

## pGL4.75[*hRluc*/CMV] 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 AY738231) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pGL4.75[*hRluc*/CMV] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
<b>AatII</b>	5	292, 345, 428, 614, 1557	<b>FokI</b>	4	1067, 3384, 3671, 4028
<b>AccI</b>	2	2066, 3746	FseI	1	1816
AflIII	1	2315	<b>HaeII</b>	2	2193, 2563
<b>Alw26I</b>	3	601, 1400, 4079	HgaI	5	693, 1736, 2426, 3004, 3763
<b>Alw44I</b>	1	2629	<b>HincII</b>	3	683, 1957, 2067
AlwNI	1	2731	<b>HindIII</b>	2	762, 825
<b>AvaI</b>	1	881	<b>HpaI</b>	1	1957
<b>AvaII</b>	1	3375	<b>MspAII</b>	5	2657, 2902, 3134, 3539, 3630
<b>BalI</b>	3	24, 78, 3471	<b>NaeI</b>	2	1814, 2185
<b>BamHI</b>	1	2059	<b>NciI</b>	3	2695, 3549, 3771
<b>BanI</b>	4	632, 2180, 3185, 4264	<b>NcoI</b>	2	527, 858
<b>BanII</b>	3	735, 925, 1065	<b>NdeI</b>	1	401
BbsI	1	2144	<b>NgoMIV</b>	2	1812, 2183
<b>BclI</b>	4	900, 1353, 3175, 3558	<b>NotI</b>	1	3135
BsaAI	1	507	<b>NruI</b>	1	1521
BsaBI	2	1619, 2058	<b>NsiI</b>	1	1079
BsaI	1	1400	NspI	1	2319
<b>BsaMI</b>	2	1878, 1971	PshAI	1	2130
BspHI	1	3035	<b>PstI</b>	1	3159
BspMI	1	4252	<b>PvuI</b>	1	3607
BsrGI	1	110	<b>SacI</b>	1	735
BssSI	1	2488	<b>SacII</b>	1	3631
<b>BstEII</b>	1	3162	<b>SalI</b>	1	2065
<b>BstXI</b>	1	3159	<b>ScaI</b>	3	3717, 3809, 4187
<b>BstZI</b>	3	1810, 1814, 3135	<b>SfiI</b>	2	9, 819
<b>Bsu36I</b>	2	1446, 3593	<b>SinI</b>	1	3375
<b>ClaI</b>	2	2052, 4180	SnaBI	1	507
<b>DraI</b>	3	2018, 3074, 3093	<b>SpeI</b>	2	166, 4066
DraIII	1	1048	<b>SspI</b>	4	19, 66, 4041, 4097
DrdI	1	2423	<b>StyI</b>	3	527, 858, 867
EarI	3	1359, 2199, 3455	TfiI	5	971, 2290, 3257, 3833, 4177
EclHKI	1	3237	<b>Tth111I</b>	1	1556
<b>Eco47III</b>	1	2191	<b>VspI</b>	1	174
<b>EcoICRI</b>	1	733	<b>XbaI</b>	1	1797
<b>EcoRV</b>	1	1345	<b>XmnI</b>	1	1394

**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 pGL4.75[*hRluc*/CMV] Vector.

<b>AccB7I</b>	AvrII	Bsp120I	<b>EcoRI</b>	<b>NarI</b>	PmlI	SgrAI	SwaI
<b>AccIII</b>	BbeI	<b>BssHII</b>	EheI	<b>NheI</b>	PpuMI	<b>SmaI</b>	XcmI
<b>Acc65I</b>	BbrPI	<b>Bst98I</b>	FspI	PacI	Psp5II	<b>SphI</b>	<b>XhoI</b>
AflIII	<b>BbuI</b>	<b>CspI</b>	<b>I-PpoI</b>	PaeR7I	PspAI	SpII	<b>XmaI</b>
<b>AgeI</b>	<b>BglII</b>	<b>Csp45I</b>	KasI	PfIMI	<b>PvuII</b>	SrfI	
<b>ApaI</b>	BlpI	Eco72I	<b>KpnI</b>	PinAI	RsrII	Sse8387I	
AscI	Bpu1102I	EcoNI	<b>MluI</b>	PmeI	<b>SgfI</b>	<b>StuI</b>	

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