

pGL4.77[hRlucP/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 AY864932) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.77[hRlucP/Hygro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	EclHKI	2	2417, 4208
Acc65I	1	15	Eco47III	2	2117, 3162
AccI	4	2186, 2386, 3037, 4717	EcoICRI	1	24
AflIII	2	2095, 3286	EcoRI	1	1033
AgeI	1	2957	EcoRV	2	42, 586
Alw26I	2	641, 5050	FseI	1	1188
Alw44I	1	3600	FspI	2	1469, 2712
AlwNI	1	3702	HincII	2	1329, 3038
ApaI	1	2323	HindIII	1	66
AvaI	5	34, 122, 1052, 2475, 2670	HpaI	1	1329
AvaII	3	2560, 2860, 4346	Hsp92I	2	795,4745
AvrII	1	1865	KpnI	1	19
BalI	2	2598, 4442	NciI	3	3666, 4520, 4742
BamHI	1	1431	NcoI	3	99, 1476, 1772
BanII	5	26, 166, 306, 2323, 2777	NheI	1	28
BbsI	2	2473, 3115	NotI	1	4106
BbuI	2	1613, 1685	NruI	1	762
BclI	4	141, 594, 4146, 4529	NsiI	3	320, 1615, 1687
BglI	3	9, 60, 101	NspI	4	1613, 1685, 2708, 3290
BglII	1	47	PmeI	1	2946
BsaAI	1	2096	PshAI	2	2290, 3101
BsaBI	2	860, 1430	PstI	1	4130
BsaI	1	641	PvuI	1	4578
BsaMI	2	1250, 1343	PvuII	1	1541
BspHI	1	4006	SacI	1	26
BspMI	1	5223	SacII	1	4602
BssHII	1	2921	SalI	1	3036
BssSI	2	2742, 3459	ScaI	3	4688, 4780, 5158
BstEII	1	4133	SfiI	2	9,60
BstXI	1	4130	SinI	3	2560, 2860, 4346
BstZI	4	1182, 1186, 2534, 4106	SpeI	1	5037
Bsu36I	2	687, 4564	SphI	2	1613,1685
ClaI	2	1424, 5151	SspI	2	5012, 5068
Csp45I	1	3022	StuI	1	1864
DraI	4	1390, 2946, 4045, 4064	Tth111I	1	797
DraIII	1	289	XbaI	2	1169, 2951
DrdI	3	1172, 2871, 3394	XhoI	1	34
EarI	5	600, 1459, 1999, 3170, 4426	XmnI	2	635,1948

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 pGL4.77[hRlucP/Hygro] Vector.

AccB7I	BbrPI	CspI	KasI	PflMI	RsrII	SplI	XcmI
AccIII	BlpI	Eco72I	MluI	PmlI	SgfI	SrfI	XmaI
AflII	Bpu 1102I	EcoNI	NarI	PpuMI	SgrAI	Sse8387I	
AscI	BsrGI	EheI	NdeI	Psp5II	SmaI	SwaI	
BbeI	Bst98I	I-PpoI	PacI	PspAI	SnaBI	VspI	

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