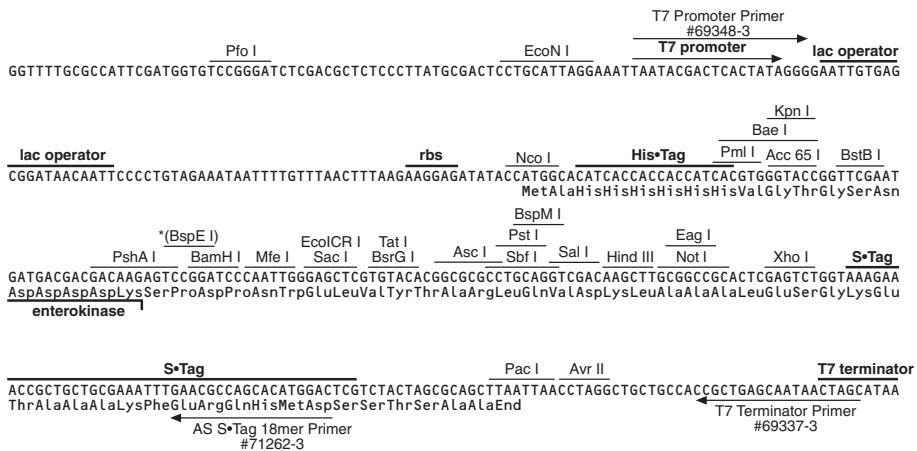
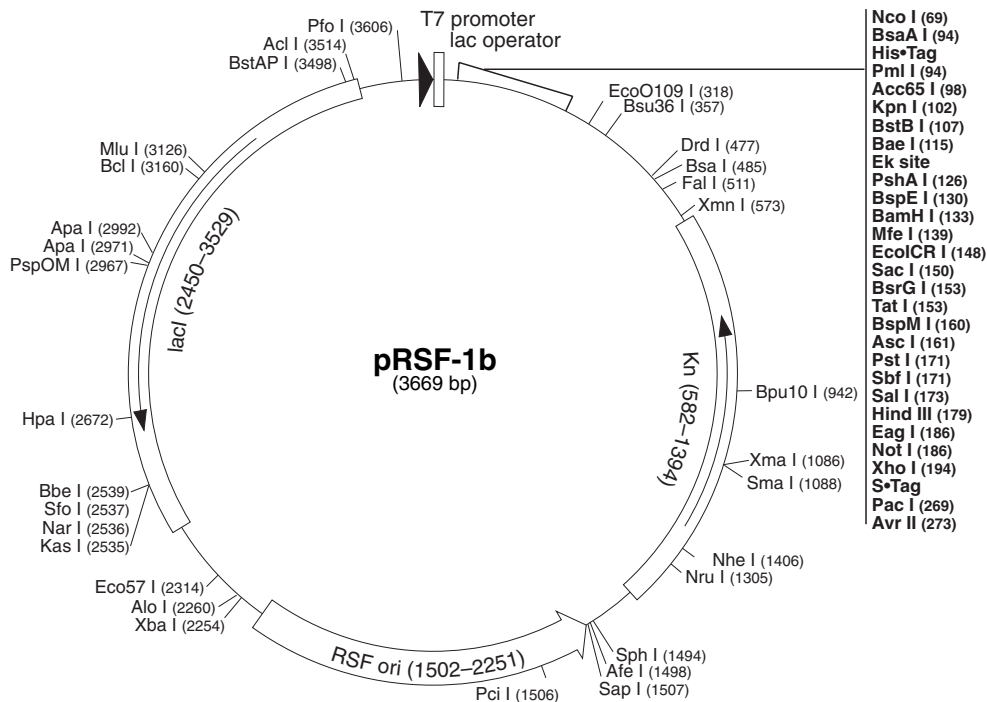


pRSF-1b Vector

	Cat. No.
pRSF-1b DNA	71363-3
pRSF-1b sequence landmarks	
T7 promoter	3653–3669
T7 transcription start	1
His•Tag [®] coding sequence	77–94
Multiple cloning sites (<i>Nco</i> I– <i>Avr</i> II)	69–278
S•Tag [™] coding sequence	206–250
T7 terminator	302–349
kan (Kn ^R) coding sequence	582–1394
RSF origin	1502–2251
<i>lac</i> I coding sequence	2450–3529

pRSF-1b carries a T7 promoter and *lac* operator to control transcription, a replication origin derived from RSF1030, and kanamycin antibiotic resistance (Kn^R). It also encodes an N-terminal His•Tag[®] sequence followed by an enterokinase (Ek) cleavage site and an optional C-terminal S•Tag[™] sequence. Unique sites are shown on the circle map. pRSF-1b is compatible with pET vectors (ColE1 origin), pCDF vectors (CloDF13 replication origin), and pACYC derived plasmids (P15A replication origin) carrying compatible antibiotic resistance markers. Sequencing can be performed using the T7 Promoter Primer (Cat. No. 69348-3) and AS S•Tag 18mer Primer (Cat. No. 71262-3) or T7 Terminator Primer (Cat. No. 69337-3).

