

## pGL4.73[*hRluc*/SV40] 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 AY738229) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pGL4.73[*hRluc*/SV40] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
<b>AatII</b>	1	1197	<b>FseI</b>	1	1456
<b>Acc65I</b>	2	15,99	<b>FspI</b>	1	53
<b>AccI</b>	2	1706, 3386	<b>HaeII</b>	2	1833, 2203
<b>AflIII</b>	1	1955	<b>HgaI</b>	4	1376, 2066, 2644, 3403
<b>Alw26I</b>	2	1040, 3719	<b>HincII</b>	2	1597, 1707
<b>Alw44I</b>	1	2269	<b>HindIII</b>	1	465
<b>AlwNI</b>	1	2371	<b>HpaI</b>	1	1597
<b>AvaI</b>	2	34, 521	<b>Hsp92I</b>	2	1194, 3414
<b>AvaII</b>	1	3015	<b>KpnI</b>	2	19, 103
<b>AvrII</b>	1	449	<b>NaeI</b>	2	1454, 1825
<b>BalI</b>	1	3111	<b>NciI</b>	3	2335, 3189, 3411
<b>BamHI</b>	1	1699	<b>NcoI</b>	3	60, 356, 498
<b>BanI</b>	5	15, 99, 1820, 2825, 3904	<b>NgoMIV</b>	2	1452, 1823
<b>BanII</b>	3	26, 565, 705	<b>NheI</b>	1	28
<b>BbsI</b>	1	1784	<b>NotI</b>	1	2775
<b>BbuI</b>	2	197, 269	<b>NruI</b>	1	1161
<b>BclI</b>	4	540, 993, 2815, 3198	<b>NsiI</b>	3	199, 271, 719
<b>BglI</b>	3	9, 402, 500	<b>NspI</b>	3	197, 269, 1959
<b>BglIII</b>	1	46	<b>PshAI</b>	1	1770
<b>BsaBI</b>	2	1259, 1698	<b>PstI</b>	1	2799
<b>BsaI</b>	1	1040	<b>PvuI</b>	1	3247
<b>BsaMI</b>	2	1518, 1611	<b>PvuII</b>	1	125
<b>BspHI</b>	1	2675	<b>SacI</b>	1	26
<b>BspMI</b>	1	3892	<b>SacII</b>	1	3271
<b>BssSI</b>	1	2128	<b>SalI</b>	1	1705
<b>BstEII</b>	1	2802	<b>ScaI</b>	3	3357, 3449, 3827
<b>BstXI</b>	1	2799	<b>SfiI</b>	2	9, 402
<b>BstZI</b>	3	1450, 1454, 2775	<b>SinI</b>	1	3015
<b>Bsu36I</b>	2	1086, 3233	<b>SpeI</b>	1	3706
<b>ClaI</b>	2	1692, 3820	<b>SphI</b>	2	197, 269
<b>DraI</b>	3	1658, 2714, 2733	<b>SspI</b>	2	3681, 3737
<b>DraIII</b>	1	688	<b>StuI</b>	1	448
<b>DrdI</b>	1	2063	<b>StyI</b>	5	60, 356, 449, 498, 507
<b>EarI</b>	3	999, 1839, 3095	<b>TfiI</b>	5	611, 1930, 2897, 3473, 3817
<b>EclHKI</b>	1	2877	<b>Tth111I</b>	1	1196
<b>Eco47III</b>	1	1831	<b>XbaI</b>	1	1437
<b>EcoICRI</b>	1	24	<b>XhoI</b>	1	34
<b>EcoRV</b>	2	42, 985	<b>XmnI</b>	1	1034
<b>FokI</b>	5	300, 707, 3024, 3311, 3668			

**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 pGL4.73[*hRluc*/SV40] Vector.

<b>AccB7I</b>	<b>BbeI</b>	<b>BsrGI</b>	<b>EcoNI</b>	<b>NarI</b>	<b>PmlI</b>	<b>SgrAI</b>	<b>SwaI</b>
<b>AccIII</b>	<b>BbrPI</b>	<b>BssHIII</b>	<b>EcoRI</b>	<b>NdeI</b>	<b>PpuMI</b>	<b>SmaI</b>	<b>VspI</b>
<b>AflII</b>	<b>BlpI</b>	<b>Bst98I</b>	<b>EheI</b>	<b>PacI</b>	<b>PspII</b>	<b>SnaBI</b>	<b>XcmI</b>
<b>AgeI</b>	<b>Bpu1102I</b>	<b>CspI</b>	<b>I-PpoI</b>	<b>PflMI</b>	<b>PspAI</b>	<b>SplI</b>	<b>XmaI</b>
<b>ApaI</b>	<b>BsaAI</b>	<b>Csp45I</b>	<b>KasI</b>	<b>PinAI</b>	<b>RsrII</b>	<b>SrfI</b>	
<b>AscI</b>	<b>Bsp120I</b>	<b>Eco72I</b>	<b>MluI</b>	<b>PmeI</b>	<b>SgfI</b>	<b>Sse838I</b>	

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