

pGL4.76[*hRluc*/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 AY864931) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.76[*hRluc*/Hygro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	EarI	5	600, 1328, 1868, 3039, 4295
Acc65I	1	15	EclHKI	2	2286, 4077
AccI	4	2055, 2255, 2906, 4586	Eco47III	2	1986, 3031
AflIII	2	1964, 3155	EcoICRI	1	24
AgeI	1	2826	EcoRV	2	42, 586
Alw26I	2	641,4919	FseI	1	1057
Alw44I	1	3469	FspI	2	1338, 2581
AlwNI	1	3571	HincII	2	1198, 2907
ApaI	1	2192	HindIII	1	66
AvaI	4	34, 122, 2344, 2539	HpaI	1	1198
AvaII	3	2429, 2729, 4215	Hsp92I	2	795, 4614
AvrII	1	1734	KpnI	1	19
BalI	2	2467, 4311	NciI	3	3535, 4389, 4611
BamHI	1	1300	NcoI	3	99, 1345, 1641
BanI	5	15, 2683, 3020, 4025, 5104	NheI	1	28
BanII	5	26, 166, 306, 2192, 2646	NotI	1	3975
BbsI	2	2342, 2984	NruI	1	762
BbuI	2	1482, 1554	NsiI	3	320, 1484, 1556
BclI	4	141, 594, 4015, 4398	NspI	4	1482, 1554, 2577, 3159
BglII	3	9, 60, 101	PmeI	1	2815
BglIII	1	47	PshAI	2	2159, 2970
BsaAI	1	1965	PstI	1	3999
BsaBI	2	860, 1299	PvuI	1	4447
BsaI	1	641	PvuII	1	1410
BsaMI	2	1119, 1212	SacI	1	26
BspHI	1	3875	SacII	1	4471
BspMI	1	5092	SalI	1	2905
BssHII	1	2790	ScaI	3	4557, 4649, 5027
BssSI	2	2611, 3328	SfiI	2	9, 60
BstEII	1	4002	SinI	3	2429, 2729, 4215
BstXI	1	3999	SpeI	1	4906
BstZI	4	1051, 1055, 2403, 3975	SphI	2	1482, 1554
Bsu36I	2	687, 4433	SspI	2	4881, 4937
ClaI	2	1293, 5020	StuI	1	1733
Csp45I	1	2891	Tth111I	1	797
DraI	4	1259, 2815, 3914, 3933	XbaI	2	1038, 2820
DraIII	1	289	XhoI	1	34
DrdI	2	2740, 3263	XmnI	2	635, 1817

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 pGL4.76[*hRluc*/Hygro] Vector.

AccB7I	BbrPI	Csp I	I-PpoI	PacI	PspAI	SnaBI	VspI
AccIII	BlpI	Eco72 I	KasI	PflMI	RsrII	SpII	XcmI
AflII	Bpu 1102I	EcoN I	MluI	PmlI	SgfI	SrfI	XmaI
AscI	BsrGI	EcoR I	NarI	PpuMI	SgrAI	Sse8387I	
BbeI	Bst98I	Ehe I	NdeI	Psp5II	SmaI	SwaI	

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