Note: the *Bsp*E I site is modified by *dam* methylation, so the plasmid must be grown in *dam*⁻ hosts to use this site for cloning.



pRSF-1b cloning/expression region

pRSF-1b Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
Acc65I	1	98	Ecl136II	1	148		
AccI	2	174 251	Eco57I	1	2314		
AccII	1	3514	Eco57MI	3	2314 2853 3342		
AfeI	1	1498	EcoICRI	1	148		
AfilIII	2	1622 3174	EcoNI	2	1049 3642		
AgeI	2	101 406	EcoO109I	1	318		
Alol	1	2260	FalI	1	511		
Apal	1	2971	HaeII	5	1500 1859 2539 2782 3563		
ApaLI	2	1925 3194	HincII	3	175 2232 2672		
AscI	1	161	HindIII	1	179		
Asel	5	572 761 2432 2491 3652	HpaI	1	2672		
AvaI	2	194 1086	KasI	1	2535		
AvrII	1	273	KpnI	1	102		
BaeI	1	115	MfeI	1	139		
BamHI	1	133	MluI	1	3174		
BanI	4	98 2405 2535 3254	MspII	3	2808 2838 3126		
BanII	3	150 1311 2971	NarI	1	2536		
BbeI	1	2539	NcoI	1	69		
BbsI	2	2689 3028	NotI	1	186		
BceAI	4	174 641 2690 3317	NruI	1	1305		
BcgI	2	198 2854	NsiI	2	850 1116		
BciVI	4	568 1444 1814 2722	NspI	2	1494 1626		
BclI	1	3160	NspV	1	107		
BfrBI	2	848 1114	Pacl	1	269		
BipI	2	291 2149	PciI	1	1622		
Bme1580I	3	1929 2971 3198	PhiMI	4	94 241 702 3599		
BmrI	3	2376 3016 3253	PfoI	1	3606		
BpmI	2	2853 3342	PinAI	2	101 406		
Bpu10I	1	942	PmlI	1	94		
BpuEI	5	355 1702 2000 2180 2366	PshAI	1	126		
BsaAI	1	94	PspOMI	1	2967		
BsaHI	2	2536 3219	PstI	1	171		
BsaI	1	485	PvuII	2	2485 2578		
BsaWI	10	101 130 391 406 823 1644 1817 1964 2352 2855	SacI	1	150		
BsaXI	3	137 495 2506	SalI	1	173		
BseYI	3	1915 2640 2775	SapI	1	1506		
BsgI	2	3129 3329	SbfI	1	171		
BsiEI	6	189 476 964 1538 1951 2395	StcI	3	29 167 3665		
BsiHKAI	3	150 1929 3198	SfoI	1	2537		
BsmAI	7	485 942 1444 2559 2946 3072 3477	SmaI	1	1088		
BsmBI	2	942 2559	SmlI	6	194 334 1717 1979 2159 2381		
BsmI	2	1003 1080	SphI	1	1494		
Bsp1286I	5	150 1311 1929 2971 3198	Sse8387I	1	171		
BspCNI	6	283 370 934 1899 2162 2595	SspI	2	1037 1411		
BspEI	1	130	StyI	4	69 273 313 2127		
BspHI	2	565 1442	TaqII	4	710 1524 2208 2381		
BspLU11I	1	1622	TatI	1	153		
BspMI	1	160	TspGWI	3	1143 1155 2236		
BsrBI	4	13 563 1448 1555	Tth111I	2	477 2049		
BsrDI	2	2767 3133	XbaI	1	2254		
BsrFI	4	101 406 1004 3488	XcmI	3	2789 2807 3323		
BsrGI	1	153	XhoI	1	194		
BssHII	2	161 2763	XmaI	1	1086		
BssSI	2	149 1784	XmnI	1	573		
BstAPI	1	3498	Enzymes that do not cut pRSF-1b:				
BstBI	1	107	AarI	AatII	AfIII	AhdI	AleI
BstEII	1	2992	AlwNI	AsiSI	BbvCI	BglI	BglII
BstXI	3	3128 3251 3380	BmgBI	BmtI	BpI	BsaBI	BseRI
BstYI	4	133 709 2398 3610	BsiWI	BsmFI	Bst1107I	BstZ17I	BtrI
Bsu36I	1	357	DraI	DraIII	EcoRI	EcoRV	FseI
BtgI	1	69	FspAI	FspI	MscI	NaeI	NdeI
BtsI	5	383 1016 1103 2447 2815	NgoMIV	NheI	PmeI	Ppil	PpuMI
Clal	1	1269	Psil	Psrl	PvuI	RsrII	SacII
DrdI	1	477	SanDI	Scal	SexAI	SfiI	SgrAI
EaeI	2	186 2500	SnaBI	SpeI	SrfI	StuI	Swal
EagI	1	186	ZraI				
EarI	5	1146 1402 1506 2282 3557					
Ecil	3	1673 1819 3389					