

pGL4.14[*luc2*/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 AY864928) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.14[*luc2*/Hygro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	EarI	5	1409, 2048, 2588, 3759, 5015
AccI	4	2775, 2975, 3626, 5306	EclHKI	2	3006, 4797
AflIII	2	2684, 3875	Eco47III	4	194, 305, 2706, 3751
AgeI	2	1232, 3546	EcoICRI	1	24
Alw26I	1	5639	EcoRV	1	42
Alw44I	1	4189	FseI	1	1777
AlwNI	1	4291	FspI	3	923, 2058, 3301
ApaI	2	131, 2912	HincII	3	1306, 1918, 3627
AvaI	5	34, 647, 1313, 3064, 3259	HindIII	1	66
AvaII	5	159, 1405, 3149, 3449, 4935	HpaI	2	1306, 1918
AvrII	1	2454	Hsp92I	4	1049, 1497, 1692, 5334
BalI	4	1082, 1607, 3187, 5031	KasI	1	1048
BamHI	1	2020	KpnI	1	19
BanII	4	26, 131, 2912, 3366	NarI	1	1049
BbeI	1	1052	NcoI	3	99, 2065, 2361
BbsI	3	159, 3062, 3704	NheI	1	28
BbuI	2	2202, 2274	NotI	1	4695
BglI	2	9, 60	NsiI	2	2204, 2276
BglII	1	47	PmeI	1	3535
BlpI	1	1055	PshAI	2	2879, 3690
BsaAI	1	2685	PstI	1	4719
BsaBI	2	548, 2019	PvuI	1	5167
BsaMI	2	1839, 1932	PvuII	2	175, 2130
BspHI	3	684, 1290, 4595	SacI	1	26
BspMI	1	5812	SacII	1	5191
BsrGI	1	591	SalI	1	3625
BssHII	1	3510	Scal	4	266, 5277, 5369, 5747
BssSI	3	1028, 3331, 4048	SfiI	2	9, 60
BstEII	1	4722	SgrAI	1	164
BstXI	1	4719	SinI	5	159, 1405, 3149, 3449, 4935
BstZI	4	1771, 1775, 3123, 4695	SpeI	1	5626
Bsu36I	1	5153	SphI	2	2202, 2274
ClaI	2	2013, 5740	SspI	2	5601, 5657
Csp45I	1	3611	StuI	1	2453
DraI	4	1979, 3535, 4634, 4653	XbaI	2	1758, 3540
DraIII	1	1247	XhoI	1	34
DrdI	3	1502, 3460, 3983	XmnI	1	2537

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 pGL4.14[*luc2*/Hygro] Vector.

AatII	BbrPI	EcoNI	NruI	Psp5II	SnaBI	Tth111I
AccB7I	BsaI	EcoRI	PacI	PspAI	SplI	VspI
AccIII	Bst98I	I-PpoI	PflMI	RsrII	SrfI	XcmI
AflIII	CspI	MluI	PmlI	SgfI	Sse8387I	XmaI
AscI	Eco72I	NdeI	PpuMI	SmaI	Swa	

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