

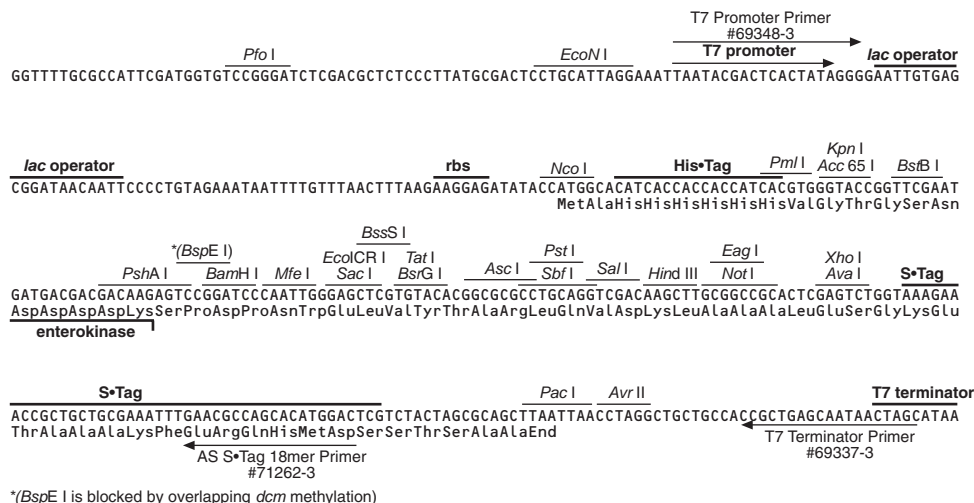
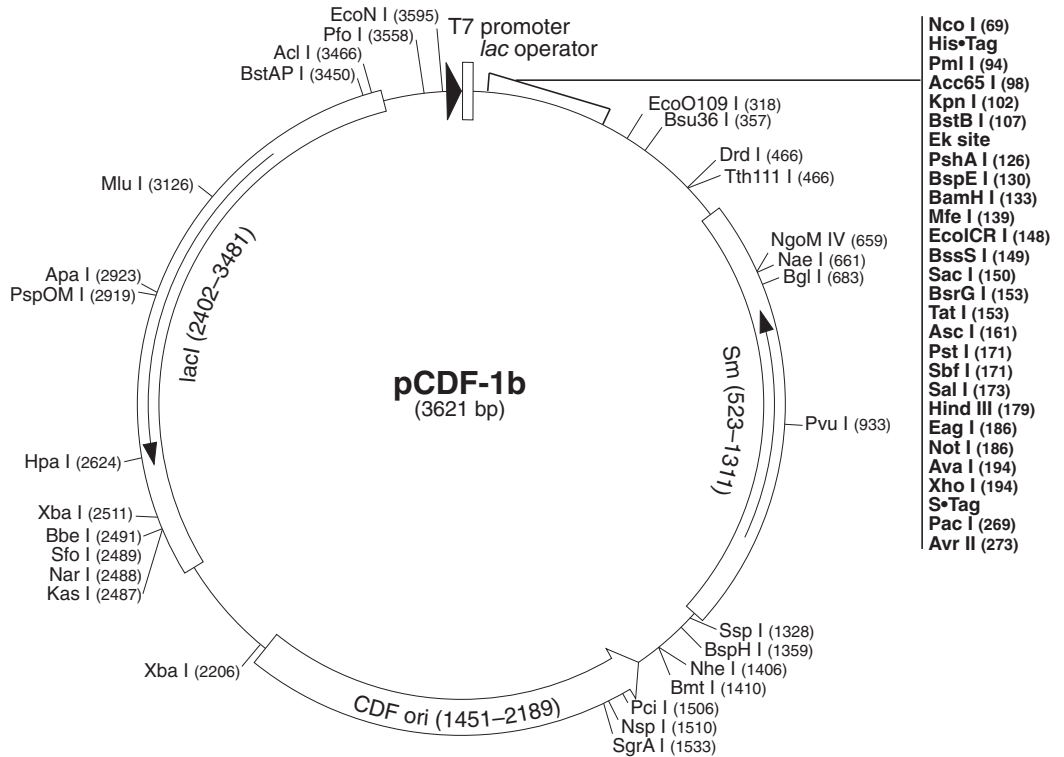
pCDF-1b Vector

TB392 0903

	Cat. No.
pCDF-1b DNA	71330-3
pCDF-1b sequence landmarks	
T7 promoter	3605-3621
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
CDF origin	1662-2411
aadA (Sm ^R) coding sequence	523-1311
lacI coding sequence	2402-3481

pCDF-1b carries a T7 promoter and *lac* operator to control transcription, a replication origin derived from CloDF13, and streptomycin/spectinomycin antibiotic resistance (Sm^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. pCDF-1b is compatible with pET vectors (ColE1 origin), pRSF vectors (RSF1030 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 *BspE* I site is modified by *dam* methylation, so the plasmid must be grown in *dam*⁻ hosts to use this site for cloning.



