

pGL4.83[*hRlucP*/Puro] 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 DQ188847) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.83[*hRlucP*/Puro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	EcoRV	2	42, 586
Acc65I	1	15	FokI	5	308, 1716, 3918, 4205, 4562
AccI	2	2600, 4280	FseI	1	1188
AflIII	1	2849	FspI	1	1469
AgeI	1	2520	HgaI	5	977, 2066, 2960, 3538, 4297
Alw26I	2	641, 4613	HincII	2	1329, 2601
Alw44I	1	3163	HindIII	1	66
AlwNI	2	2063, 3265	HpaI	1	1329
AvaI	5	34, 122, 1052, 2145, 2463	Hsp92I	4	795, 2164, 2503, 4308
AvaII	4	2021, 2469, 2476, 3909	KasI	2	2163, 2502
AvrII	2	1865, 2472	KpnI	1	19
BalI	3	1941, 2286, 4005	NaeI	4	1075, 1186, 2162, 2719
BamHI	1	1431	NarI	2	2164, 2503
BanII	4	26, 166, 306, 2304	NciI	5	2009, 2347, 3229, 4083, 4305
BbeI	2	2167, 2506	NcoI	3	99, 1476, 1772
BbsI	1	2678	NgoMIV	4	1073, 1184, 2160, 2717
BbuI	2	1613, 1685	NheI	1	28
BclI	4	141, 594, 3709, 4092	NotI	1	3669
BglI	3	9, 60, 101	NruI	2	762, 1948
BglII	1	47	NsiI	3	320, 1615, 1687
BlpI	1	2202	NspI	3	1613, 1685, 2853
BsaBI	2	860, 1430	PpuMI	2	2469, 2476
BsaI	1	641	PshAI	2	2076, 2664
BsaMI	2	1250, 1343	PstI	1	3693
BspHI	1	3569	PvuI	1	4141
BspMI	1	4786	PvuII	1	1541
BssSI	1	3022	SacI	1	26
BstEII	1	3696	SacII	1	4165
BstXI	1	3693	Sall	1	2599
BstZI	4	1182, 1186, 2122, 3669	ScaI	3	4251, 4343, 4721
Bsu36I	2	687, 4127	SfiI	2	9, 60
Clal	2	1424, 4714	SgrAI	1	2370
Csp45I	1	2585	SinI	4	2021, 2469, 2476, 3909
DraI	3	1390, 3608, 3627	SpeI	1	4600
DraIII	2	289, 2018	SphI	2	1613, 1685
DrdI	2	1172, 2957	SspI	2	4575, 4631
EarI	4	600, 1459, 2733, 3989	StuI	1	1864
EclHKI	1	3771	Tth111I	1	797
Eco47III	1	2725	XbaI	2	1169, 2514
EcoCRI	1	24	XhoI	1	34
EcoRI	1	1033	XmnI	1	635

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 pGL4.83[*hRlucP*/Puro] Vector.

AccB7I	BbrPI	Bst98I	MluI	PmlI	SnaBI	VspI
AccIII	BsaAI	CspI	NdeI	PspAI	SplI	XcmI
AflII	Bsp120I	Eco72I	PacI	RsrII	SrfI	XmaI
ApaI	BsrGI	EcoNI	PflMI	Sgfl	Sse8387I	
AscI	BssHII	I-PpoI	PmeI	SmaI	SwaI	

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