

pGL4.81[*hRlucCP/Neo*] 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 DQ188845) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.81[*hRlucCP/Neo*] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	EcoICRI	1	24
Acc65I	1	15	EcoNI	1	2436
AccI	2	2846, 4526	EcoRI	1	1033
AflIII	1	3095	EcoRV	2	42, 586
AgeI	1	2766	FseI	3	1239, 2093, 2413
Alw26I	2	641, 4859	FspI	1	1520
Alw44I	1	3409	HgaI	4	977, 3206, 3784, 4543
AlwNI	3	2264, 2453, 3511	HincII	2	1380, 2847
ApaI	1	2036	HindIII	1	66
AvaI	3	34, 122, 1106	HpaI	1	1380
AvrII	2	1916, 2594	Hsp92I	4	795, 2125, 2324, 4554
BalI	4	2172, 2257, 2631, 4251	KasI	2	2124, 2323
BamHI	1	1482	KpnI	1	19
BanII	4	26,166,306,2036	NarI	2	2125, 2324
BbeI	2	2128, 2327	NciI	4	2226, 3475, 4329, 4551
BbsI	2	2115, 2924	NcoI	3	99, 1527, 1823
BbuI	3	1664, 1736, 2495	NheI	1	28
BclI	4	141, 594, 3955, 4338	NotI	1	3915
BglI	3	9, 60, 101	NruI	1	762
BglII	1	47	NsiI	3	320, 1666, 1738
BlpI	2	2268, 2371	NspI	4	1664, 1736, 2495, 3099
BsaBI	2	860, 1481	PpuMI	2	2249, 2507
BsaI	1	641	PshAI	2	2587, 2910
BsaMI	2	1301, 1394	PstI	2	1218, 3939
BspHI	1	3815	PvuI	1	4387
BspMI	2	2358, 5032	PvuII	1	1592
BssSI	1	3268	SacI	1	26
Bst98I	1	1063	SacII	1	4411
BstEII	1	3942	SalI	1	2845
BstXI	1	3939	ScaI	3	4497, 4589, 4967
BstZI	5	1233, 1237, 2062, 2222, 3915	SfiI	2	9, 60
Bsu36I	2	687, 4373	SgrAI	1	2178
ClaI	2	1475, 4960	SpeI	1	4846
Csp45I	1	2831	SphI	3	1664, 1736, 2495
DraI	3	1441, 3854, 3873	SspI	2	4821, 4877
DraIII	2	289, 1075	StuI	1	1915
DrdI	2	2514, 3203	Tth111I	1	797
EarI	4	600,1510,2979,4235	XbaI	2	1220, 2760
EclHKI	1	4017	XhoI	1	34
Eco47III	2	2081, 2971	XmnI	2	635, 1052

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 pGL4.81[*hRlucCP/Neo*] Vector.

AccB7I	BsaAI	Eco72I	PacI	PspAI	SnaBI	Swal
AccIII	BsrGI	I-PpoI	PfiMI	RsrII	SpII	VspI
AscI	BssHIII	MluI	PmeI	SgfI	SrfI	XcmI
BbrPI	CspI	NdeI	PmlI	SmaI	Sse8387I	XmaI

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