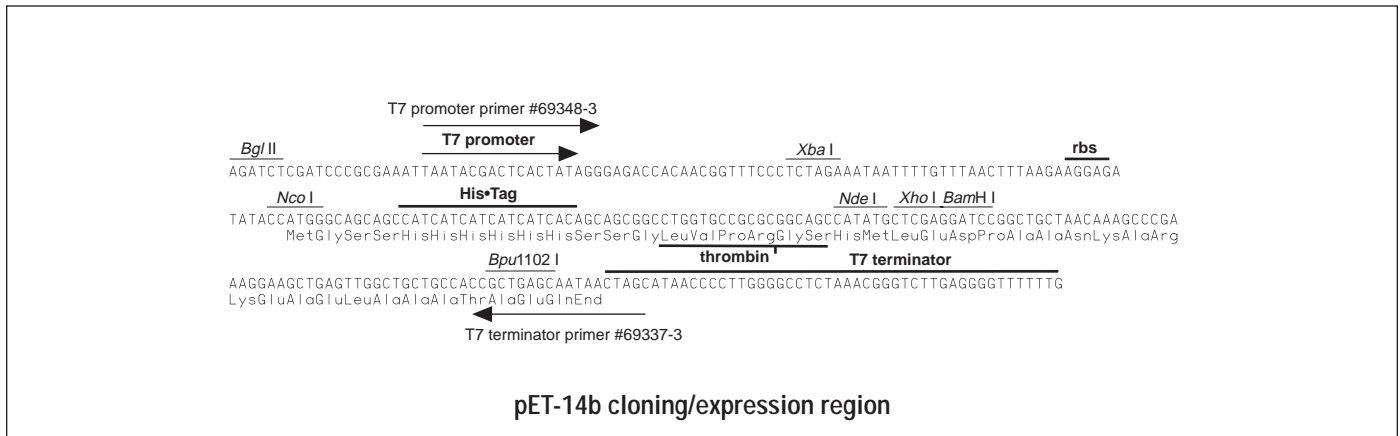
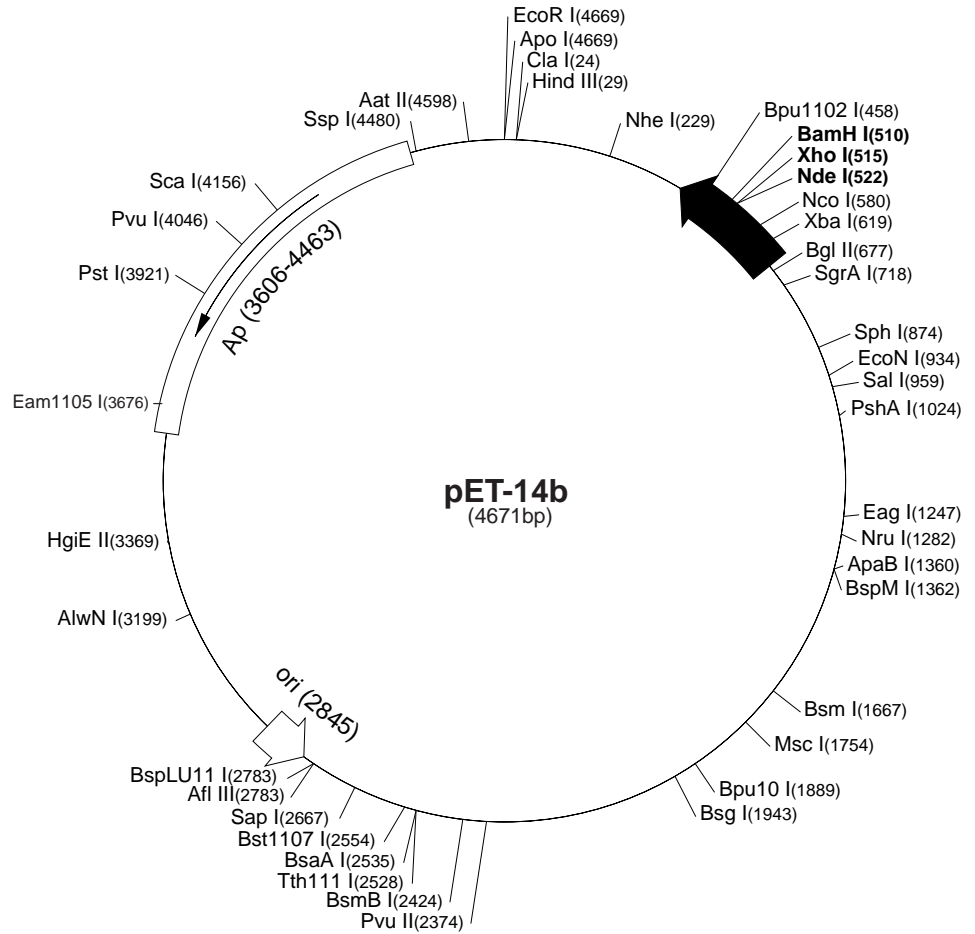


pET-14b Vector

The pET-14b vector (Cat. No. 69660-3) carries an N-terminal His•Tag® sequence followed by a thrombin site and three cloning sites. 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.

