

pBACsurf-1 Transfer Plasmid

TB148 06/00

b	Locus	polh
	Promoter	polh
	N-terminal fusion	gp64 signal sequence
	C-terminal fusion option	gp64 mature domain
	Cloning options	polylinker

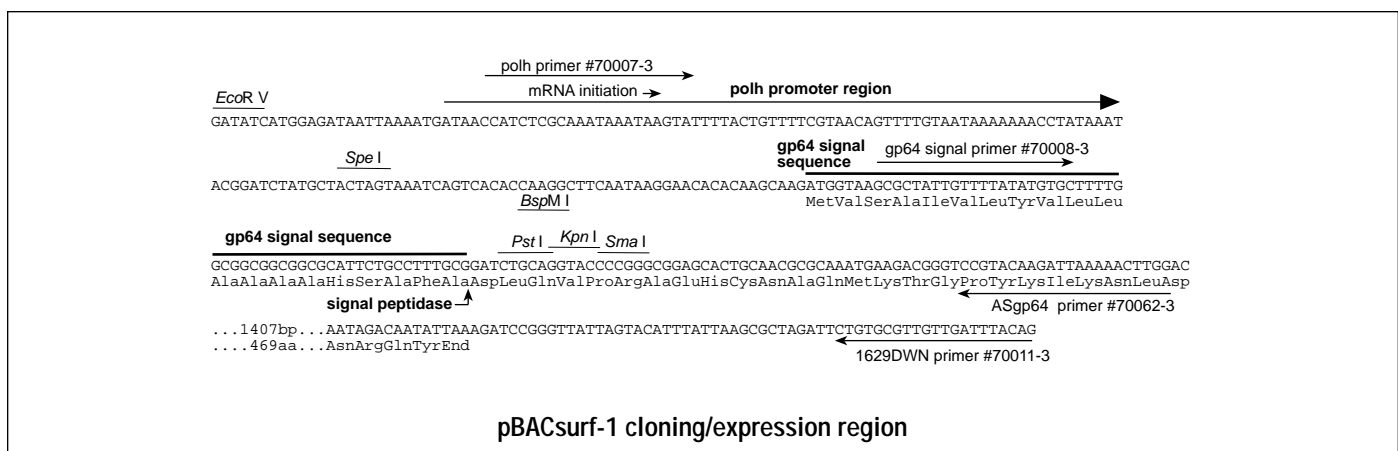
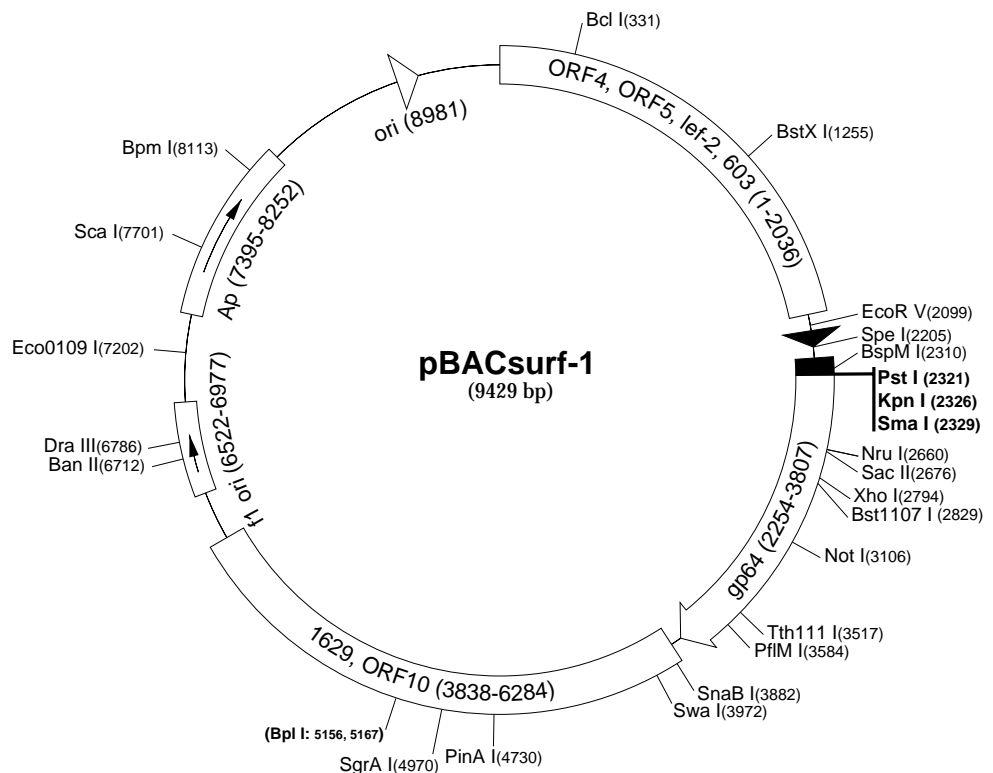
pBACsurf-1 is a baculovirus transfer plasmid (Cat. No. 70055-3) designed for expression of target proteins on the virion surface (1). For virus display, inserts are cloned in-frame between the upstream AcNPV gp64 signal sequence and downstream gp64 mature domain. The transmembrane region of the mature domain anchors the fusion protein in the membrane of the baculovirus nucleocapsid. By including a translation termination codon in the insert, target proteins can be secreted from the cell. The plasmid is compatible with BacVector™-1000, -2000 or -3000 Triple Cut Virus DNA for low background transfection and efficient utilization of the polh promoter. Unique restriction sites are indicated on the circle map. The cloning/expression region of the coding strand transcribed from the polh promoter is shown below. The f1 origin is oriented so that infection with helper phage will produce virions containing single stranded DNA that corresponds to the non-coding strand. Single stranded sequencing of phage-derived DNA can be performed using the gp64 signal primer (Cat. No. 70008-3).

