

pGL4.15[*luc2P*/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 AY864929) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.15[*luc2P*/Hygro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	EclHKI	2	3134, 4925
AccI	4	2903, 3103, 3754, 5434	Eco47III	4	194, 305, 2834, 3879
AflIII	2	2812, 4003	EcoICRI	1	24
AgeI	2	1232, 3674	EcoRI	1	1750
Alw26I	1	5767	EcoRV	1	42
Alw44I	1	4317	FseI	1	1905
AlwNI	1	4419	FspI	3	923, 2186, 3429
ApaI	2	131, 3040	HincII		1306, 2046, 3755
AvaII	5	159, 1405, 3277, 3577, 5063	HindIII	1	66
AvrII	1	2582	HpaI	2	1306, 2046
BalI	4	1082, 1607, 3315, 5159	Hsp92I	4	1049, 1497, 1692, 5462
BamHI	1	2148	KasI	1	1048
BanII	4	26, 131, 3040, 3494	KpnI	1	19
BbeI	1	1052	NarI	1	1049
BbsI	3	159, 3190, 3832	NcoI	3	99, 2193, 2489
BbuI	2	2330, 2402	NheI	1	28
BglI	2	9, 60	NotI	1	4823
BglII	1	47	NsiI	2	2332, 2404
BlpI	1	1055	PmeI	1	3663
BsaAI	1	2813	PshAI	2	3007, 3818
BsaBI	2	548, 2147	PstI	1	4847
BsaMI	2	1967, 2060	PvuI	1	5295
BspHI	3	684, 1290, 4723	PvuII	2	175, 2258
BspMI		5940	SacI	1	26
BsrGI	1	591	SacII	1	5319
BssHII	1	3638	SalI	1	3753
BssSI	3	1028, 3459, 4176	ScaI	4	266, 5405, 5497, 5875
BstEII	1	4850	SfiI	2	9, 60
BstXI	1	4847	SgrAI	1	164
BstZI	4	1899, 1903, 3251, 4823	SinI	5	159, 1405, 3277, 3577, 5063
Bsu36I	1	5281	SpeI	1	5754
ClaI	2	2141, 5868	SphI	2	2330, 2402
Csp45I	1	3739	SspI	2	5729, 5785
DraI	4	2107, 3663, 4762, 4781	StuI	1	2581
DraIII	1	1247	XbaI	2	1886, 3668
DrdI	4	1502, 1889, 3588, 4111	XhoI	1	34
EarI	5	1409, 2176, 2716, 3887, 5143	XmnI	1	2665

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 pGL4.15[*luc2P*/Hygro] Vector.

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

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