

pGL4.71[*hRlucP*] 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 AY738227) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.71[*hRlucP*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	FokI	4	308, 2756, 3043, 3400
Acc65I	1	15	FseI	1	1188
AccI	2	1438, 3118	HaellI	3	1143, 1565, 1935
AflIII	1	1687	HgaI	4	977, 1798, 2376, 3135
Alw26I	2	641, 3451	HincII	2	1329, 1439
Alw44I	1	2001	HindIII	1	66
AlwNI	1	2103	HpaI	1	1329
AvaI	3	34, 122, 1052	Hsp92I	2	795, 3146
AvaII	1	2747	KpnI	1	19
BalI	1	2843	MspA1I	5	2029, 2274, 2506, 2911, 3002
BamHI	1	1431	NaeI	3	1075, 1186, 1557
BanI	5	15, 1076, 1552, 2557, 3636	NciI	3	2067, 2921, 3143
BanII	3	26, 166, 306	NcoI	1	99
BbsI	1	1516	NgoMIV	3	1073, 1184, 1555
BclI	4	141, 594, 2547, 2930	NheI	1	28
BglI	3	9, 60, 101	NotI	1	2507
BglII	1	47	NruI	1	762
BsaBI	2	860, 1430	NsiI	1	320
BsaI	1	641	NspI	1	1691
BsaMI	2	1250, 1343	PshAI	1	1502
BspHI	1	2407	PstI	1	2531
BspMI	1	3624	PvuI	1	2979
BssSI	1	1860	SacI	1	26
BstEII	1	2534	SacII	1	3003
BstXI	1	2531	SalI	1	1437
BstZI	3	1182, 1186, 2507	ScaI	3	3089, 3181, 3559
Bsu36I	2	687, 2965	SfiI	2	9, 60
Clal	2	1424, 3552	SinI	1	2747
DraI	3	1390, 2446, 2465	SpeI	1	3438
DraIII	1	289	SspI	2	3413, 3469
DrdI	2	1172, 1795	StyI	2	99, 108
EarI	3	600, 1571, 2827	TfiI	5	212, 1662, 2629, 3205, 3549
EclHKI	1	2609	Tth111I	1	797
Eco47III	1	1563	XbaI	1	1169
EcoICRI	1	24	XhoI	1	34
EcoRI	1	1033	XmnI	1	635
EcoRV	2	42, 586			

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 pGL4.71[*hRlucP*] Vector.

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

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