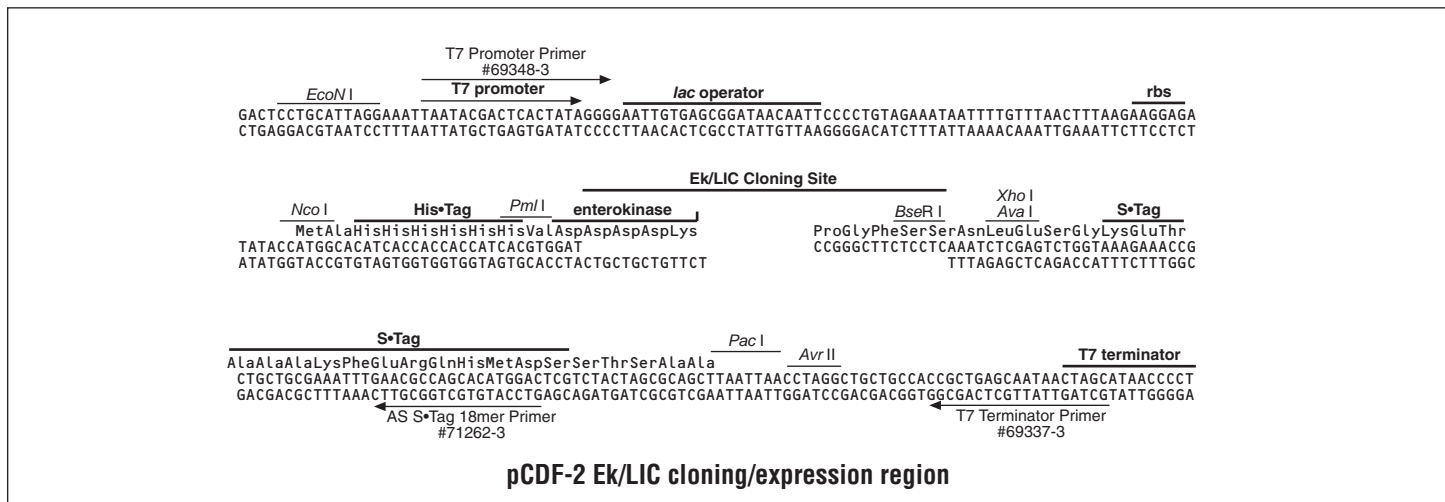
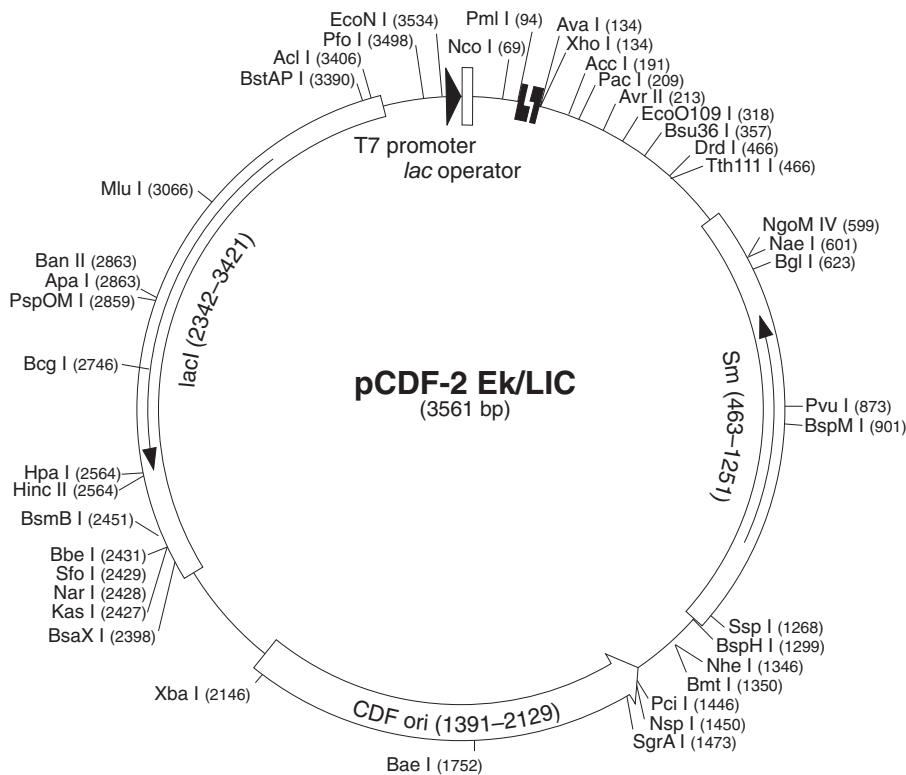


# pCDF-2 Ek/LIC Vector

TB394 0903

	Cat. No.
pCDF-2 Ek/LIC Cloning Kit	71337-3
<b>pCDF-2 Ek/LIC sequence landmarks</b>	
T7 promoter	3545-3561
T7 transcription start	1
His•Tag® coding sequence	77-94
Ek/LIC cloning site	101-129
S•Tag™ coding sequence	146-190
T7 terminator	242-289
<i>aadA</i> (SmR) coding sequence	463-1251
CDF origin	1391-2129
<i>lacI</i> coding sequence	2342-3421