Note: Before transfection of insect cells with pBACsurf-1 derived transfer vectors, it is highly recommended to linearize the vector with the enzyme BpI (Fermentas). This treatment prevents a frame-shift mutation from occurring within the ORF 1629 open reading frame.

pBACsurf-1 sequence landmarks

polh promoter region	2121-2191
polh transcription start	2142-2143
gp64 signal sequence	2254-2325
Multiple cloning sites (Pst I - Sma I)	2321-2329
gp64 mature domain	2332-3807

1. Boublik, Y., Di Bonito, P. and Jones, I.M. (1995) *BioTechnology* **13**, 1079-1084.



pBACsurf-1 Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
Enzyme	# Sites	Locations	Dsal	5	2673 2867 3290 3626 3680	Sfcl	6	2317 5215 6560 7936 8614		
AatII	3	3521 4163 7263	EaeI	11		SgrAI	1	8805 4970		
AccI	6	1047 1332 2828 2920 5250 6239	EagI	2	3106 6083	SmaI	1	2329		
Acil	92		EarI	5	1139 2609 6419 7383 9187	SnaBI	1	3882		
AflIII	15		Eco47III	4	2262 3840 5023 5839	SpeI	1	2205		
AhdI	3	721 1587 8182	Eco57I	3	3707 7516 8528	SphI	2	234 3244		
AluI	37		EcoO109I	1	7202	Sspl	11			
AlwI	19		EcoRII	8	2486 2801 2822 6360 8909 8922 9043 9331	StyI	5	2223 2399 3296 3626 3680		
Alw26I	14		EcoRV	1	2099	Swal	1	3972		
AlwNI	2	3584 8661	FauI	16		TaiI	44			
ApaLI	7	553 564 3384 5970 7013 7510 8756	Fnu4HI	60		TaqI	32			
ApoI	15		FokI	7	267 693 6380 7100 7743 8030 8211	TfiI	8	727 1534 2904 3845 5448 5461 9096 9236		
AvaI	5	365 2327 2794 2983 6890	FspI	7	184 1747 2732 3245 5402 6460 7959	ThaI	32			
Avall	4	733 2362 7818 8040	HaeII	14		TseI	27			
BamHI	2	2517 3530	HaeIII	24		Tsp45I	12			
BanI	12		HgaI	17		Tsp509I	71			
BanII	1	6712	HhaI	52		TspRI	20			
BbsI	4	1586 2362 4341 6115	HincII	8	14 942 1048 1333 3738 5251 6061 6191	Tth111I	1	3517		
BbvI	27		HindIII	2	4245 5291	VspI	7	960 1068 3754 4367 8007 9242 9301		
Bcgl	3	208 4697 7678	Hinfl	22		XhoI	1	2794		
Bcgl'	3	174 4663 7644	HpaI	4	942 3738 6061 6191	XmnI	3	4423 6007 7582		
BclI	1	331	HphI	17		Enzymes that do not cut pBACsurf-1:				
Bfal	12		KpnI	1	2326	AflIII	Apal	ApalBI	AscI	AvrII
