

pGL2-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 X65324) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL2-Control Vector 1-5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccB7I	1	2236	EheI	1	302
AccI	1	3195	HaeII	5	304, 3322, 3692, 5331, 5339
AccIII	2	963, 1479	HincII	4	1572, 2800, 3196, 5900
Acc65I	1	8	HindII	4	1572, 2800, 3196, 5900
AcyI	4	275, 301, 1694, 4874	HindIII	1	239
AflIII	3	22, 761, 3444	HpaI	2	2800, 5900
Alw26I	4	1291, 1523, 4398, 5174	Hsp92I	4	275, 301, 1694, 4874
Alw44I	2	3758, 5004	KasI	1	300
AlwNI	1	3860	KpnI	1	12
AspHI	5	18, 1733, 3762, 4923, 5008	MluI	1	22
AvaI	3	1, 33, 1324	MspAII	5	331, 1621, 3786, 4031, 4972
AvaII	3	1447, 4475, 4697	NaeI	2	3314, 5383
AvrII	1	223	NarI	1	301
BamHI	1	3188	NcoI	2	130, 2940
BanII	3	18, 1292, 5415	NgoMIV	2	3312, 5381
BbeI	1	304	NheI	1	28
BbsI	5	278, 1556, 1672, 2384, 3273	NsiI	2	3033, 3105
BbuI	3	931, 3035, 3107	NspI	4	931, 3035, 3107, 3448
BglI	3	176, 4457, 5725	Pacl	1	1592
BglII	1	37	PaeR7I	1	33
BsaI	1	4398	PflMI	1	2236
BsaAI	3	422, 5486, 5819	Ppu10I	2	3029, 3101
BsaBI	2	42, 2699	PpuMI	1	1447
BsaHI	4	275, 301, 1694, 4874	PshAI	1	3259
BspHI	2	4164, 5172	Psp5II	1	1447
BspMI	1	1666	PspAI	1	1
BsrGI	1	758	PvuI	2	4707, 5753
BssSI	2	3617, 5001	SacI	1	18
BstEII	1	875	SalI	1	3194
Bsu36I	1	881	ScaI	1	4817
Cfr10I	5	537, 1696, 3312, 4417, 5381	SfiI	1	176
ClaI	1	1633	SgrAI	1	1696
Csp45I	2	437, 1225	SinI	3	1447, 4475, 4697
DraII	1	1447	SmaI	1	3
DraIII	1	5489	SphI	3	931, 3035, 3107
DrdI	2	3552, 5533	SpII	1	422
DsaI	2	130, 2940	SspI	5	2515, 2668, 5141, 5694, 5809
EaeI	1	4725	StuI	1	222
EclHKI	1	4337	StyI	4	130, 223, 2384, 2940
Eco47III	1	3320	VspI	2	2603, 4509
Eco81I	1	881	XbaI	1	315
EcoICRI	1	16	XcmI	1	1003
EcoNI	1	1885	XhoI	1	33
EcoRI	1	855	XmaI	1	1
EcoRV	1	1606	XmnI	1	4936

Restriction Enzymes That Do Not Cut the pGL2-Control Vector.

AatII	BclI	BstXI	I-PpoI	PstI	SrfI
AflIII	BlpI	BstZI	NdeI	PvuII	Sse8387I
AgeI	Bpu1102I	CspI	NotI	RsrII	Swal
ApaI	Bsp120I	EagI	NruI	SacII	Tth111I
AscI	BssHII	Eco52I	PinAI	SgfI	
BalI	Bst1107I	Eco72I	PmeI	SnaBI	
BbrPI	Bst98I	FseI	PmlI	SpeI	

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