

pGL4.25[luc2CP/minP] 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 5' end of the cut DNA (the base to the right of the cut site). Vector sequences are also available in the GenBank® database (GenBank®/EMBL Accession Number DQ904457) and on the Internet at: www.promega.com/vectors/.

Restriction Enzymes That Cut the pGL4.25[luc2CP/minP] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	EcoICRI	1	24
AccI	2	2247, 3927	EcoRI	1	1791
AflIII	1	2946	EcoRV	1	42
AgeI	1	1273	FseI	1	1997
Alw26I	1	4260	FspI	1	964
Alw44I	1	2810	HincII	3	1347, 2138, 2248
AvaI	4	34, 688, 1354, 1864	HindIII	1	66
AvaII	3	200, 1446, 3556	HpaI	2	1347, 2138
BalI	3	1123, 1648, 3652	Hsp92I	4	1090, 1538, 1733, 3955
BamHI	1	2240	KasI	1	1089
BanII	2	26, 172	KpnI	1	19
BbeI	1	1093	NarI	1	1090
BbsI	2	200, 2325	NcoI	1	140
BglI	2	9, 60	NheI	1	28
BglII	1	47	NotI	1	3316
BlpI	1	1096	NspI	3	633, 888, 2500
BsaBI	2	589, 2239	PleI	5	1990, 2390, 2875, 3407, 4030
BsaMI	2	2059, 2152	PshAI	1	2311
BspHI	3	725, 1331, 3 216	PstI	2	1976, 3340
BspMI	1	4433	PvuI	1	3788
BsrGI	1	632	PvuII	1	216
BssSI	2	1069, 2669	SacI	1	26
Bst98I	1	1821	SacII	1	3812
BstEII	1	3343	SalI	1	2246
BstXI	1	3340	ScaI	4	307, 3898, 3990, 4368
BstZI	3	1991, 1995, 3316	SfiI	2	9, 60
Bsu36I	1	3774	SgrAI	1	205
ClaI	2	2233, 4361	SinI	3	200, 1446, 3556
DraI	3	2199, 3255, 3274	SpeI	1	4247
DraIII	2	1288, 1833	SspI	2	4222, 4278
DrdI	2	1543, 2604	StyI	2	140, 755
EarI	3	1450, 2380, 3636	XbaI	1	1978
EclHKI	1	3418	XhoI	1	34
Eco47III	3	235, 346, 2372	XmnI	1	1810

Note: XbaI cuts the vector once in *dam+* strains, as a second XbaI site downstream of the selectable marker is protected by methylation. XbaI cuts twice in *dam-* strains.

Restriction Enzymes That Do Not Cut the pGL4.25[luc2CP/minP] Vector.

AatII	AvrII	BssHII	I-PpoI	NsiI	PpuMI	SphI	VspI
AccB7I	BbuI	Csp45I	MluI	PacI	SgfI	StuI	XcmI
AccIII	BsaAI	CspI	NdeI	PmeI	SmaI	SwaI	XmaI
AscI	BsaI	EcoNI	NruI	PmlI	SnaBI	Tth111I	

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