

pGL4.18[*luc2P*/Neo] 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 DQ188838) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.18[*luc2P*/Neo] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	Eco47III	4	194, 305, 2747, 3637
AccI	2	3512, 5192	EcoICRI	1	24
AflIII	1	3761	EcoNI	1	3102
AgeI	2	1232, 3432	EcoRI	1	1750
Alw26I	1	5525	EcoRV	1	42
Alw44I	1	4075	FseI	3	1905, 2759, 3079
AlwNI	3	2930, 3119, 4177	FspI	2	923, 2186
ApaI	2	131, 2702	HincII	3	1306, 2046, 3513
AvaI	4	34, 647, 1313, 1769	HindIII	1	66
AvrII	2	2582, 3260	HpaI	2	1306, 2046
BamHI	1	2148	KasI	3	1048, 2790, 2989
BanII	3	26, 131, 2702	KpnI	1	19
BbeI	3	1052, 2794, 2993	NarI	3	1049, 2791, 2990
BbsI	3	159, 2781, 3590	NcoI	3	99, 2193, 2489
BbuI	3	2330, 2402, 3161	NheI	1	28
BglI	2	9, 60	NotI	1	4581
BglII	1	47	NsiI	2	2332, 2404
BlpI	3	1055, 2934, 3037	PpuMI	2	2915, 3173
BsaBI	2	548, 2147	PshAI	2	3253, 3576
BsaMI	2	1967, 2060	PstI	1	4605
BspHI	3	684, 1290, 4481	PvuI	1	5053
BspMI	2	3024, 5698	PvuII	2	175, 2258
BsrGI	1	591	SacI	1	26
BssSI	2	1028, 3934	SacII	1	5077
BstEII	1	4608	SalI	1	3511
BstXI	1	4605	ScaI	4	266, 5163, 5255, 5633
BstZI	5	1899, 1903, 2728, 2888, 4581	SfiI	2	9, 60
Bsu36I	1	5039	SgrAI	2	164, 2844
Clal	2	2141, 5626	SpeI	1	5512
Csp45I	1	3497	SphI	3	2330, 2402, 3161
DraI	3	2107, 4520, 4539	SspI	2	5487, 5543
DraIII	1	1247	StuI	1	2581
DrdI	4	1502, 1889, 3180, 3869	XbaI	2	1886, 3426
EarI	4	1409, 2176, 3645, 4901	XhoI	1	34
EclHKI	1	4683			

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 pGL4.18[*luc2P*/Neo] Vector.

AatII	BbrPI	CspI	NruI	PspAI	SpII	VspI
AccB7I	BsaI	Eco72I	PacI	RsrII	SrfI	XcmI
AccIII	BsaAI	I-PpoI	PfiMI	SgfI	Sse8387I	XmaI
AflII	BssHIII	MluI	PmeI	SmaI	SwaI	XmnI
AscI	Bst98I	NdeI	PmlI	SnaBI	Tth111I	

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