

pTK-neo Plasmid

Cat. No. 71284-3

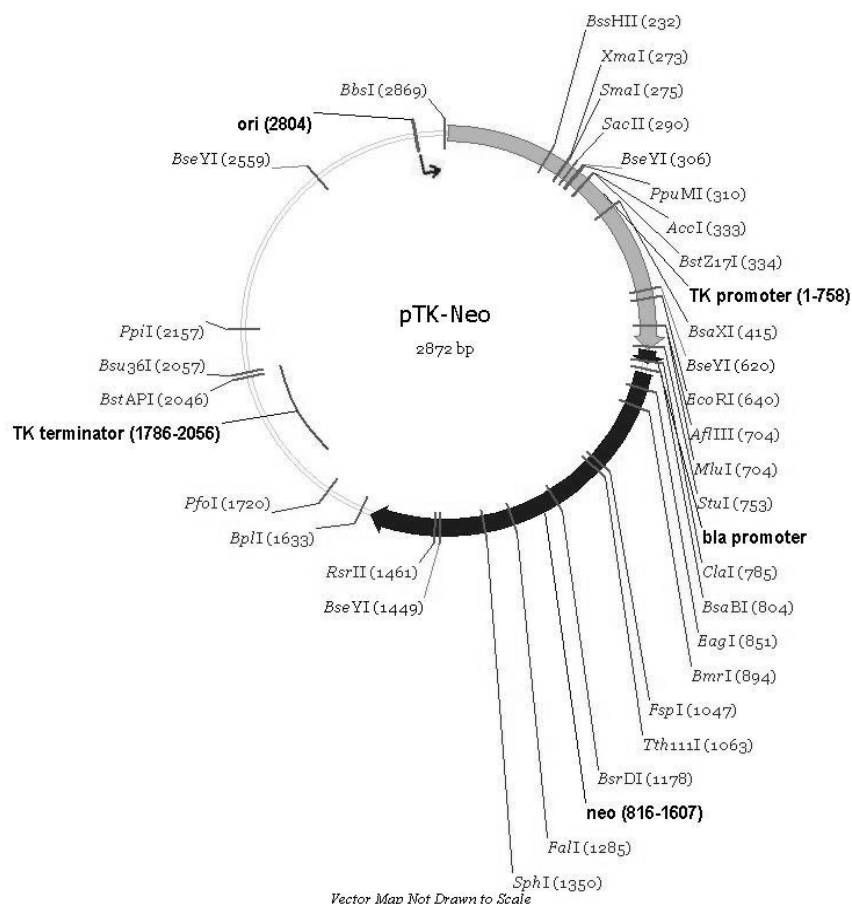
pTK-neo sequence landmarks

TK promoter	1-758
<i>bla</i> promoter	762-790
<i>neo</i> (Kan ^R) coding sequence	816-1607
TK terminator	1786-2056
pUC origin	2804

pTK-neo (Cat. No. 71284-3) can be used to select stably transformed mammalian cell lines using G418 (Cat. No. 345812). A minimal thymidine kinase (TK) promoter controls expression of the neomycin resistance gene. This promoter facilitates selection for stable integration of both the selection plasmid and a contranected expression plasmid into transcriptionally active sites in the genome. In bacteria, the beta-lactamase (*bla*) promoter drives expression of the *neo* gene, allowing growth in the presence of kanamycin.

Notes

All corresponding maps and tables are generated using VectorNTI® software. Restriction sites are numbered in accordance with the conventions of the respective software.



USA and Canada

Tel (800) 628-8470
bioscienceshelp@
emdchemicals.com

France

Freephone
0800 126 461

Germany

Freecall
0800 100 3496

Europe

Ireland

Toll Free
1800 409 445

United Kingdom

Freephone
0800 622 935

All other

European Countries
+44 115 943 0840

All Other Countries

Contact Your Local Distributor

www.merck4biosciences.com
bioscienceshelp@
emdchemicals.com

techservice@merckbio.eu

www.merck4biosciences.com

FOR RESEARCH USE ONLY. NOT FOR HUMAN OR DIAGNOSTIC USE.

