

## pGL4.70[*hRluc*] 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 AY738226) and on the Internet at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

### Restriction Enzymes That Cut the pGL4.70[*hRluc*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
<b>AatII</b>	1	798	<b>FokI</b>	4	308, 2625, 2912, 3269
<b>Acc65I</b>	1	15	FseI	1	1057
<b>AccI</b>	2	1307, 2987	<b>HaellI</b>	2	1434, 1804
AflIII	1	1556	HgaI	4	977, 1667, 2245, 3004
<b>Alw26I</b>	2	641, 3320	<b>HincII</b>	2	1198, 1308
<b>Alw44I</b>	1	1870	<b>HindIII</b>	1	66
AlwNI	1	1972	<b>HpaI</b>	1	1198
<b>AvaI</b>	2	34, 122	<b>Hsp92I</b>	2	795, 3015
<b>AvaII</b>	1	2616	<b>KpnI</b>	1	19
<b>BalI</b>	1	2712	<b>MspA1I</b>	5	1898, 2143, 2375, 2780, 2871
<b>BamHI</b>	1	1300	<b>NaeI</b>	2	1055, 1426
<b>BanI</b>	4	15, 1421, 2426, 3505	<b>NciI</b>	3	1936, 2790, 3012
<b>BanII</b>	3	26, 166, 306	<b>NcoI</b>	1	99
BbsI	1	1385	<b>NgoMIV</b>	2	1053, 1424
<b>BclI</b>	4	141, 594, 2416, 2799	<b>NheI</b>	1	28
<b>BglI</b>	3	9, 60, 101	<b>NotI</b>	1	2376
<b>BglII</b>	1	47	<b>NruI</b>	1	762
BsaBI	2	860, 1299	<b>NsiI</b>	1	320
BsaI	1	641	NspI	1	1560
<b>BsaMI</b>	2	1119, 1212	PshAI	1	1371
BspHI	1	2276	<b>PstI</b>	1	2400
BspMI	1	3493	<b>PvuI</b>	1	2848
BssSI	1	1729	<b>SacI</b>	1	26
<b>BstEII</b>	1	2403	<b>SacII</b>	1	2872
<b>BstXI</b>	1	2400	<b>SalI</b>	1	1306
<b>BstZI</b>	3	1051, 1055, 2376	<b>ScaI</b>	3	2958, 3050, 3428
<b>Bsu36I</b>	2	687, 2834	<b>SfiI</b>	2	9, 60
<b>ClaI</b>	2	1293, 3421	<b>SinI</b>	1	2616
<b>DraI</b>	3	1259, 2315, 2334	<b>SpeI</b>	1	3307
DraIII	1	289	<b>SspI</b>	2	3282, 3338
DrdI	1	1664	<b>StyI</b>	2	99, 108
EarI	3	600, 1440, 2696	TfiI	5	212, 1531, 2498, 3074, 3418
<b>EclHKI</b>	1	2478	<b>Tth111I</b>	1	797
<b>Eco47III</b>	1	1432	<b>XbaI</b>	1	1038
<b>EcoICRI</b>	1	24	<b>XhoI</b>	1	34
<b>EcoRV</b>	2	42, 586	<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 pGL4.70[*hRluc*] Vector.

<b>AccB7I</b>	BbeI	BsrGI	<b>EcoRI</b>	<b>NdeI</b>	Psp5II	<b>SnaBI</b>	<b>VspI</b>
<b>AccIII</b>	BbrPI	<b>BssHII</b>	EheI	PacI	PspAI	<b>SphI</b>	XcmI
AflII	<b>BbuI</b>	<b>Bst98I</b>	FspI	PflMI	<b>PvuII</b>	SplI	<b>XmaI</b>
<b>AgeI</b>	BlpI	<b>CspI</b>	<b>I-PpoI</b>	PinAI	RsrII	SrfI	
<b>ApaI</b>	Bpu1102I	<b>Csp45I</b>	KasI	PmeI	<b>SgfI</b>	Sse8387I	
AscI	BsaAI	Eco72I	<b>MluI</b>	PmlI	SgrAI	<b>StuI</b>	
AvrII	Bsp120I	EcoNI	<b>NarI</b>	PpuMI	<b>SmaI</b>	Swal	

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