

pGL4.24[luc2P/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 DQ904456) and on the Internet at: www.promega.com/vectors/.

Restriction Enzymes That Cut the pGL4.24[luc2P/minP] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	Eco47III	3	235, 346, 2321
AccI	2	2196, 3876	EcoICRI	1	24
AflIII	1	2445	EcoRI	1	1791
AgeI	1	1273	EcoRV	1	42
Alw26I	1	4209	FseI	1	1946
Alw44I	1	2759	FspI	1	964
AlwNI	1	2861	HincII	3	1347, 2087, 2197
ApaI	1	172	HindIII	1	66
AvaI	4	34, 688, 1354, 1810	HpaI	2	1347, 2087
AvaII	3	200, 1446, 3505	Hsp92I	4	1090, 1538, 1733, 3904
BalI	3	1123, 1648, 3601	KasI	1	1089
BamHI	1	2189	KpnI	1	19
BanII	2	26, 172	NarI	1	1090
BbeI	1	1093	NcoI	1	140
BbsI	2	200, 2274	NheI	1	28
BglI	2	9, 60	NotI	1	3265
BglIII	1	47	NspI	3	633, 888, 2449
BlpI	1	1096	PshAI	1	2260
BsaBI	2	589, 2188	PstI	1	3289
BsaMI	2	2008, 2101	PvuI	1	3737
BspHI	3	725, 1331, 3165	PvuII	1	216
BspMI	1	4382	SacI	1	26
BsrGI	1	632	SacII	1	3761
BssSI	2	1069, 2618	SalI	1	2195
BstEII	1	3292	ScaI	4	307, 3847, 3939, 4317
BstXI	1	3289	SfiI	2	9, 60
BstZI	3	1940, 1944, 3265	SgrAI	1	205
Bsu36I	1	3723	SinI	3	200, 1446, 3505
ClaI	2	2182, 4310	SpeI	1	4196
DraI	3	2148, 3204, 3223	SspI	2	4171, 4227
DraIII	1	1288	StyI	2	140, 755
DrdI	3	1543, 1930, 2553	XbaI	1	1927
EarI	3	1450, 2329, 3585	XhoI	1	34
EclHKI	1	3367			

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 pGL4.24[luc2P/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.