

pGL2-Enhancer 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 X65325) and on the Internet at: www.promega.com/vectors/

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

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccB7I	1	2044	EheI	1	110
AccI	1	3003	FspI	2	4367, 5540
AccIII	2	771, 1287	HaeII	5	112, 3130, 3500, 5139, 5147
Acc65I	1	8	HincII	4	1380, 2608, 3004, 5708
AcyI	4	83,109,1502, 4682	HindIII	4	1380, 2608, 3004, 5708
AflIII	3	22, 569, 3252	HindIII	1	47
Alw26I	4	1099, 1331, 4206, 4982	HpaI	2	2608, 5708
Alw44I	2	3566, 4812	Hsp92I	4	83, 109, 1502, 4682
AlwNI	1	3668	KasI	1	108
AspHI	5	18, 1541, 3570, 4731, 4816	KpnI	1	12
AvaI	3	1, 33, 1132	MluI	1	22
AvaII	3	1255, 4283, 4505	MspAII	5	139, 1429, 3594, 3839, 4780
BamHI	1	2996	NaeI	2	3122, 5191
BanII	3	18, 1100, 5223	NarI	1	109
BbeI	1	112	NcoI	1	2748
BbsI	5	86, 1364, 1480, 2192, 3081	NgoMIV	2	3120, 5189
BbuI	3	739, 2843, 2915	NheI	1	28
BglI	2	4265, 5533	NsiI	2	2841, 2913
BglII	1	37	NspI	4	739, 2843, 2915, 3256
BsaI	1	4206	Pacl	1	1400
BsaAI	3	230, 5294, 5627	PaeR7I	1	33
BsaBI	1	2507	PflMI	1	2044
BsaHI	4	83, 109,1502, 4682	Ppu10I	2	2837, 2909
BspHI	2	3972, 4980	PpuMI	1	1255
BspMI	1	1474	PshAI	1	3067
BsrGI	1	566	Psp5II	1	1255
BssSI	2	3425, 4809	PspAI	1	1
BstEII	1	683	PvuI	2	4515, 5561
Bsu36I	1	689	SacI	1	18
Cfr10I	5	345, 1504, 3120, 4225, 5189	SalI	1	3002
ClaI	1	1441	Scal	1	4625
Csp45I	2	245, 1033	SgrAI	1	1504
DraII	1	1255	SinI	3	1255, 4283, 4505
DraIII	1	5297	SmaI	1	3
DrdI	2	3360, 5341	SphI	3	739, 2843, 2915
DsaI	1	2748	SpII	1	230
EaeI	1	4533	SspI	5	2323, 2476, 4949 5502, 5617
EclHKI	1	4145	StyI	2	2192, 2748
Eco47III	1	3128	VspI	2	2411, 4317
Eco81I	1	689	XbaI	1	123
EcoICRI	1	16	XcmI	1	811
EcoNI	1	1693	XhoI	1	33
EcoRI	1	663	XmaI	1	1
EcoRV	1	1414	XmnI	1	4744

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

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

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