

## pGL2-Basic 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 X65323) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pGL2-Basic Vector 1-5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
<b>AccB7I</b>	1	2044	<b>EcoRV</b>	1	1414
<b>AccI</b>	1	2746	EheI	1	110
<b>AccIII</b>	2	771, 1287	FspI	2	4110, 5283
<b>Acc65I</b>	1	8	<b>HaellI</b>	5	112, 2873, 3243, 4882, 4890
AcyI	4	83, 109, 1502, 4425	<b>HincII</b>	4	1380, 2608, 2747, 5451
AflIII	3	22, 569, 2995	HindIII	4	1380, 2608, 2747, 5451
<b>Alw26I</b>	4	1099, 1331, 3949, 4725	<b>HindIII</b>	1	47
<b>Alw44I</b>	2	3309, 4555	<b>HpaI</b>	2	2608, 5451
AlwNI	1	3411	<b>Hsp92I</b>	4	83, 109, 1502, 4425
AspHI	5	18, 1541, 3313, 4474, 4559	KasI	1	108
<b>AvaI</b>	3	1, 33, 1132	<b>KpnI</b>	1	12
<b>AvaII</b>	3	1255, 4026, 4248	<b>MluI</b>	1	22
<b>BamHI</b>	1	2739	<b>MspA1I</b>	5	139, 1429, 3337, 3582, 4523
<b>BanII</b>	3	18, 1100, 4966	<b>NaeI</b>	2	2865, 4934
BbeI	1	112	<b>NarI</b>	1	109
BbsI	5	86, 1364, 1480, 2192, 2824	<b>NgoMIV</b>	2	2863, 4932
<b>BbuI</b>	1	739	<b>NheI</b>	1	28
<b>BglI</b>	2	4008, 5276	NspI	2	739, 2999
<b>BglII</b>	1	37	Pacl	1	1400
BsaI	1	3949	PaeR7I	1	33
BsaAI	3	230, 5037, 5370	PfIMI	1	2044
BsaBI	1	2507	PpuMI	1	1255
BsaHI	4	83, 109, 1502, 4425	PshAI	1	2810
BspHI	2	3715, 4723	Psp5II	1	1255
BspMI	1	1474	PspAI	1	1
BsrGI	1	566	<b>PvuI</b>	2	4258, 5304
BssSI	2	3168, 4552	<b>SacI</b>	1	18
<b>BstEII</b>	1	683	<b>SalI</b>	1	2745
<b>Bsu36I</b>	1	689	<b>ScaI</b>	1	4368
Cfr10I	5	345, 1504, 2863, 3968, 4932	SgrAI	1	1504
<b>ClaI</b>	1	1441	<b>SinI</b>	3	1255, 4026, 4248
<b>Csp45I</b>	2	245, 1033	<b>SmaI</b>	1	3
DraII	1	1255	<b>SphI</b>	1	739
DraIII	1	5040	SpII	1	230
DrdI	2	3103, 5084	<b>SspI</b>	5	2323, 2476, 4692, 5245, 5360
EaeI	1	4276	<b>StyI</b>	1	2192
<b>EclHKI</b>	1	3888	<b>VspI</b>	2	2411, 4060
<b>Eco47III</b>	1	2871	<b>XbaI</b>	1	123
Eco81I	1	689	XcmI	1	811
<b>EcoICRI</b>	1	16	<b>XhoI</b>	1	33
EcoNI	1	1693	<b>XmaI</b>	1	1
<b>EcoRI</b>	1	663	<b>XmnI</b>	1	4487

### Restriction Enzymes That Do Not Cut the pGL2-Basic Vector.

<b>AatII</b>	BbrPI	<b>Bst98I</b>	Eco72I	<b>NsiI</b>	RsrII	Sse8387I
AflII	<b>BclI</b>	<b>BstXI</b>	FseI	PinAI	<b>SacII</b>	<b>StuI</b>
<b>AgeI</b>	BlpI	<b>BstZI</b>	<b>I-PpoI</b>	PmeI	<b>SfiI</b>	Swal
<b>ApaI</b>	Bpu1102I	<b>CspI</b>	NcoI	PmlI	<b>Sgfl</b>	<b>Tth111I</b>
AscI	Bsp120I	DsaI	<b>NdeI</b>	Ppu10I	<b>SnaBI</b>	
AvrII	<b>BssHIII</b>	EagI	<b>NotI</b>	<b>PstI</b>	<b>SpeI</b>	
<b>Ball</b>	Bst1107I	Eco52I	<b>NruI</b>	<b>PvuII</b>	SrfI	

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