The pCDF-2 Ek/LIC vector is prepared for rapid, directional cloning of PCR-amplified DNA for high-level expression of polypeptides. The vector carries a T7lac promoter to control transcription, a replication origin derived from CloDF13, and streptomycin/spectinomycin antibiotic resistance. Using specifically designed primers for amplification and the pCDF-2 Ek/LIC Cloning Kit (71337-3), inserts can be efficiently cloned without the need for restriction digestion or ligation. Fusion proteins contain an N-terminal cleavable His•Tag® sequence. Fusion to an optional C-terminal S•Tag™ sequence can also be created for detection, purification and quantification of fusion proteins. pCDF-2 is compatible with pET vectors (ColE1 origin), with pRSF vectors (RSF1030 replication origin), and pACYC plasmids (P15A replication origin) carrying compatible antibiotic resistance markers. Unique sites are shown on the circle map. Sequencing can be performed using the T7 Promoter Primer (Cat. No. 69348-3) and AS S•Tag 18 mer Primer (Cat. No. 71262-3) or T7 Terminator Primer (Cat. No. 69337-3).



## pCDF-2 Ek/LIC Restriction Sites

TB394 0903

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations
AccI	1	191	MslI	5	96 980 2700 2730 3018
AccII	1	3406	NaeI	1	601
AflIII	2	1446 3066	NarI	1	2428
AgeI	1	346	NcoI	1	69
ApaI	1	2863	NgoMIV	1	599
ApaLI	2	973 3086	NheI	1	1346
AseI	3	2324 2383 3544	NspI	1	1450
AvaI	1	134	PacI	1	209
AvrII	1	213	PciI	1	1446
BaeI	1	1752	PfIMI	3	94 181 3491
BanI	3	2297 2427 3146	PfoI	1	3498
BanII	1	2863	PmlI	1	94
BbeI	1	2431	PshAI	1	111
BbsI	2	2581 2920	PspOMI	1	2859
BceAI	6	1132 1176 1537 1648 2582	PvuI	1	873
		3209	PvuII	2	2377 2470
BcgI	1	2746	SfiI	3	29 993 3557
BciVI	3	1301 1667 2614	SfoI	1	2429
BclI	2	1050 3052	SgrAI	1	1473
BglI	1	623	SmlI	5	134 274 1833 2110 2273
BlpI	2	231 940	SspI	1	1268
Bme1580I	3	977 2863 3090	StyI	5	69 213 253 474 816
BmrI	7	604 613 709 1898 2268	TaqII	2	1346 2273
		2908 3145	TspGWI	2	972 1414
BmtI	1	1350	Tth111I	1	406
BpmI	3	1002 2745 3234	XbaI	1	2146
BpuEI	4	295 1854 2095 2258	XcmI	4	1423 2681 2699 3215
BsaAI	2	94 495	XhoI	1	134
BsaHI	3	1196 2428 3111			
BsaWI	8	331 346 782 1534 1670			
		1818 2244 2747			
BsaXI	1	2398			
BseRI	1	115			
BseYI	3	1769 2532 2667			
BsgI	2	3021 3221			
BsiEI	4	405 873 1360 2287			
BsiHKAI	3	977 1989 3090			
BsmAI	6	1301 1372 2451 2838 2964			
		3369			
BsmBI	1	2451			
BsmFI	2	692 1453			
Bsp1286I	4	977 1989 2863 3090			
BspCNI	3	223 310 2487			
BspHI	1	1299			
BspMI	1	901			
BsrBI	3	13 1305 2144			
BsrDI	4	612 905 2659 3025			
BsrFI	5	346 599 652 1473 3380			
BssHII	2	524 2655			
BstAPI	1	3390			
BstEII	2	1087 2884			
BstXI	3	3020 3143 3272			
BstYI	7	421 767 1607 2106 2117			
		2290 3502			
Bsu36I	1	297			
BtgI	4	69 955 1425 2006			
BtsI	4	323 1137 2339 2707			
DrallI	2	498 1118			
DrdI	1	406			
EaeI	3	416 1161 2392			
EarI	3	1259 2174 3449			
Ecil	4	790 1486 1673 3281			
Eco57I	2	1108 2206			
Eco57MI	5	1002 1108 2206 2745 3234			
EcoNI	1	3534			
EcoO109I	1	258			
HaeII	8	610 759 1189 1394 1714			
		2431 2674 3455			
HincII	1	2564			
HpaI	1	2564			
KasI	1	2427			
MluI	1	3066			

## Enzymes that do not cut pCDF-2 Ek/LIC:

AarI	AatII	Acc65I	AfeI	AflIII
AhdI	AleI	Alol	AlwNI	AscI
AsiSI	BamHI	BbvCI	BfrBI	BglII
BmgBI	BpII	Bpu10I	BsaBI	Bsal
BsiWI	BsmI	BspEI	BsrGI	BssSI
BstBI	BstZ17I	Clal	DraI	EagI
EcoICRI	EcoRI	EcoRV	FalI	FseI
FspAI	FspI	HindIII	KpnI	MfeI
MscI	NdeI	Nott	NruI	NsiI
PmeI	PpiI	PpuMI	PsiI	PsiI
PstI	RsrII	SacI	SacII	SalI
SanDI	SapI	SbfI	Scal	SexAI
SfiI	SmaI	SnaBI	SpeI	SphI
SrfI	StuI	Swal	TatI	XmaI
XmnI	ZraI			