

pGL4.10[*luc2*] 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 AY738222) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.10[*luc2*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65 I	1	15	EcoICRI	1	24
AccI	2	2027, 3707	EcoRV	1	42
AflIII	1	2276	FseI	1	1777
AgeI	1	1232	FspI	1	923
Alw26I	1	4040	HincII	3	1306, 1918, 2028
Alw44I	1	2590	HindIII	1	66
AlwNI	1	2692	HpaI	2	1306, 1918
ApaI	1	131	Hsp92I	4	1049, 1497, 1692, 3735
AvaI	3	34, 647, 1313	KasI	1	1048
AvaII	3	159, 1405, 3336	KpnI	1	19
BalI	3	1082, 1607, 3432	NaeI	5	166, 1510, 1531, 1775, 2146
BamHI	1	2020	NarI	1	1049
BanII	2	26, 131	NcoI	1	99
BbeI	1	1052	NgoMIV	5	164, 1508, 1529, 1773, 2144
BbsI	2	159, 2105	NheI	1	28
BglII	2	9, 60	NotI	1	3096
BglIII	1	47	NspI	3	592, 847, 2280
BlpI	1	1055	PleI	5	1770, 2170, 2655, 3187, 3810
BsaBI	2	548, 2019	PshAI	1	2091
BsaMI	2	1839, 1932	PstI	1	3120
BspHI	3	684, 1290, 2996	PvuI	1	3568
BspMI	1	4213	PvuII	1	175
BsrGI	1	591	SacI	1	26
BssSI	2	1028, 2449	SacII	1	3592
BstEII	1	3123	SalI	1	2026
BstXI	1	3120	ScaI	4	266, 3678, 3770, 4148
BstZI	3	1771, 1775, 3096	SfiI	2	9, 60
Bsu36I	1	3554	SgrAI	1	164
ClaI	2	2013, 4141	SinI	3	159, 1405, 3336
DraI	3	1979, 3035, 3054	SpeI	1	4027
DraIII	1	1247	SspI	2	4002, 4058
DrdI	2	1502, 2384	StyI	2	99, 714
EarI	3	1409, 2160, 3416	XbaI	1	1758
EclHKI	1	3198	XhoI	1	34
Eco47III	3	194, 305, 2152			

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 pGL4.10[*luc2*] Vector.

AatII	BbrPI	CspI	MluI	PmeI	RsrII	SrfI	XcmI
AccB7I	BbuI	Csp45I	NdeI	PmlI	SgfI	Sse 8387I	XmaI
AccIII	BsaI	Eco72I	NruI	Ppu10I	SmaI	StuI	XmnI
AflII	BsaAI	EcoNI	NsiI	PpuMI	SnaBI	SwaI	
AscI	BssHIII	EcoRI	PacI	Psp5II	SphI	Tth111I	
AvrII	Bst98I	I-PpoI	PflMI	PspAI	SplI	VspI	

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