

pGL3-Enhancer 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 U47297) and on the Internet at: www.promega.com/vectors/

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

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AccI	1	2257	EheI	1	122
AccIII	2	783,1299	FseI	1	1761
Acc65I	1	1	FspI	2	3621, 4794
AcyI	4	95, 121, 1514, 3936	HincII	3	1392, 1902, 2258
AflIII	3	15, 581, 2506	HindII	3	1392, 1902, 2258
Alw26I	5	1111, 1343, 1409, 3460, 4236	HindIII	1	53
Alw44I	2	2820, 4066	HpaI	1	1902
AlwNI	1	2922	Hsp92I	4	95, 121, 1514, 3936
AspHI	5	11, 1553, 2824, 3985, 4070	KasI	1	120
AvaI	3	26, 32, 1144	KpnI	1	5
AvaII	3	1267, 3537, 3759	MluI	1	15
BamHI	1	2250	NaeI	3	1759, 2376, 4445
BanII	4	11, 33, 1112, 4477	NarI	1	121
BbeI	1	124	NcoI	1	86
BbsI	4	98, 1376, 1492, 2335	NgoMIV	3	1757, 2374, 4443
BbuI	3	751, 2108, 2180	NheI	1	21
BclI	1	668	NotI	1	4897
BglI	2	3519, 4787	NsiI	2	2106, 2178
BglII	1	36	NspI	4	751, 2108, 2180, 2510
BsaI	1	3460	PaeR7I	2	1675, 4512
BsaAI	1	4548	Ppu10I	2	2102, 2174
BsaBI	1	2003	PpuMI	1	1267
BsaHI	4	95, 121, 1514, 3936	PshAI	1	2321
BsaMI	3	60, 1823, 1916	Psp5II	1	1267
BsmI	3	60, 1823, 1916	PspAI	1	26
BspHI	3	671, 3226, 4234	PvuI	2	3769, 4815
BspMI	3	1477, 1486, 5027	SacI	1	11
BsrGI	1	578	SalI	1	2256
BssSI	2	2679, 4063	ScaI	3	253, 3879, 4962
BstZI	3	1755, 1759, 4897	SgrAI	1	1516
ClaI	3	1997, 4955, 5059	SinI	3	1267, 3537, 3759
Csp45I	1	257	SmaI	1	28
DraI	4	1963, 3265, 3284, 3976	SphI	3	751, 2108, 2180
DraII	1	1267	SrfI	1	28
DraIII	1	4551	SspI	3	4203, 4756, 4871
DrdI	3	1489, 2614, 4595	StyI	1	86
DsaI	2	86, 458	VspI	1	3571
EaeI	4	1755, 1759, 3787, 4897	XbaI	1	1742
EagI	3	1755, 1759, 4897	XcmI	1	823
EclHKI	1	3399	XhoI*	1	32
Eco47III	1	2382	XmaI	1	26
Eco52I	3	1755, 1759, 4897	XmnI	1	3998
EcoICRI	1	9			
EcoNI	3	645, 1045, 1705			

*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-Enhancer Vector.

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

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