

pGL4.22[*luc2CP/Puro*] 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 DQ188842) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.22[*luc2CP/Puro*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	EcoICRI	1	24
AccI	2	3368, 5048	EcoRI	1	1750
AflIII	1	3617	EcoRV	1	42
AgeI	2	1232, 3288	FseI	1	1956
Alw26I	1	5381	FspI	2	923, 2237
Alw44I	1	3931	HincII	3	1306, 2097, 3369
AlwNI	2	2831, 4033	HindIII	1	66
ApaI	1	131	HpaI	2	1306, 2097
AvrII	2	2633, 3240	KasI	3	1048, 2931, 3270
BalI	5	1082, 1607, 2709, 3054, 4773	KpnI	1	19
BamHI	1	2199	NarI	3	1049, 2932, 3271
BanII	3	26, 131, 3072	NcoI	3	99, 2244, 2540
BbeI	3	1052, 2935, 3274	NheI	1	28
BbsI	2	159, 3446	NotI	1	4437
BbuI	2	2381, 2453	NruI	1	2716
BglI	2	9, 60	NsiI	2	2383, 2455
BglII	1	47	NspI	5	592, 847, 2381, 2453, 3621
BlpI	2	1055, 2970	PleI	5	1949, 3511, 3996, 4528, 5151
BsaBI	2	548, 2198	PpuMI	2	3237, 3244
BsaMI	2	2018, 2111	PshAI	2	2844, 3432
BspHI	3	684, 1290, 4337	PstI	2	1935, 4461
BspMI	1	5554	PvuI	1	4909
BsrGI	1	591	PvuII	2	175, 2309
BssSI	2	1028, 3790	SacI	1	26
Bst98I	1	1780	SacII	1	4933
BstEII	1	4464	SalI	1	3367
BstXI	1	4461	ScaI	4	266, 5019, 5111, 5489
BstZI	4	1950, 1954, 2890, 4437	SfiI	2	9, 60
Bsu36I	1	4895	SgrAI	2	164, 3138
Clal	2	2192, 5482	SpeI	1	5368
Csp45I	1	3353	SphI	2	2381, 2453
DraI	3	2158, 4376, 4395	SspI	2	5343, 5399
DraIII	3	1247, 1792, 2786	StuI	1	2632
DrdI	2	1502, 3725	XbaI	2	1937, 3282
EarI	4	1409, 2227, 3501, 4757	XhoI	1	34
EclHKI	1	4539	XmnI	1	1769
Eco47III	3	194, 305, 3493			

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 pGL4.22[*luc2CP/Puro*] Vector.

AatII	BbrPI	CspI	MluI	PmeI	SgfI	SrfI	VspI
AccB7I	BsaI	Eco72I	NdeI	PmlI	SmaI	Sse8387I	XcmI
AccIII	BsaAI	EcoNI	PacI	PspAI	SnaBI	Swal	XmaI
AscI	BssHII	I-PpoI	PflMI	RsrII	SplI	Tth111I	

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