

pGL4.80[*hRlucP/Neo*] 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 DQ188844) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.80[*hRlucP/Neo*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	EcoNI	1	2385
Acc65I	1	15	EcoRI	1	1033
AccI	2	2795, 4475	EcoRV	2	42, 586
AflIII	1	3044	FseI	3	1188, 2042, 2362
AgeI	1	2715	FspI	1	1469
Alw26I	2	641, 4808	HgaI	4	977, 3155, 3733, 4492
Alw44I	1	3358	HincII	2	1329, 2796
AlwNI	3	2213, 2402, 3460	HindIII	1	66
ApaI	1	1985	HpaI	1	1329
AvaI	3	34, 122, 1052	Hsp92I	4	795, 2074, 2273, 4503
AvrII	2	1865, 2543	KasI	2	2073, 2272
BalI	4	2121, 2206, 2580, 4200	KpnI	1	19
BamHI	1	1431	NarI	2	2074, 2273
BanII	4	26, 166, 306, 1985	NciI	4	2175, 3424, 4278, 4500
BbeI	2	2077, 2276	NcoI	3	99, 1476, 1772
BbsI	2	2064, 2873	NheI	1	28
BbuI	3	1613, 1685, 2444	NotI	1	3864
BclI	4	141, 594, 3904, 4287	NruI	1	762
BglI	3	9, 60, 101	NsiI	3	320, 1615, 1687
BglII	1	47	NspI	4	1613, 1685, 2444, 3048
BlpI	2	2217, 2320	PpuMI	2	2198, 2456
BsaBI	2	860, 1430	PshAI	2	2536, 2859
BsaI	1	641	PstI	1	3888
BsaMI	2	1250, 1343	PvuI	1	4336
BspHI	1	3764	PvuII	1	1541
BspMI	2	2307, 4981	SacI	1	26
BssSI	1	3217	SacII	1	4360
BstEII	1	3891	SalI	1	2794
BstXI	1	3888	ScaI	3	4446, 4538, 4916
BstZI	5	1182, 1186, 2011, 2171, 3864	SfiI	2	9, 60
Bsu36I	2	687, 4322	SgrAI	1	2127
ClaI	2	1424, 4909	SpeI	1	4795
Csp45I	1	2780	SphI	3	1613, 1685, 2444
DraI	3	1390, 3803, 3822	SspI	2	4770, 4826
DraIII	1	289	StuI	1	1864
DrdI	3	1172, 2463, 3152	Tth111I	1	797
EarI	4	600, 1459, 2928, 4184	XbaI	2	1169, 2709
EclHKI	1	3966	XhoI	1	34
Eco47III	2	2030, 2920	XmnI	1	635
EcoICRI	1	24			

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 pGL4.80[*hRlucP/Neo*] Vector.

AccB7I	BbrPI	Bst98I	MluI	PmeI	SgfI	SrfI	XcmI
AccIII	BsaAI	CspI	NdeI	PmlI	SmaI	Sse8387I	XmaI
AflIII	BsrGI	Eco72I	PacI	PspAI	SnaBI	Swal	
AscI	BssHIII	I-PpoI	PflMI	RsrII	SpII	VspI	

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