

pGL4.27[*luc2P*/minP/Hygro] 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 DQ904459) and on the Internet at: www.promega.com/vectors/.

Restriction Enzymes That Cut the pGL4.27[*luc2P*/minP/Hygro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	EclHKI	2	3175, 4966
AccI	4	2944, 3144, 3795, 5475	Eco47III	4	235, 346, 2875, 3920
AflIII	2	2853, 4044	EcoICRI	1	24
AgeI	2	1273, 3715	EcoRI	1	1791
Alw26I	1	5808	EcoRV	1	42
Alw44I	1	4358	FseI	1	1946
AlwNI	1	4460	FspI	3	964, 2227, 3470
ApaI	2	172, 3081	HincII	3	1347, 2087, 3796
AvaII	5	200, 1446, 3318, 3618, 5104	HindIII	1	66
AvrII	1	2623	HpaI	2	1347, 2087
BalI	4	1123, 1648, 3356, 5200	Hsp92I	4	1090, 1538, 1733, 5503
BamHI	1	2189	KasI	1	1089
BanII	4	26, 172, 3081, 3535	KpnI	1	19
BbeI	1	1093	NarI	1	1090
BbsI	3	200, 3231, 3873	NcoI	3	140, 2234, 2530
BbuI	2	2371, 2443	NheI	1	28
BglI	2	9, 60	NotI	1	4864
BglII	1	47	NsiI	2	2373, 2445
BlpI	1	1096	Pme I	1	3704
BsaAI	1	2854	PshAI	2	3048, 3859
BsaBI	2	589, 2188	PstI	1	4888
BsaMI	2	2008, 2101	PvuI	1	5336
BspHI	3	725, 1331, 4764	PvuII	2	216, 2299
BspMI	1	5981	SacI	1	26
BsrGI	1	632	SacII	1	5360
BssHII	1	3679	SalI	1	3794
BssSI	3	1069, 3500, 4217	Scal	4	307, 5446, 5538, 5916
BstEII	1	4891	SfiI	2	9, 60
BstXI	1	4888	SgrAI	1	205
BstZI	4	1940, 1944, 3292, 4864	SinI	5	200, 1446, 3318, 3618, 5104
Bsu36I	1	5322	SpeI	1	5795
ClaI	2	2182, 5909	SphI	2	2371, 2443
Csp45I	1	3780	SspI	2	5770, 5826
DraI	4	2148, 3704, 4803, 4822	StuI	1	2622
DraIII	1	1288	XbaI	2	1927, 3709
DrdI	4	1543, 1930, 3629, 4152	XhoI	1	34
EarI	5	1450, 2217, 2757, 3928, 5184	XmnI	1	2706

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 pGL4.27[*luc2P*/minP/Hygro] Vector.

AatII	AscI	CspI	MluI	PacI	SgfI	Swal	XcmI
AccB7I	BsaI	EcoNI	NdeI	PmlI	SmaI	Tth111I	XmaI
AccIII	Bst98I	I-PpoI	NruI	PpuMI	SnaBI	VspI	

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