

pGL4.74[*hRluc*/TK] 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 AY738230) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.74[*hRluc*/TK] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	1513	EcoRI	1	670
Acc65I	1	15	EcoRV	1	1301
AccI	3	363, 2022, 3702	FseI	1	1772
AflIII	2	734, 2271	HaeII	4	171, 314, 2149, 2519
Alw26I	3	407, 1356, 4035	HincII	2	1913, 2023
Alw44I	1	2585	HindIII	1	781
AlwNI	2	378, 2687	HpaI	1	1913
AvaI	4	129, 250, 303, 837	Hsp92I	3	311, 1510, 3730
AvrII	1	343	KasI	1	310
BalI	2	149, 3427	KpnI	1	19
BamHI	1	2015	MluI	1	734
BanI	5	15, 310, 2136, 3141, 4220	NaeI	2	1770, 2141
BanII	3	334, 881, 1021	NarI	1	311
BbeI	1	314	NcoI	2	351, 814
BbsI	2	27, 2100	NgoMIV	2	1768, 2139
BclI	4	856, 1309, 3131, 3514	NotI	1	3091
BglI	4	9, 319, 322, 816	NruI	1	1477
BsaAI	1	367	NsiI	1	1035
BsaBI	2	1575, 2014	NspI	2	620, 2275
BsaI	2	407, 1356	PpuMI	1	340
BsaMI	2	1834, 1927	PshAI	1	2086
BspHI	1	2991	PstI	3	530, 767, 3115
BspMI	1	4208	PvuI	1	3563
BssHII	1	262	PvuII	1	552
BssSI	1	2444	SacII	2	320, 3587
Bst98I	2	57, 181	SalI	1	2021
BstEII	1	3118	ScaI	3	3673, 3765, 4143
BstXI	1	3115	SfiI	1	9
BstZI	3	1766, 1770, 3091	SmaI	1	305
Bsu36I	2	1402, 3549	SpeI	1	4022
ClaI	2	2008, 4136	SspI	2	3997, 4053
Csp45I	1	674	StyI	4	343, 351, 814, 823
DraI	3	1974, 3030, 3049	TfiI	5	927, 2246, 3213, 3789, 4133
DraIII	1	1004	Tth111I	1	1512
DrdI	1	2379	XbaI	1	1753
EarI	4	158, 1315, 2155, 3411	XmaI	1	303
EclHKI	1	3193	XmnI	1	1350
Eco47III	1	2147			

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 pGL4.74[*hRluc*/TK] Vector.

AccB7I	BbrPI	Bsp120I	EcoNI	PacI	PmlI	SnaBI	StuI
AccIII	BbuI	BsrGI	FspI	Paer7I	RsrII	SphI	Swal
AgeI	BglII	CspI	I-PpoI	PflMI	SacI	SpII	VspI
ApaI	BlpI	Eco72I	NdeI	PinAI	SgfI	SrfI	XcmI
AscI	Bpu1102I	EcoCRI	NheI	PmeI	SgrAI	Sse8387I	XhoI

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