

pGL4.82[*hRluc*/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 DQ188846) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pGL4.82[*hRluc*/Puro] Vector Between 1 and 5 Times.

Enzyme	No. of Sites	Location	Enzyme	No. of Sites	Location
AatII	1	798	FokI	5	308, 1585, 3787, 4074, 4431
Acc65I	1	15	FseI	1	1057
AccI	2	2469, 4149	FspI	1	1338
AflIII	1	2718	HgaI	5	977, 1935, 2829, 3407, 4166
AgeI	1	2389	HincII	2	1198, 2470
Alw26I	2	641, 4482	HindIII	1	66
Alw44I	1	3032	HpaI	1	1198
AlwNI	2	1932, 3134	Hsp92I	4	795, 2033, 2372, 4177
AvaI	4	34, 122, 2014, 2332	KasI	2	2032, 2371
AvaII	4	1890, 2338, 2345, 3778	KpnI	1	19
AvrII	2	1734, 2341	NaeI	3	1055, 2031, 2588
BalI	3	1810, 2155, 3874	NarI	2	2033, 2372
BamHI	1	1300	NciI	5	1878, 2216, 3098, 3952, 4174
BanII	4	26, 166, 306, 2173	NcoI	3	99, 1345, 1641
BbeI	2	2036, 2375	NgoMIV	3	1053, 2029, 2586
BbsI	1	2547	NheI	1	28
BbuI	2	1482, 1554	NotI	1	3538
BclI	4	141, 594, 3578, 3961	NruI	2	762, 1817
BglI	3	9, 60, 101	NsiI	3	320, 1484, 1556
BglII	1	47	NspI	3	1482, 1554, 2722
BlpI	1	2071	PpuMI	2	2338, 2345
BsaBI	2	860, 1299	PshAI	2	1945, 2533
BsaI	1	641	PstI	1	3562
BsaMI	2	1119, 1212	PvuI	1	4010
BspHI	1	3438	PvuII	1	1410
BspMI	1	4655	SacI	1	26
BssSI	1	2891	SacII	1	4034
BstEII	1	3565	SalI	1	2468
BstXI	1	3562	ScaI	3	4120, 4212, 4590
BstZI	4	1051, 1055, 1991, 3538	SfiI	2	9, 60
Bsu36I	2	687, 3996	SgrAI	1	2239
Clal	2	1293, 4583	SinI	4	1890, 2338, 2345, 3778
Csp45I	1	2454	SpeI	1	4469
DraI	3	1259, 3477, 3496	SphI	2	1482, 1554
DraIII	2	289, 1887	SspI	2	4444, 4500
DrdI	1	2826	StuI	1	1733
EarI	4	600, 1328, 2602, 3858	Tth111I	1	797
EclHKI	1	3640	XbaI	2	1038, 2383
Eco47III	1	2594	XhoI	1	34
EcoICRI	1	24	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 pGL4.82[*hRluc*/Puro] Vector.

AccB7I	BbrPI	Bst98I	I-PpoI	PmeI	SmaI	Swal
AccIII	BsaAI	CspI	MluI	PmlI	SnaBI	VspI
AflII	Bsp120I	Eco72I	NdeI	PspAI	SpII	XcmI
ApaI	BsrGI	EcoNI	PacI	RsrII	SrfI	XmaI
AscI	BssHIII	EcoRI	PfiMI	SgfI	Sse8387I	

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