

pGL4.23[*luc2*/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 DQ904455) and on the Internet at: www.promega.com/vectors/.

Restriction Enzymes That Cut the pGL4.23[*luc2*/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	2068, 3748	EcoRV	1	42
AflIII	1	2317	FseI	1	1818
AgeI	1	1273	FspI	1	964
Alw26I	1	4081	HincII	3	1347, 1959, 2069
Alw44I	1	2631	HindIII	1	66
AlwNI	1	2733	HpaI	2	1347, 1959
ApaI	1	172	Hsp92I	4	1090, 1538, 1733, 3776
AvaI	3	34, 688, 1354	KasI	1	1089
AvaII	3	200, 1446, 3377	KpnI	1	19
BalI	3	1123, 1648, 3473	NaeI	5	207, 1551, 1572, 1816, 2187
BamHI	1	2061	NarI	1	1090
BanII	2	26, 172	NcoI	1	140
BbeI	1	1093	NgoMIV	5	205, 1549, 1570, 1814, 2185
BbsI	2	200, 2146	NheI	1	28
BglI	2	9, 60	NotI	1	3137
BglII	1	47	NspI	3	633, 888, 2321
BlpI	1	1096	PleI	5	1811, 2211, 2696, 3228, 3851
BsaBI	2	589, 2060	PshAI	1	2132
BsaMI	2	1880, 1973	PstI	1	3161
BspHI	3	725, 1331, 3037	PvuI	1	3609
BspMI	1	4254	PvuII	1	216
BsrGI	1	632	SacI	1	26
BssSI	2	1069, 2490	SacII	1	3633
BstEII	1	3164	SalI	1	2067
BstXI	1	3161	ScaI	4	307, 3719, 3811, 4189
BstZI	3	1812, 1816, 3137	SfiI	2	9, 60
Bsu36I	1	3595	SgrAI	1	205
ClaI	2	2054, 4182	SinI	3	200, 1446, 3377
DraI	3	2020, 3076, 3095	SpeI	1	4068
DraIII	1	1288	SspI	2	4043, 4099
DrdI	2	1543, 2425	StyI	2	140, 755
EarI	3	1450, 2201, 3457	XbaI	1	1799
EclHKI	1	3239	XhoI	1	34
Eco47III	3	235, 346, 2193			

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 pGL4.23[*luc2*/minP] Vector.

AatII	BbuI	Csp45I	NdeI	PmlI	SphI	XcmI
AccB7I	BsaAI	CspI	NruI	PpuMI	StuI	XmaI
AccIII	BsaI	EcoNI	NsiI	SgfI	Swal	XmnI
AscI	BssHII	I-PpoI	Pacl	SmaI	Tth111I	
AvrII	Bst98I	MluI	PmeI	SnaBI	VspI	

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