

pGL4.28[*luc2CP/minP/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 DQ904460) and on the Internet at: www.promega.com/vectors/.

Restriction Enzymes That Cut the pGL4.28[*luc2CP/minP/Hygro*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	EclHKI	2	3226, 5017
AccI	4	2995, 3195, 3846, 5526	Eco47III	4	235, 346, 2926, 3971
AflIII	2	2904, 4095	EcoICRI	1	24
AgeI	2	1273, 3766	EcoRI	1	1791
Alw26I	1	5859	EcoRV	1	42
Alw44I	1	4409	FseI	1	1997
AlwNI	1	4511	FspI	3	964, 2278, 3521
ApaI	2	172, 3132	HincII	3	1347, 2138, 3847
AvaII	5	200, 1446, 3369, 3669, 5155	HindIII	1	66
AvrII	1	2674	HpaI	2	1347, 2138
BalI	4	1123, 1648, 3407, 5251	Hsp92I	4	1090, 1538, 1733, 5554
BamHI	1	2240	KasI	1	1089
BanII	4	26, 172, 3132, 3586	KpnI	1	19
BbeI	1	1093	NarI	1	1090
BbsI	3	200, 3282, 3924	NcoI	3	140, 2285, 2581
BbuI	2	2422, 2494	NheI	1	28
BglI	2	9, 60	NotI	1	4915
BglII	1	47	NsiI	2	2424, 2496
BlpI	1	1096	PmeI	1	3755
BsaAI	1	2905	PshAI	2	3099, 3910
BsaBI	2	589, 2239	PstI	2	1976, 4939
BsaMI	2	2059, 2152	PvuI	1	5387
BspHI	3	725, 1331, 4815	PvuII	2	216, 2350
BspMI	1	6032	SacI	1	26
BsrGI	1	632	SacII	1	5411
BssHII	1	3730	SalI	1	3845
BssSI	3	1069, 3551, 4268	ScaI	4	307, 5497, 5589, 5967
Bst98I	1	1821	SfiI	2	9, 60
BstEII	1	4942	SgrAI	1	205
BstXI	1	4939	SinI	5	200, 1446, 3369, 3669, 5155
BstZI	4	1991, 1995, 3343, 4915	SpeI	1	5846
Bsu36I	1	5373	SphI	2	2422, 2494
ClaI	2	2233, 5960	SspI	2	5821, 5877
Csp45I	1	3831	StuI	1	2673
DraI	4	2199, 3755, 4854, 4873	XbaI	2	1978, 3760
DraIII	2	1288, 1833	XhoI	1	34
DrdI	3	1543, 3680, 4203	XmnI	2	1810, 2757
EarI	5	1450, 2268, 2808, 3979, 5235			

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 pGL4.28[*luc2CP/minP/Hygro*] Vector.

AatII	AscI	EcoNI	NdeI	PmlI	SmaI	Tth111I	XmaI
AccB7I	BsaI	I-PpoI	NruI	PpuMI	SnaBI	VspI	
AccIII	CspI	MluI	Pacl	SgfI	Swal	Xcml	

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