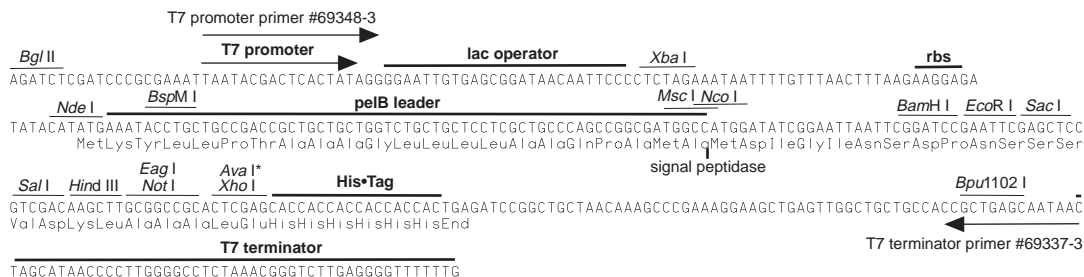
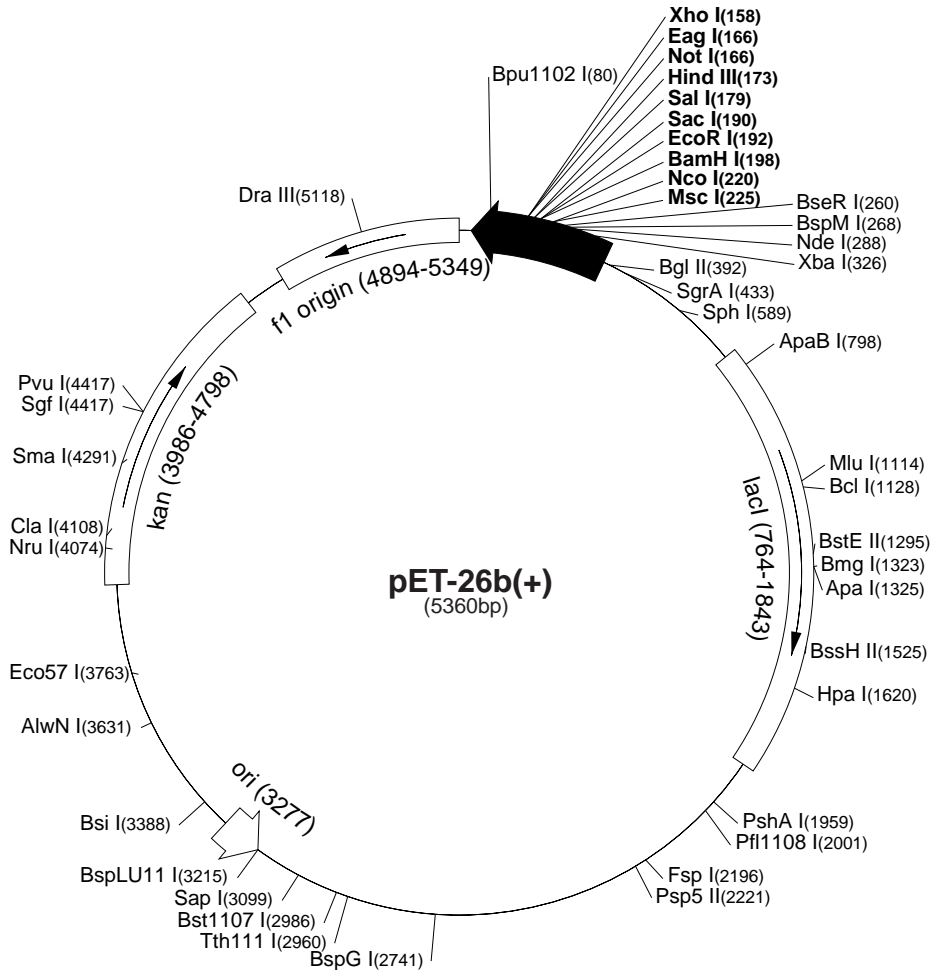


The pET-26b(+)⁺ vector (Cat. No. 69862-3) carries an N-terminal *pelB* signal sequence for potential periplasmic localization, plus optional C-terminal His•Tag[®] sequence. Unique sites are shown on the circle map. Note that the sequence is numbered by the pBR322 convention, so the T7 expression region is reversed on the circular map. The cloning/expression region of the coding strand transcribed by T7 RNA polymerase 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 coding strand. Therefore, single-stranded sequencing should be performed using the T7 terminator primer (Cat. No. 69337-3).

pET-26b(+)⁺ sequence landmarks

T7 promoter	361-377
T7 transcription start	360
<i>pelB</i> coding sequence	224-289
Multiple cloning sites	
(<i>Nco</i> I - <i>Xho</i> I)	158-225
His•Tag coding sequence	140-157
T7 terminator	26-72
<i>lacI</i> coding sequence	764-1843
pBR322 origin	3277
Kan coding sequence	3986-4798
f1 origin	4894-5349



