

pGL4.11[*luc2P*] 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 AY738223) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.11[*luc2P*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	Eco47III	3	194, 305, 2280
AccI	2	2155, 3835	EcoICRI	1	24
AflIII	1	2404	EcoRI	1	1750
AgeI	1	1232	EcoRV	1	42
Alw26I	1	4168	FseI	1	1905
Alw44I	1	2718	FspI	1	923
AlwNI	1	2820	HincII	3	1306, 2046, 2156
ApaI	1	131	HindIII	1	66
AvaI	4	34, 647, 1313, 1769	HpaI	2	1306, 2046
AvaII	3	159, 1405, 3464	Hsp92I	4	1049, 1497, 1692, 3863
Ball	3	1082, 1607, 3560	KasI	1	1048
BamHI	1	2148	KpnI	1	19
BanII	2	26, 131	NarI	1	1049
BbeI	1	1052	NcoI	1	99
BbsI	2	159, 2233	NheI	1	28
BglI	2	9, 60	NotI	1	3224
BglII	1	47	NspI	3	592, 847, 2408
BlpI	1	1055	PshAI	1	2219
BsaBI	2	548, 2147	PstI	1	3248
BsaMI	2	1967, 2060	PvuI	1	3696
BspHI	3	684, 1290, 3124	PvuII	1	175
BspMI	1	4341	SacI	1	26
BsrGI	1	591	SacII	1	3720
BssSI	2	1028, 2577	SalI	1	2154
BstEII	1	3251	ScaI	4	266, 3806, 3898, 4276
BstXI	1	3248	SfiI	2	9, 60
BstZI	3	1899, 1903, 3224	SgrAI	1	164
Bsu36I	1	3682	SinI	3	159, 1405, 3464
Clal	2	2141, 4269	SpeI	1	4155
DraI	3	2107, 3163, 3182	SspI	2	4130, 4186
DraIII	1	1247	StyI	2	99, 714
DrdI	3	1502, 1889, 2512	XbaI	1	1886
EarI	3	1409, 2288, 3544	XhoI	1	34
EclHKI	1	3326			

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 pGL4.11[*luc2P*] Vector.

AatII	BbrPI	CspI	NdeI	PmlI	SgfI	Sse8387I	XmaI
AccB7I	BbuI	Csp45 I	NruI	Ppu10I	SmaI	StuI	XmnI
AccIII	BsaI	Eco72I	NsiI	PpuMI	SnaBI	Swal	
AflII	BsaAI	EcoNI	PacI	Psp5 II	SphI	Tth111I	
AscI	BssHIII	I-PpoI	PflMI	PspAI	SpII	VspI	
AvrII	Bst98I	MluI	PmeI	RsrII	SrfI	XcmI	

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