859, 3192, 3968	HincII	4	399, 1635, 1880, 1990
Alw44I	2	2552, 3798	HindII	4	399, 1635, 1880, 1990
AlwNI	1	2654	HindIII	1	53
ApaI	1	1228	HpaI	1	1880
AspHI	5	11, 177, 2556, 3717, 3802	Hsp92I	2	311, 3668
AvaI	3	26, 32, 1135	KpnI	1	5
AvaII	2	3269, 3491	MluI	1	15
BamHI	1	1982	NciI	5	27, 28, 2618, 3314, 3665
BanII	5	11, 33, 973, 1228, 4209	NcoI	1	86
BbsI	3	453, 989, 2067	NheI	1	21
BclI	2	882, 1400	NotI	1	4629
BglI	2	3251, 4519	NruI	2	1205, 1472
BglII	1	36	NsiI	1	1468
BsaI	2	859, 3192	NspI	1	2242
BsaAI	1	4280	PaeR7I	1	32
BsaBI	1	1981	PfiMI	1	138
BsaHI	2	311, 3668	Ppu10I	1	1464
Bsp120I	1	1224	PshAI	1	2053
BspHI	2	2958, 3966	PspAI	1	26
BspMI	1	4759	PvuI	2	3501, 4547
BsrBRI	1	1981	PvuII	2	768, 1432
BssSI	3	976, 2411, 3795	SacI	1	11
BstXI	1	292	SalI	1	1988
BstZI	3	1733, 1737, 4629	ScaI	3	1609, 3611, 4694
ClaI	3	1975, 4687, 4791	SinI	2	3269, 3491
DraI	5	621, 1941, 2997, 3016, 3708	SmaI	1	28
DraII	1	1258	SrfI	1	28
DraIII	2	1006, 4283	SspI	5	477, 555, 3935, 4488, 4603
DrdI	2	2346, 4327	StyI	3	86, 206, 1301
DsaI	1	86	VspI	1	3303
EaeI	5	1055, 1733, 1737, 3519, 4629	XbaI	1	1720
EagI	3	1733, 1737, 4629	XhoI	1	32
EarI	3	2122, 3926, 4564	XmaI	1	26
EclHKI	1	3131	XmnI	1	3730
Eco47III	1	2114			

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

AccIII	BbrPI	Bst98I	EcoNI	NdeI	PstI	SpeI	XcmI
AflIII	BbuI	BstEII	EcoRI	Pacl	RsrII	SphI	
AgeI	BlpI	Bsu36I	EcoRV	PinAI	SacII	SpII	
AscI	Bpu1102I	CspI	EheI	Pmel	SfiI	Sse8387I	
AvrII	BsrGI	Csp45I	I-PpoI	PmlI	SgfI	StuI	
BalI	BssHIII	Eco72I	KasI	PpuMI	SgrAI	Swal	
BbeI	Bst1107I	Eco81I	NarI	Psp5II	SnaBI	Tth111I	

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