

pGL4.78[*hRlucCP/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 AY864933) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.78[*hRlucCP/Hygro*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	EcIHKI	2	2468, 4259
Acc65I	1	15	Eco47III	2	2168, 3213
AccI	4	2237, 2437, 3088, 4768	EcoICRI	1	24
AflIII	2	2146, 3337	EcoRI	1	1033
AgeI	1	3008	EcoRV	2	42, 586
Alw26I	2	641, 5101	FseI	1	1239
Alw44I	1	3651	FspI	2	1520, 2763
AlwNI	1	3753	HincII	2	1380, 3089
ApaI	1	2374	HindIII	1	66
AvaI	5	34, 122, 1106, 2526, 2721	HpaI	1	1380
AvaII	3	2611, 2911, 4397	Hsp92I	2	795, 4796
AvrII	1	1916	KpnI	1	19
BalI	2	2649, 4493	NciI	3	3717, 4571, 4793
BamHI	1	1482	NcoI	3	99, 1527, 1823
BanII	5	26, 166, 306, 2374, 2828	NheI	1	28
BbsI	2	2524, 3166	NotI	1	4157
BbuI	2	1664, 1736	NruI	1	762
BclI	4	141, 594, 4197, 4580	NsiI	3	320, 1666, 1738
BglI	3	9, 60, 101	NspI	4	1664, 1736, 2759, 3341
BglII	1	47	PmeI	1	2997
BsaAI	1	2147	PshAI	2	2341, 3152
BsaBI	2	860, 1481	PstI	2	1218, 4181
BsaI	1	641	PvuI	1	4629
BsaMI	2	1301, 1394	PvuII	1	1592
BspHI	1	4057	SacI	1	26
BspMI	1	5274	SacII	1	4653
BssHII	1	2972	SalI	1	3087
BssSI	2	2793, 3510	ScaI	3	4739, 4831, 5209
Bst98I	1	1063	SfiI	2	9, 60
BstEII	1	4184	SinI	3	2611, 2911, 4397
BstXI	1	4181	SpeI	1	5088
BstZI	4	1233, 1237, 2585, 4157	SphI	2	1664, 1736
Bsu36I	2	687, 4615	SspI	2	5063, 5119
ClaI	2	1475, 5202	StuI	1	1915
Csp45I	1	3073	Tth111I	1	797
DraI	4	1441, 2997, 4096, 4115	XbaI	2	1220, 3002
DraIII	2	289, 1075	XhoI	1	34
DrdI	2	2922, 3445	XmnI	3	635, 1052, 1999
EarI	5	600, 1510, 2050, 3221, 4477			

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 pGL4.78[*hRlucCP/Hygro*] Vector.

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

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