Enzyme	#	Locations					Enzyme	#	Locations					Enzymes that do not cut pTK-neo:									
AccI	1	332					EcoO109I	2	309	2042				AatII	Acc65I	AccI	AfeI	AgeI					
AflII	2	26	150				EcoRI	1	639					AhdI	AleI	Alol	Apal	AscI					
AflIII	1	703					FalI	1	1284					AseI	AsiSI	BaeI	BamHI	BbvCI					
AlwNI	3	347	2045	2453			FspI	1	1046					BcgI	BclI	BfrBI	BglII	BlpI					
ApaLI	1	2548					HaeII	4	140	283	947	2622		BmgBI	BmtI	Bpu10I	BseRI	BsgI					
AvaI	3	98	219	272			KasI	2	279	943				BsiWI	BsmBI	BsmI	BspEI	BspLU11I					
AvrII	2	312	1781				MluI	1	703					BsrGI	BstEII	BstXI	BtrI	BtsI					
BanI	3	279	943	978			MscI	2	118	1026				DrallI	Ecl136II	EcoCRI	EcoNI	EcoRV					
BanII	3	303	1309	2019			MslI	3	532	652	1381			FseI	FspAI	HincII	HindIII	HpaI					
BbeI	2	283	947				NaeI	2	1446	1729				KpnI	MfeI	NcoI	NdeI	NheI					
BbsI	1	2868					NarI	2	280	944				NotI	NruI	NsiI	PacI	PciI					
BceAI	4	903	1370	1857	2362		NgoMIV	2	1444	1727				PflMI	PinAI	PmeI	PmlI	PshAI					
BciVI	3	776	1163	2664			NspI	3	589	1349	2866			Psil	PspOMI	Psrl	PvuI	SacI					
BglI	2	288	291				NspV	2	643	1626				Sall	SanDI	SbfI	Scal	SexAI					
Bme1580I	5	127	458	890	983	2552	PfoI	1	1719					Sfil	SgrAI	SnaBI	SpeI	SrfI					
BmrI	1	893					Ppil	1	2156					Sse8387I	Sspl	SwaI	TatI	XbaI					
BpII	1	1632					PpuMI	1	309					XcmI	XhoI	XmnI	ZraI						
Bpml	3	369	1727	1784			PstI	2	499	736													
BpuEI	4	1320	2238	2479	2777		PvuII	2	521	1050													
BsaAI	2	336	1248				RsrII	1	1460														
BsaBI	1	803					SacII	1	289														
BsaHI	4	280	944	1646	1725		SapI	3	127	1288	1498												
BsaI	2	376	1933				Sfcl	4	495	732	2406	2597											
BsaWI	5	378	975	1819	2509	2656	SfoI	2	281	945													
BsaXI	1	414					SmaI	1	274														
BseYI	4	305	619	1448	2558		SmlI	6	26	150	1335	2217	2494										
BsiEI	2	853	2528						2756														
BsiHKAI	3	1057	1247	2552			SphI	1	1349														
BsmAI	5	376	771	786	1805	1933	StuI	1	752														
BsmFI	9	339	408	505	549	659	StyI	2	312	1781													
		1095	1627	1736	1891		TaqII	4	244	981	1649	1893											
Bsp1286I	10	127	303	458	890	983	TspGWI	3	159	418	1819												
		1057	1247	1309	2019	2552	Tth111I	1	1062														
BspCNI	4	1599	2069	2171	2580		XmaI	1	272														
BspHI	2	773	2142																				
BspMI	3	831	1212	1662																			
BsrBI	3	771	1557	1611																			
BsrDI	1	1177																					
BsrFI	3	1263	1444	1727																			
BssHII	1	231																					
BssSI	3	1536	1662	2689																			
Bst1107I	1	333																					
BstAPI	1	2045																					
BstBI	2	643	1626																				
BstYI	7	1115	1361	1753	2112	2124																	
		2210	2221																				
BstZ17I	1	333																					
Bsu36I	1	2056																					
Btgl	5	119	170	286	320	533																	
Clal	1	784																					
DraI	2	2088	2107																				
DrdI	2	971	2760																				
EaeI	5	116	850	1024	1415	1442																	
EagI	1	850																					
EarI	3	127	1288	1498																			
Ecil	3	1473	2660	2806																			
Eco57I	4	489	1090	1522	2320																		
Eco57MI	7	369	489	1090	1522	1727																	
		1784	2320																				

© 2010 EMD Chemicals Inc., an Affiliate of Merck KGaA, Darmstadt, Germany. All rights reserved. The Novagen® name and Novagen® logo are registered trademarks of EMD Chemicals Inc. in the United States and in certain other jurisdictions. VectorNTI® is a registered trademark of Life Technologies Corporation in the US and/or certain other countries.

The terms and conditions governing EMD Chemicals' sale of this product are as indicated on our website (www.emdbiosciences.com).

USA and Canada
Tel (800) 628-8470
bioscienceshelp@emdchemicals.com

Germany
Tel 0800 100 3496
techservice@merckbiosciences.de

United Kingdom and Ireland
UK Freephone 0800 622935
Ireland Toll Free 1800 409445
customer.service@merckbiosciences.co.uk

All Other Countries
www.merck4biosciences.com
bioscienceshelp@emdchemicals.com