BglI	3	3117 6470 8064	MaeIII	24		BaeI	BglIII	Bpu1102I	BssHIII	BstEII
BpII	2	5156 5167	MbolI	20		Bsu36I	EcoNI	EcoRI	FseI	NdeI
Bpml	1	8113	MluI	3	568 1487 2529	NheI	Pacl	Pmel	PshAI	Psp5II
Bpu10I	2	1504 2843	MnlI	38		RsrII	Sacl	SexAI	Sfil	SgfI
BsaI	4	2741 3047 3882 6783	MscI	3	3295 3454 5472	SrfI	Sse8387I	StuI	SunI	XbaI
BsaBI	4	512 2124 5402 5452	MseI	75		XcmI				
BsaHI	12		MslI	13						
BsaJI	13		MspI	25						
BsaWI	7	917 1921 4730 5999 7886 8717 8864	MspA1I	14						
BseRI	2	1620 3205	MunI	5	3506 3871 4216 4777 5142					
BsgI	4	3256 4527 4821 4975	MwoI	36						
BsiEI	9	718 2628 3109 6086 6441 7664 7813	NarI	4	3120 4713 4980 6480					
BsiHKAI	10	557 568 2339 3092 3388 5974 7017 7514 7599 8760	NciI	8	2328 2329 3816 7110 7145 7646 7997 8693					
BsII	18		NcoI	2	3626 3680					
BsmI	3	772 1025 2298	NgoAIV	3	1869 3396 6678					
BsmBI	6	2037 2434 2882 4732 7145 7187	NlaIII	30						
BsmFI	2	4288 5740	NlaIV	26						
Bsp1286I	13		NotI	1	3106					
BspEI	2	917 5999	NruI	1	2660					
BspLU11I	2	5257 9070	Nsil	6	772 1272 3073 3182 3242 3478					
BspMI	1	2310	NspI	8	210 234 937 1054 3244 5261 7157 9074					
BsrI	19		NspV	3	156 2776 5905					
BsrBI	8	47 793 4835 5130 6639 7340 9141 9382	PfIMI	1	3584					
BsrDI	5	250 1167 3478 7948 8122	PinAI	1	4730					
BsrFI	8	1869 3396 4730 4922 4970 6078 6678 8097	PleI	14						
BsrGI	8	95 277 930 1137 1856 2783 4212 5442	PmlI	2	2741 3047					
BssSI	5	579 6196 7206 7513 8897	Psp1406I	7	220 273 3128 6187 6996 7580 7953					
Bst1107I	1	2829	PstI	1	2321					
BstXI	1	1255	PvuI	2	6441 7813					
BstYI	11		Pvull	3	5828 6410 9250					
Cac8I	39		RcaI	4	3326 7237 7342 8350					
Clal	2	4854 5177	RsaI	31						
CviJI	104		SacII	1	2676					
Ddel	11		Sall	3	1046 1331 5249					
Dpnl	33		SapI	2	2609 9187					
Dral	9	985 1517 3972 5297 5306 6172 7604 8296 8315	Sau3AI	33						
DrallI	1	6786	Sau96I	10	733 2362 2670 6429 6789 7202 7818 8040 8057 8136					
DrdI	5	3283 3517 6830 7099 8968	Scal	1	7701					
			ScrFI	16						
			SfaNI	22						