

pGL4.84[*hRlucCP/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 DQ188848) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.84[*hRlucCP/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, 1767, 3969, 4256, 4613
AccI	2	2651, 4331	FseI	1	1239
AflIII	1	2900	FspI	1	1520
AgeI	1	2571	HgaI	5	977, 2117, 3011, 3589, 4348
Alw26I	2	641, 4664	HincII	2	1380, 2652
Alw44I	1	3214	HindIII	1	66
AlwNI	2	2114, 3316	HpaI	1	1380
AvaI	5	34, 122, 1106, 2196, 2514	Hsp92I	4	795, 2215, 2554, 4359
AvaII	4	2072, 2520, 2527, 3960	KasI	2	2214, 2553
AvrII	2	1916, 2523	KpnI	1	19
BalI	3	1992, 2337, 4056	NaeI	4	1129, 1237, 2213, 2770
BamHI	1	1482	NarI	2	2215, 2554
BanII	4	26, 166, 306, 2355	NciI	5	2060, 2398, 3280, 4134, 4356
BbeI	2	2218, 2557	NcoI	3	99, 1527, 1823
BbsI	1	2729	NgoMIV	4	1127, 1235, 2211, 2768
BbuI	2	1664, 1736	NheI	1	28
BclI	4	141, 594, 3760, 4143	NotI	1	3720
BglI	3	9, 60, 101	NruI	2	762, 1999
BglII	1	47	NsiI	3	320, 1666, 1738
BlpI	1	2253	NspI	3	1664, 1736, 2904
BsaBI	2	860, 1481	PpuMI	2	2520, 2527
BsaI	1	641	PshAI	2	2127, 2715
BsaMI	2	1301, 1394	PstI	2	1218, 3744
BspHI	1	3620	PvuI	1	4192
BspMI	1	4837	PvuII	1	1592
BssSI	1	3073	SacI	1	26
Bst98I	1	1063	SacII	1	4216
BstEII	1	3747	SalI	1	2650
BstXI	1	3744	ScaI	3	4302, 4394, 4772
BstZI	4	1233, 1237, 2173, 3720	SfiI	2	9, 60
Bsu36I	2	687, 4178	SgrAI	1	2421
ClaI	2	1475, 4765	SinI	4	2072, 2520, 2527, 3960
Csp45I	1	2636	SpeI	1	4651
DraI	3	1441, 3659, 3678	SphI	2	1664, 1736
DraIII	3	289, 1075, 2069	SspI	2	4626, 4682
DrdI	1	3008	StuI	1	1915
EarI	4	600, 1510, 2784, 4040	Tth111I	1	797
EclHKI	1	3822	XbaI	2	1220, 2565
Eco47III	1	2776	XhoI	1	34
EcoCRI	1	24	XmnI	2	635, 1052
EcoRI	1	1033			

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 pGL4.84[*hRlucCP/Puro*] Vector.

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

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