pET-26b(+)⁺ cloning/expression region

pET-26b(+) Restriction Map

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AccI	2	180 2985	BsrFI	8	231 424 433 800 2012	NotI	1	166		
AccIII	7	881 1609 1940 2724 2865				NruI	1	4074		
		3167 4958	BssHII	1	1525	NsiI	2	4267 4533		
Acil	74		Bst1107I	1	2986	NspI	4	589 2560 2852 3219		
AfIII	2	1114 3215	BstEII	1	1295	Pfi1108I	1	2001		
AluI	22		BstXI	3	916 1045 1168	PfiMI	2	696 4680		
AlwI	13		BstYI	9	132 198 392 678 1890	PleI	9	375 663 750 1546 3109		
Alw21I	7	159 190 614 1098 2209						3594 4649 5053 5061		
		3033 3533	Cac8I	40		PshAI	1	1959		
Alw44I	3	1094 3029 3529	CjeI	24		Psp5II	1	2221		
AlwNI	1	3631	CjePI	18		Psp1406I	4	776 2144 2540 4903		
Apal	1	1325	Clal	1	4108	PvuI	1	4417		
ApaBI	1	798	CviJI	84		PvuII	3	1714 1807 2806		
ApoI	6	192 1389 4030 4214 4920	CviRI	22		RcaI	3	512 3935 4810		
		4931	DdeI	11		RsaI	3	1261 3021 4252		
AvaI	2	158 4289	DpnI	21		SacI	1	190		
AvaII	5	1666 2042 2130 2221 2500	DraIII	1	5118	SalI	1	179		
BamHI	1	198	DrdI	3	2908 3323 5073	SapI	1	3099		
BanI	8	436 457 571 1034 1753	DrdII	2	837 5123	Sau96I	14			
		1883 2009 5155	Dsal	3	220 551 2187	Sau3AI	21			
BanII	6	190 498 512 1325 4072	EaeI	5	166 223 422 554 1788	ScrFI	21			
		5193	EagI	1	166	SfaNI	23			
BbsI	4	1260 1599 1973 2333	EarI	3	732 3099 4230	SfiI	4	360 3480 3671 5337		
BbvI	28		Ecil	3	891 3289 3435	Sgfl	1	4417		
BccI	14		Eco47III	3	519 2020 2469	SgrAI	1	433		
Bce83I	6	21 1928 2098 3306 3604	Eco57I	1	3763	SmaI	1	4291		
		3845	EcoNI	2	649 4329	SphI	1	589		
Bcefl	6	633 974 1601 3717 4736	EcoO109I	3	53 547 2221	Sspl	2	4342 4910		
		5144	EcoRI	1	192	StyI	2	57 220		
Bcgl	8	160 194 1406 1440 1940	EcoRII	9	837 1152 1692 1749 3241	TaqI	15			
		1974 2792 2826				TaqII	6	1022 1240 1913 3117 4671		
BclI	1	1128	EcoRV	2	217 1564			5022		
Bfal	6	70 327 2229 3710 4017	FauI	17		TfiI	9	1793 2095 2265 2769 3190		
		5269	FokI	9	1160 1169 2434 2496 2574			4328 4384 4556 4647		
BglI	2	237 2178				Thal	35			
BglII	1	392	FspI	1	2196	Tsel	28			
BmgI	1	1323	GdiII	4	166 422 554 1788	Tsp45I	7	1295 2123 2654 2867 2962		
BpmI	4	952 1441 2075 2742	HaeI	7	225 842 2163 3230 3241			4564 5291		
Bpu10I	2	2321 4434				Tsp509I	22			
Bpu1102I	1	80	HaeII	14		Tth111I	1	2960		
BsaAI	2	2967 5118	HaeIII	24		Tth111II	8	953 1646 2676 3805 3812		
BsaBI	3	391 397 2412	Hgal	11				3844 4253 4380		
BsaHI	5	437 458 572 1071 1754	HgiEII	2	712 3801	UbaII	18			
BsaJI	10	57 220 551 557 1749	Hhal	46		VspI	6	207 375 1799 1858 4616		
		2187 3375 4288 4289 4690	Hin4I	3	1013 4103 4645			4805		
BsaWI	7	2 1433 1936 2404 3421	HincII	2	181 1620	XbaI	1	326		
		3568 4552	HindIII	1	173	XcmI	3	970 1486 1504		
BsaXI	2	1773 5066	Hinfl	18		XhoI	1	158		
Bsbl	2	2931 5025	HpaI	1	1620	XmnI	3	208 2773 4806		
BscGI	11		HphI	16						
BseRI	1	260	MaeII	14		Enzymes that do not cut pET-26b(+):				
BsgI	3	965 1165 2375	MaeIII	16		AatII	AfIII	AgeI	AscI	AvrII
Bsil	1	3388	MbolI	12		BaeI	BsaI	BsrGI	Bsu36I	DraI
BsiEI	6	169 271 1899 3131 3555	MluI	1	1114	Eam1105I	FseI	KpnI	MunI	NheI
		4417	MmeI	7	3430 3614 4059 4253 4615	NspV	Pacl	PmeI	PmlI	PstI
BsII	23					RleAI	RsrII	SacII	Scal	SexAI
Bsml	2	4301 4378	MnII	26		SfiI	SnaBI	SpeI	SrfI	Sse8387I
BsmAI	6	811 1216 1342 1729 2856	MscI	1	225	StuI	SunI	Swal		
		4433	MseI	26						
BsmBI	3	1729 2856 4433	MslI	6	1166 1454 1484 2202 2397					
BsmFI	4	575 2116 2486 5333								
BsoFI	46		MspI	30						
Bsp24I	10	404 436 955 987 1257	MspA1I	9	84 267 1144 1714 1807					
		1289 3708 3740 3886 3918								
Bsp1286I	12		MwoI	40						
BspEI	2	2 2404	NarI	4	437 458 572 1754					
BspGI	1	2741	NciI	12						
BspLU11I	1	3215	NcoI	1	220					
BspMI	1	268	NdeI	1	288					
BsrI	21		NgoAIV	5	231 424 2012 2172 5219					
BsrBI	4	347 3148 4816 5262	NlaIII	25						
BsrDI	2	1161 1527	NlaIV	21						