

pGL4.13[*luc2*/SV40] 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 AY738225) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.13[*luc2*/SV40] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	2	15,99	EcoICRI	1	24
AccI	2	2426,4106	EcoRV	1	42
AflIII	1	2675	FseI	1	2176
AgeI	1	1631	FspI	2	53, 1322
Alw26I	1	4439	HincII	3	1705, 2317, 2427
Alw44I	1	2989	HindIII	1	465
AlwNI	1	3091	HpaI	2	1705, 2317
ApaI	1	530	Hsp92I	4	1448, 1896, 2091, 4134
AvaI	3	34, 1046, 1712	KasI	1	1447
AvaII	3	558, 1804, 3735	KpnI	2	19, 103
AvrII	1	449	NaeI	5	565, 1909, 1930, 2174, 2545
BalI	3	1481, 2006, 3831	NarI	1	1448
BamHI	1	2419	NcoI	3	60, 356, 498
BanII	2	26, 530	NgoMIV	5	563, 1907, 1928, 2172, 2543
BbeI	1	1451	NheI	1	28
BbsI	2	558, 2504	NotI	1	3495
BbuI	2	197, 269	NsiI	2	199, 271
BglI	2	9, 402	NspI	5	197, 269, 991, 1246, 2679
BglII	1	46	PleI	5	2169, 2569, 3054, 3586, 4209
BlpI	1	1454	PshAI	1	2490
BsaBI	2	947, 2418	PstI	1	3519
BsaMI	2	2238, 2331	PvuI	1	3967
BspHI	3	1083, 1689, 3395	PvuII	2	125,574
BspMI	1	4612	SacI	1	26
BsrGI	1	990	SacII	1	3991
BssSI	2	1427, 2848	SalI	1	2425
BstEII	1	3522	ScaI	4	665, 4077, 4169, 4547
BstXI	1	3519	SfiI	2	9, 402
BstZI	3	2170, 2174, 3495	SgrAI	1	563
Bsu36I	1	3953	SinI	3	558, 1804, 3735
ClaI	2	2412,4540	SpeI	1	4426
DraI	3	2378, 3434, 3453	SphI	2	197, 269
DraIII	1	1646	SspI	2	4401, 4457
DrdI	2	1901, 2783	StuI	1	448
EarI	3	1808, 2559, 3815	StyI	5	60, 356, 449, 498, 1113
EclHKI	1	3597	XbaI	1	2157
Eco47III	3	593,704,2551	XhoI	1	34

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 pGL4.13[*luc2*/SV40] Vector.

AatII	BbrPI	CspI	I-PpoI	PflMI	PspAI	SplI	VspI
AccB7I	BsaI	Csp45I	MluI	PmeI	RsrII	SrfI	XcmI
AccIII	BsaAI	Eco72I	NdeI	PmlI	SgfI	Sse8387I	XmaI
AflII	BssHIII	EcoNI	NruI	PpuMI	SmaI	Swal	XmnI
AscI	Bst98I	EcoRI	PacI	Psp5II	SnaBI	Tth111I	

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