

pGL4.12[*luc2CP*] 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 AY738224) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.12[*luc2CP*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	Eco47III	3	194, 305, 2331
AccI	2	2206, 3886	EcoICRI	1	24
AflIII	1	2455	EcoRI	1	1750
AgeI	1	1232	EcoRV	1	42
Alw26I	1	4219	FseI	1	1956
Alw44I	1	2769	FspI	1	923
AlwNI	1	2871	HincII	3	1306, 2097, 2207
ApaI	1	131	HindIII	1	66
AvaI	4	34, 647, 1313, 1823	HpaI	2	1306, 2097
AvaII	3	159, 1405, 3515	Hsp92I	4	1049, 1497, 1692, 3914
BalI	3	1082, 1607, 3611	KasI	1	1048
BamHI	1	2199	KpnI	1	19
BanII	2	26, 131	NarI	1	1049
BbeI	1	1052	NcoI	1	99
BbsI	2	159, 2284	NheI	1	28
BglI	2	9, 60	NotI	1	3275
BglIII	1	47	NspI	3	592, 847, 2459
BlpI	1	1055	PleI	5	1949, 2349, 2834, 3366, 3989
BsaBI	2	548, 2198	PshAI	1	2270
BsaMI	2	2018, 2111	PstI	2	1935, 3299
BspHI	3	684, 1290, 3175	PvuI	1	3747
BspMI	1	4392	PvuII	1	175
BsrGI	1	591	SacI	1	26
BssSI	2	1028, 2628	SacII	1	3771
Bst98I	1	1780	SalI	1	2205
BstEII	1	3302	ScaI	4	266, 3857, 3949, 4327
BstXI	1	3299	SfiI	2	9, 60
BstZI	3	1950, 1954, 3275	SgrAI	1	164
Bsu36I	1	3733	SinI	3	159, 1405, 3515
ClaI	2	2192, 4320	SpeI	1	4206
DraI	3	2158, 3214, 3233	SspI	2	4181, 4237
DraIII	2	1247, 1792	StyI	2	99, 714
DrdI	2	1502, 2563	XbaI	1	1937
EarI	3	1409, 2339, 3595	XhoI	1	34
EclHKI	1	3377	XmnI	1	1769

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 pGL4.12[*luc2CP*] Vector.

AatII	BbuI	Eco72I	NsiI	PpuMI	SnaBI	Swal	XcmI
AccB7I	BsaI	EcoNI	PacI	Psp5II	SphI	Tth111I	XmaI
AccIII	BsaAI	I-PpoI	PflMI	PspI	SpII	VspI	
AscI	BssHII	MluI	PmeI	RsrII	SrfI		
AvrII	CspI	NdeI	PmlI	SgfI	Sse8387I		
BbrPI	Csp45I	NruI	Ppu10I	SmaI	StuI		

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