

## pGL4.79[*hRluc*/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 DQ188843) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pGL4.79[*hRluc*/Neo] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
<b>AatII</b>	1	798	EcoNI	1	2254
<b>Acc65I</b>	1	15	<b>EcoRV</b>	2	42, 586
<b>AccI</b>	2	2664, 4344	FseI	3	1057, 1911, 2231
AflIII	1	2913	FspI	1	1338
<b>AgeI</b>	1	2584	HgaI	4	977, 3024, 3602, 4361
<b>Alw26I</b>	2	641, 4677	HincII	2	1198, 2665
<b>Alw44I</b>	1	3227	<b>HindIII</b>	1	66
AlwNI	3	2082, 2271, 3329	<b>HpaI</b>	1	1198
<b>ApaI</b>	1	1854	<b>Hsp92I</b>	4	795, 1943, 2142, 4372
<b>AvaI</b>	2	34, 122	KasI	2	1942, 2141
AvrII	2	1734, 2412	<b>KpnI</b>	1	19
<b>BalI</b>	4	1990, 2075, 2449, 4069	<b>NaeI</b>	5	1055, 1909, 2040, 2229, 2783
<b>BamHI</b>	1	1300	<b>NarI</b>	2	1943, 2142
<b>BanII</b>	4	26, 166, 306, 1854	<b>NciI</b>	4	2044, 3293, 4147, 4369
BbeI	2	1946, 2145	<b>NcoI</b>	3	99, 1345, 1641
BbsI	2	1933, 2742	<b>NgoMIV</b>	5	1053, 1907, 2038, 2227, 2781
<b>BbuI</b>	3	1482, 1554, 2313	<b>NheI</b>	1	28
<b>BclI</b>	4	141, 594, 3773, 4156	<b>NotI</b>	1	3733
<b>BglI</b>	3	9, 60, 101	<b>NruI</b>	1	762
<b>BglIII</b>	1	47	<b>NsiI</b>	3	320, 1484, 1556
BlpI	2	2086, 2189	NspI	4	1482, 1554, 2313, 2917
BsaBI	2	860, 1299	PpuMI	2	2067, 2325
BsaI	1	641	PshAI	2	2405, 2728
<b>BsaMI</b>	2	1119, 1212	<b>PstI</b>	1	3757
BspHI	1	3633	<b>PvuI</b>	1	4205
BspMI	2	2176, 4850	<b>PvuII</b>	1	1410
BssSI	1	3086	<b>SacI</b>	1	26
<b>BstEII</b>	1	3760	<b>SacII</b>	1	4229
<b>BstXI</b>	1	3757	<b>SalI</b>	1	2663
<b>BstZI</b>	5	1051, 1055, 1880, 2040, 3733	<b>ScaI</b>	3	4315, 4407, 4785
<b>Bsu36I</b>	2	687, 4191	<b>SfiI</b>	2	9, 60
<b>ClaI</b>	2	1293, 4778	SgrAI	1	1996
<b>Csp45I</b>	1	2649	<b>SpeI</b>	1	4664
<b>DraI</b>	3	1259, 3672, 3691	<b>SphI</b>	3	1482, 1554, 2313
DraIII	1	289	<b>SspI</b>	2	4639, 4695
DrdI	2	2332, 3021	<b>StuI</b>	1	1733
EarI	4	600, 1328, 2797, 4053	<b>Tth111I</b>	1	797
<b>EclHKI</b>	1	3835	<b>XbaI</b>	2	1038, 2578
<b>Eco47III</b>	2	1899, 2789	<b>XhoI</b>	1	34
<b>EcoICRI</b>	1	24	<b>XmnI</b>	1	635

**Note:** XbaI cuts the vector once in *dam*<sup>+</sup> strains, as a second XbaI site downstream of the selectable marker is protected by methylation. XbaI cuts twice in *dam*<sup>-</sup> strains.

### Restriction Enzymes That Do Not Cut the pGL4.79[*hRluc*/Neo] Vector.

<b>AccB7I</b>	BbrPI	<b>Bst98I</b>	<b>I-PpoI</b>	PflMI	RsrII	SpII	<b>VspI</b>
<b>AccIII</b>	BsaAI	<b>CspI</b>	<b>MluI</b>	PmeI	<b>SgfI</b>	SrfI	XcmI
AflII	BsrGI	Eco72I	<b>NdeI</b>	PmlI	<b>SmaI</b>	Sse8387I	<b>XmaI</b>
AscI	<b>BssHIII</b>	<b>EcoRI</b>	PaCI	PspAI	<b>SnaBI</b>	Swal	

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