pCDF-1b cloning/expression region

pCDF-1b Restriction Sites

TB392 0903

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations
Acc65I	1	98	Eco57I	2	1168 2266
AccI	2	174 251	Eco57MI	5	1062 1168 2266 2805 3294
AcII	1	3466	EcoICRI	1	148
AflIII	2	1506 3126	EcoNI	1	3594
AgeI	2	101 406	EcoO109I	1	318
ApaI	1	2923	HaeII	8	670 819 1249 1454 1774
ApaLI	2	1033 3146			2491 2734 3515
AscI	1	161	HincII	2	175 2624
Asel	3	2384 2443 3604	HindIII	1	179
AvaI	1	194	HpaI	1	2624
AvrII	1	273	KasI	1	2487
BaeI	2	115 1812	KpnI	1	102
BamHI	1	133	MfeI	1	139
BanI	4	98 2357 2487 3206	MluI	1	3126
BanII	2	150 2923	MslI	4	1040 2760 2790 3078
BbeI	1	2491	NaeI	1	661
BbsI	2	2641 2980	NarI	1	2488
BceAI	7	174 1192 1236 1597 1708	NcoI	1	69
		2642 3269	NgoMIV	1	659
BcgI	2	198 2806	NheI	1	1406
BciVI	3	1361 1727 2674	NottI	1	186
BclI	2	1110 3112	NspI	1	1510
BglI	1	683	NspV	1	107
BlnI	2	291 1000	PacI	1	269
Bme1580I	3	1037 2923 3150	PciI	1	1506
Bmrl	7	664 673 769 1958 2328	PfIMI	3	94 241 3551
		2968 3205	PfoI	1	3558
BmtI	1	1410	PinAI	2	101 406
BpmI	3	1062 2805 3294	PmlI	1	94
BpuEI	4	355 1914 2155 2318	PshAI	1	126
BsaAI	2	94 555	PspOMI	1	2919
BsaHI	3	1256 2488 3171	PstI	1	171
BsaWI	10	101 130 391 406 842	PvuI	1	933
		1594 1730 1878 2304 2807	PvuII	2	2437 2530
BsaXI	2	137 2458	SacI	1	150
BseYI	3	1829 2592 2727	SalI	1	173
BsgI	2	3081 3281	SbfI	1	171
BsiEI	5	189 465 933 1420 2347	SfiI	4	29 167 1053 3617
BsiHKAII	4	150 1037 2049 3150	SfoI	1	2489
BsmAI	6	1361 1432 2511 2898 3024	SgrAI	1	1533
		3429	SmlI	5	194 334 1893 2170 2333
BsmBI	1	2511	Sse8387I	1	171
BsmFI	2	752 1513	Sspl	1	1328
Bsp1286I	5	150 1037 2049 2923 3150	StyI	5	69 273 313 534 876
BspCNI	3	283 370 2547	TaqII	2	1406 2333
BspEI	1	130	TatI	1	153
BspHI	1	1359	TspGWI	2	1032 1474
BspLU11I	1	1506	Tth111I	1	466
BspMI	2	160 961	XbaI	1	2206
BsrBI	3	13 1365 2204	XcmI	4	1483 2741 2759 3275
BsrDI	4	672 965 2719 3085	XhoI	1	194
BsrFI	6	101 406 659 712 1533			
		3440			
BsrGI	1	153			
BssHII	3	161 584 2715			
BssSI	1	149			
BstAPI	1	3450			
BstBI	1	107			
BstEII	2	1147 2944			
BstXI	3	3080 3203 3332			
BstYI	8	133 481 827 1667 2166			
		2177 2350 3562			
Bsu36I	1	357			
BtgI	4	69 1015 1485 2066			
BtsI	4	383 1197 2399 2767			
DrallI	2	558 1178			
DrdI	1	466			
EaeI	4	186 476 1221 2452			
EagI	1	186			
EarI	3	1319 2234 3509			
Ecil	4	850 1546 1733 3341			
Ecl136II	1	148			

Enzymes that do not cut pCDF-1b:

AarI	AatII	AfeI	AflIII	AhdI
AleI	Alol	AlwNI	AsiSI	BbvCI
BfrBI	BglIII	BmgBI	BpII	Bpu10I
BsaBI	Bsal	BseRI	BsiWI	BsmI
Bst1107I	BstZ17I	BtrI	Clal	Dral
EcoRI	EcoRV	FalI	FseI	FspAI
FspI	MscI	NdeI	NruI	NsiI
PmeI	Ppil	PpuMI	PsiI	PsiI
RsrII	SacII	SanDI	SapI	Scal
SexAI	SfiI	SmaI	SnaBI	SpeI
SphI	SrfI	StuI	Swal	XmaI
XmnI	ZraI			