

pGL4.26[luc2P/minP/Hygro] 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 DQ904458) and on the Internet at: www.promega.com/vectors/.

Restriction Enzymes That Cut the pGL4.26[luc2P/minP/Hygro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	EarI	5	1450, 2089, 2629, 3800, 5056
AccI	4	2816, 3016, 3667, 5347	EclHKI	2	3047, 4838
AflIII	2	2725, 3916	Eco47III	4	235, 346, 2747, 3792
AgeI	2	1273, 3587	EcoICRI	1	24
Alw26I	1	5680	EcoRV	1	42
Alw44I	1	4230	FseI	1	1818
AlwNI	1	4332	FspI	3	964, 2099, 3342
ApaI	2	172, 2953	HincII	3	1347, 1959, 3668
AvaI	5	34, 688, 1354, 3105, 3300	HindIII	1	66
AvaII	5	200, 1446, 3190, 3490, 4976	HpaI	2	1347, 1959
AvrII	1	2495	Hsp92I	4	1090, 1538, 1733, 5375
BalI	4	1123, 1648, 3228, 5072	KasI	1	1089
BamHI	1	2061	KpnI	1	19
BanII	4	26, 172, 2953, 3407	NarI	1	1090
BbeI	1	1093	NcoI	3	140, 2106, 2402
BbsI	3	200, 3103, 3745	NheI	1	28
BbuI	2	2243, 2315	NotI	1	4736
BglI	2	9, 60	NsiI	2	2245, 2317
BglII	1	47	PmeI	1	3576
BlpI	1	1096	PshAI	2	2920, 3731
BsaAI	1	2726	PstI	1	4760
BsaBI	2	589, 2060	PvuI	1	5208
BsaMI	2	1880, 1973	PvuII	2	216, 2171
BspHI	3	725, 1331, 4636	SacI	1	26
BspMI	1	5853	SacII	1	5232
BsrGI	1	632	SalI	1	3666
BssHII	1	3551	ScaI	4	307, 5318, 5410, 5788
BssSI	3	1069, 3372, 4089	SfiI	2	9, 60
BstEII	1	4763	SgrAI	1	205
BstXI	1	4760	SinI	5	200, 1446, 3190, 3490, 4976
BstZI	4	1812, 1816, 3164, 4736	SpeI	1	5667
Bsu36I	1	5194	SphI	2	2243, 2315
ClaI	2	2054, 5781	SspI	2	5642, 5698
Csp45I	1	3652	StuI	1	2494
DraI	4	2020, 3576, 4675, 4694	XbaI	2	1799, 3581
DraIII	1	1288	XhoI	1	34
DrdI	3	1543, 3501, 4024	XmnI	1	2578

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 pGL4.26[luc2P/minP/Hygro] Vector.

AatII	AscI	CspI	I-PpoI	NruI	PpuMI	SnaBI	VspI
AccB7I	BsaI	EcoNI	MluI	PacI	SgfI	Swal	XcmI
AccIII	Bst98I	EcoRI	NdeI	PmlI	SmaI	Tth111I	XmaI

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