

pGL2-Promoter 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 X65326) and on the Internet at: www.promega.com/vectors/

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

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccB7I	1	2236	EheI	1	302
AccI	1	2938	FspI	2	4302, 5475
AccIII	2	963, 1479	HaeII	5	304, 3065, 3435, 5074, 5082
Acc65I	1	8	HincII	4	1572, 2800, 2939, 5643
AcyI	4	275, 301, 1694, 4617	HindII	4	1572, 2800, 2939, 5643
AflIII	3	22, 761, 3187	HindIII	1	239
Alw26I	4	1291, 1523, 4141, 4917	HpaI	2	2800, 5643
Alw44I	2	3501, 4747	Hsp92I	4	275, 301, 1694, 4617
AlwNI	1	3603	KasI	1	300
AspHI	5	18, 1733, 3505, 4666, 4751	KpnI	1	12
AvaI	3	1, 33, 1324	MluI	1	22
AvaII	3	1447, 4218, 4440	MspAII	5	331, 1621, 3529, 3774, 4715
AvrII	1	223	NaeI	2	3057, 5126
BamHI	1	2931	NarI	1	301
BanII	3	18, 1292, 5158	NcoI	1	130
BbeI	1	304	NgoMIV	2	3055, 5124
BbsI	5	278, 1556, 1672, 2384, 3016	NheI	1	28
BbuI	1	931	NspI	2	931, 3191
BglI	3	176, 4200, 5468	Pacl	1	1592
BglIII	1	37	PaeR7I	1	33
BsaI	1	4141	PflMI	1	2236
BsaAI	3	422, 5229, 5562	PpuMI	1	1447
BsaBI	2	42, 2699	PshAI	1	3002
BsaHI	4	275, 301, 1694, 4617	Psp5II	1	1447
BspHI	2	3907, 4915	PspAI	1	1
BspMI	1	1666	PvuI	2	4450, 5496
BsrGI	1	758	SacI	1	18
BssSI	2	3360, 4744	SalI	1	2937
BstEII	1	875	Scal	1	4560
Bsu36I	1	881	SfiI	1	176
Cfr10I	5	537, 1696, 3055, 4160, 5124	SgrAI	1	1696
ClaI	1	1633	SinI	3	1447, 4218, 4440
Csp45I	2	437, 1225	SmaI	1	3
DraII	1	1447	SphI	1	931
DraIII	1	5232	SpII	1	422
DrdI	2	3295, 5276	SspI	5	2515, 2668, 4884, 5437, 5552
DsaI	1	130	StuI	1	222
EaeI	1	4468	StyI	3	130, 223, 2384
EclHKI	1	4080	VspI	2	2603, 4252
Eco47III	1	3063	XbaI	1	315
Eco81I	1	881	XcmI	1	1003
EcoICRI	1	16	XhoI	1	33
EcoNI	1	1885	XmaI	1	1
EcoRI	1	855	XmnI	1	4679
EcoRV	1	1606			

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

AatII	BclI	BstXI	I-PpoI	PmlI	SnaBI
AflIII	BlpI	BstZI	NdeI	Ppu10I	SpeI
AgeI	Bpu1102I	CspI	NotI	PstI	SrfI
ApaI	Bsp120I	EagI	NruI	PvuII	Sse8387I
AscI	BssHII	Eco52I	NsiI	RsrII	Swal
BalI	Bst1107I	Eco72I	PinAI	SacII	Tth111I
BbrPI	Bst98I	FseI	PmeI	SgfI	

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