

pGL3-Promoter 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 3'-end of the cut DNA (the base to the left of the cut site). Vector sequences are also available in the GenBank® database (GenBank®/EMBL Accession Number U47298) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL3-Promoter Vector 1-5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccI	1	2203	EcoICRI	1	9
AccIII	2	975, 1491	EcoNI	3	837, 1237, 1897
Acc65I	1	1	EheI	1	314
AcyI	4	28, 313, 1706, 3882	FseI	1	1953
AflIII	3	15, 773, 2452	FspI	2	3567, 4740
Alw26I	5	1303, 1535, 1601, 3406, 4182	HincII	3	1584, 2094, 2204
Alw44I	2	2766, 4012	HindII	3	1584, 2094, 2204
AlwNI	1	2868	HindIII	1	245
AspHI	5	11, 1745, 2770, 3931, 4016	HpaI	1	2094
AvaI	3	26, 32, 1336	Hsp92I	4	287, 313, 1706, 3882
AvaII	3	1459, 3483, 3705	KasI	1	312
AvrII	1	229	KpnI	1	5
BamHI	1	2196	MluI	1	15
BanII	4	11, 33, 1304, 4423	NaeI	3	1951, 2322, 4391
BbeI	1	316	NarI	1	313
BbsI	4	290, 1568, 1684, 2281	NcoI	1	278
BbuI	1	943	NgoMIV	3	1949, 2320, 4389
BclI	1	860	NheI	1	21
BglI	3	182, 3465, 4733	NotI	1	4843
BglII	1	36	NspI	2	943, 2456
BsaI	1	3406	PaeR7I	2	1867, 4458
BsaAI	1	4494	PpuMI	1	1459
BsaBI	2	48, 2195	PshAI	1	2267
BsaHI	4	287, 313, 1706, 3882	Psp5II	1	1459
BsaMI	3	252, 2015, 2108	PspAI	1	26
BsmI	3	252, 2015, 2108	PvuI	2	3715, 4761
BspHI	3	863, 3172, 4180	SacI	1	11
BspMI	3	1669, 1678, 4973	SalI	1	2202
BsrGI	1	770	ScaI	3	445, 3825, 4908
BssSI	2	2625, 4009	SfiI	1	182
BstZI	3	1947, 1951, 4843	SgrAI	1	1708
ClaI	3	2189, 4901, 5005	SinI	3	1459, 3483, 3705
Csp45I	1	449	SmaI	1	28
DraI	4	2155, 3211, 3230, 3922	SphI	1	943
DraII	1	1459	SrfI	1	28
DraIII	1	4497	SspI	3	4149, 4702, 4817
DrdI	3	1681, 2560, 4541	StuI	1	228
DsaI	2	278, 650	StyI	2	229, 278
EaeI	4	1947, 1951, 3733, 4843	VspI	1	3517
EagI	3	1947, 1951, 4843	XbaI	1	1934
EclHKI	1	3345	XcmI	1	1015
Eco47III	1	2328	XhoI*	1	32
Eco52I	3	1947, 1951, 4843	XmaI	1	26
			XmnI	1	3944

*Due to the extent of supercoiling in this vector, the XhoI site has proven difficult to cut to completion under standard restriction digest conditions. For single XhoI digests, we recommend digesting the vector for a minimum of 2 hours using 20 units of enzyme per microgram of DNA at 37°C to ensure complete digestion. If performing a double digest with XhoI and another enzyme, linearize the vector using the companion enzyme prior to carrying out the XhoI digest. Under these conditions, XhoI will cut the vector following standard reactions conditions.

Restriction Enzymes That Do Not Cut the pGL3-Promoter Vector.

AatII	BalI	Bst1107I	Eco72I	NruI	PmlI	SgfI	Tth111I
AccB7I	BbrPI	Bst98I	Eco81I	NsiI	Ppu10I	SnaBI	
AflIII	BlpI	BstEII	EcoRI	PacI	PstI	SpeI	
AgeI	Bpu1102I	BstXI	EcoRV	PfIMI	PvuII	SpII	
ApaI	Bsp120I	Bsu36I	I-PpoI	PinAI	RsrII	Sse8387I	
AscI	BssHII	CspI	NdeI	PmeI	SacII	Swal	

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