

pGL3-Control 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 U47296) 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	EcoNI	3	837, 1237, 1897
AccIII	2	975, 1491	EheI	1	314
Acc65I	1	1	FseI	1	1953
AcyI	4	28, 313, 1706, 3882	FspI	2	3567, 4740
AflIII	3	15, 773, 2452	HincII	3	1584, 2094, 2204
Alw26I	5	1303, 1535, 1601, 3406, 4182	HindIII	3	1584, 2094, 2204
Alw44I	2	2766, 4012	HindIII	1	245
AlwNI	1	2868	HpaI	1	2094
AspHI	5	11, 1745, 2770, 3931, 4016	Hsp92I	4	287, 313, 1706, 3882
AvaI	3	26, 32, 1336	KasI	1	312
AvaII	3	1459, 3483, 3705	KpnI	1	5
AvrII	1	229	MluI	1	15
BamHI	1	2196	NaeI	3	1951, 2322, 4391
BanII	4	11, 33, 1304, 4423	NarI	1	313
BbeI	1	316	NcoI	1	278
BbsI	4	290, 1568, 1684, 2281	NgoMIV	3	1949, 2320, 4389
BbuI	1	943	NheI	1	21
BclI	1	860	NotI	1	4843
BglI	3	182, 3465, 4733	NspI	2	943, 2456
BglII	1	36	PaeR7I	2	1867, 4458
BsaI	1	3406	PpuMI	1	1459
BsaAI	1	4494	PshAI	1	2267
BsaBI	2	48, 2195	Psp5II	1	1459
BsaHI	4	287, 313, 1706, 3882	PspAI	1	26
BsaMI	3	252, 2015, 2108	PvuI	2	3715, 4761
BsmI	3	252, 2015, 2108	SacI	1	11
BspHI	3	863, 3172, 4180	SalI	1	2202
BspMI	3	1669, 1678, 4973	ScaI	3	445, 3825, 4908
BsrGI	1	770	SfiI	1	182
BssSI	2	2625, 4009	SgrAI	1	1708
BstZI	3	1947, 1951, 4843	SinI	3	1459, 3483, 3705
ClaI	3	2189, 4901, 5005	SmaI	1	28
Csp45I	1	449	SphI	1	943
DraI	4	2155, 3211, 3230, 3922	SrfI	1	28
DraII	1	1459	SspI	3	4149, 4702, 4817
DraIII	1	4497	StuI	1	228
DrdI	3	1681, 2560, 4541	StyI	2	229, 278
DsaI	2	278, 650	VspI	1	3517
EaeI	4	1947, 1951, 3733, 4843	XbaI	1	1934
EagI	3	1947, 1951, 4843	XcmI	1	1015
EclHKI	1	3345	XhoI*	1	32
Eco47III	1	2328	XmaI	1	26
Eco52I	3	1947, 1951, 4843	XmnI	1	3944
EcoICRI	1	9			

*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-Control 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	BssHIII	CspI	NdeI	PmeI	SacII	Swal	

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