

pGL4.21[*luc2P/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 DQ188841) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.21[*luc2P/Puro*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	Eco47III	3	194, 305, 3442
AccI	2	3317, 4997	EcoICRI	1	24
AflIII	1	3566	EcoRI	1	1750
AgeI	2	1232, 3237	EcoRV	1	42
Alw26I	1	5330	FseI	1	1905
Alw44I	1	3880	FspI	2	923, 2186
AlwNI	2	2780, 3982	HincII	3	1306, 2046, 3318
ApaI	1	131	HindIII	1	66
AvrII	2	2582, 3189	HpaI	2	1306, 2046
BalI	5	1082, 1607, 2658, 3003, 4722	KasI	3	1048, 2880, 3219
BamHI	1	2148	KpnI	1	19
BanII	3	26, 131, 3021	NarI	3	1049, 2881, 3220
BbeI	3	1052, 2884, 3223	NcoI	3	99, 2193, 2489
BbsI	2	159, 3395	NheI	1	28
BbuI	2	2330, 2402	NotI	1	4386
BglI	2	9, 60	NruI	1	2665
BglII	1	47	NsiI	2	2332, 2404
BlpI	2	1055, 2919	NspI	5	592, 847, 2330, 2402, 3570
BsaBI	2	548, 2147	PpuMI	2	3186, 3193
BsaMI	2	1967, 2060	PshAI	2	2793, 3381
BspHI	3	684, 1290, 4286	PstI	1	410
BspMI	1	5503	PvuI	1	4858
BsrGI	1	591	PvuII	2	175, 2258
BssSI	2	1028, 3739	SacI	1	26
BstEII	1	4413	SacII	1	4882
BstXI	1	4410	SalI	1	3316
BstZI	4	1899, 1903, 2839, 4386	ScaI	4	266, 4968, 5060, 5438
Bsu36I	1	4844	SfiI	2	9, 60
Clal	2	2141, 5431	SgrAI	2	164, 3087
Csp45I	1	3302	SpeI	1	5317
DraI	3	2107, 4325, 4344	SphI	2	2330, 2402
DraIII	2	1247, 2735	SspI	2	5292, 5348
DrdI	3	1502, 1889, 3674	StuI	1	2581
EarI	4	1409, 2176, 3450, 4706	XbaI	2	1886, 3231
EclHKI	1	4488	XhoI	1	34

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 pGL4.21[*luc2P/Puro*] Vector.

AatII	BbrPI	CspI	NdeI	PspAI	SpII	VspI
AccB7I	BsaI	Eco72I	PacI	RsrII	SrfI	XcmI
AccIII	BsaAI	EcoNI	PflMI	SgfI	Sse8387I	XmaI
AflII	BssHII	I-PpoI	PmeI	Sma	Swal	XmnI
AscI	Bst98I	MluI	PmlI	SnaBI	Tth111I	

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