

pGL4.20[*luc2/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 DQ188840) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.20[*luc2/Puro*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
Acc65I	1	15	Eco47III	3	194, 305, 3314
AccI	2	3189, 4869	EcoICRI	1	24
AflIII	1	3438	EcoRV	1	42
AgeI	2	1232, 3109	FseI	1	1777
Alw26I	1	5202	FspI	2	923, 2058
Alw44I	1	3752	HincII	3	1306, 1918, 3190
AlwNI	2	2652, 3854	HindIII	1	66
ApaI	1	131	HpaI	2	1306, 1918
AvaI	5	34, 647, 1313, 2734, 3052	KasI	3	1048, 2752, 3091
AvrII	2	2454, 3061	KpnI	1	19
BalI	5	1082, 1607, 2530, 2875, 4594	NarI	3	1049, 2753, 3092
BamHI	1	2020	NcoI	3	99, 2065, 2361
BanII	3	26, 131, 2893	NheI	1	28
BbeI	3	1052, 2756, 3095	NotI	1	4258
BbsI	2	159, 3267	NruI	1	2537
BbuI	2	2202, 2274	NsiI	2	2204, 2276
BglI	2	9, 60	NspI	5	592, 847, 2202, 2274, 3442
BglII	1	47	PleI	5	1770, 3332, 3817, 4349, 4972
BlpI	2	1055, 2791	PpuMI	2	3058, 3065
BsaBI	2	548, 2019	PshAI	2	2665, 3253
BsaMI	2	1839, 1932	PstI	1	4282
BspHI	3	684, 1290, 4158	PvuI	1	4730
BspMI	1	5375	PvuII	2	175, 2130
BsrGI	1	591	SacI	1	26
BssSI	2	1028, 3611	SacII	1	4754
BstEII	1	4285	SalI	1	3188
BstXI	1	4282	ScaI	4	266, 4840, 4932, 5310
BstZI	4	1771, 1775, 2711, 4258	SfiI	2	9, 60
Bsu36I	1	4716	SgrAI	2	164, 2959
ClaI	2	2013, 5303	SpeI	1	5189
Csp45I	1	3174	SphI	2	2202, 2274
DraI	3	1979, 4197, 4216	SspI	2	5164, 5220
DraIII	2	1247, 2607	StuI	1	2453
DrdI	2	1502, 3546	XbaI	2	1758, 3103
EarI	4	1409, 2048, 3322, 4578	XhoI	1	34
EclHKI	1	4360			

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 pGL4.20[*luc2/Puro*] Vector.

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

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