pET-14b sequence landmarks	
T7 promoter	646-662
T7 transcription start	645
His•Tag coding sequence	554-571
Multiple cloning sites (<i>Nde</i> I - <i>Bam</i> H I)	510-526
T7 terminator	404-450
pBR322 origin	2845
<i>bla</i> coding sequence	3606-4463



pET-14b Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AatII	1	4598	BstYI	9	510 677 1975 3424 3435 3521 3533 4301 4318	PshAI	1	1024		
AccI	2	960 2553	Cac8I	31		Psp5II	2	1747 1789		
AceIII	5	1005 2292 2433 2735 3975	CjeI	16		Psp1406I	4	1209 2108 3902 4275		
Acil	72		CjePI	22		PstI	1	3921		
AflIII	1	2783	Clal	1	24	PvuI	1	4046		
AluI	18		CviJI	81		PvuII	1	2374		
AlwI	14		CvIRI	21		RcaI	4	797 3503 4511 4616		
Alw21I	8	280 899 1486 1777 2601 3101 4262 4347	DdeI	10	458 479 1889 2051 2591 3058 3467 3633 4173 4599	RsaI	3	165 2589 4156		
Alw44I	3	2597 3097 4343	DpnI	25		Sall	1	959		
AlwNI	1	3199	DraI	3	3542 3561 4253	SapI	1	2667		
ApaBI	1	1360	DrdI	2	2476 2891	Sau96I	16			
ApoI	1	4669	Dsal	3	580 836 1755	Sau3AI	25			
AvaI	2	515 1733	EaeI	6	295 707 839 1247 1752 4064	Scal	1	4156		
Avall	8	1107 1195 1444 1747 1789 2068 3814 4036	EagI	1	1247	ScrFI	17			
BamHI	1	510	Eam1105I	1	3676	SfaNI	22			
BanI	10	76 119 537 721 742 856 1074 1513 1597 3624	EarI	2	2667 4471	Sfcl	5	138 645 3048 3239 3917		
BanII	2	783 797	Ecil	4	1703 2857 3003 3831	SgrAI	1	718		
BbsI	3	1038 1901 4654	Eco47III	4	234 804 1085 2037	SphI	1	874		
BbvI	28		Eco57I	2	3331 4343	Sspl	1	4480		
BccI	10	570 768 861 1298 1387 1694 1706 3713 3837 4124	EcoNI	1	934	StyI	3	435 580 1677		
Bce83I	7	399 993 1163 2874 3172 3413 4281	EcoO109I	5	431 832 1747 1789 4652	TaqI	10	24 339 516 674 682 960 1435 1576 2883 4327		
Bcefl	3	918 1475 3285	EcoRI	1	4669	TaqII	6	978 2685 4024 4209 4362 4379		
Bcgl	6	1005 1039 2360 2394 4181 4215	EcoRII	7	129 540 1366 1749 2809 2930 2943	TfiI	6	1160 1314 1612 1833 2337 2758		
Bfal	7	230 448 620 1797 3278 3531 3866	EcoRV	2	187 378	Thal	26			
BglI	3	1243 1477 3796	FauI	11		Tsel	28			
BglII	1	677	FokI	12		Tsp45I	9	124 212 1188 1455 2222 2435 2530 3932 4143		
Bpml	4	1140 1694 2310 3746	Fspl	4	262 1666 1764 3898	Tsp509I	10	58 251 611 661 1627 1641 3543 3849 4104 4669		
Bpu10I	1	1889	GdIII	5	295 707 839 1247 4064	Tth111I	1	2528		
Bpu1102I	1	458	HaeI	7	1228 1300 1357 1754 2798 2809 3261\	Tth111III	5	2244 3373 3380 3412 4668		
BsaI	2	644 3737	HaeII	11		UbaJI	24			
BsaAI	1	2535	HaeIII	24		VspI	2	660 3848		
BsaBI	3	676 682 1980	Hgal	11		XbaI	1	619		
BsaHI	6	722 743 857 1514 4213 4595	HgiEI	1	3369	XhoI	1	515		
BsaJI	10	115 129 435 580 836 842 1475 1677 1755 2943	HhaI	32		XmnI	2	2341 4275		
BsaWI	6	380 1001 1972 2989 3136 3967	Hin4I	5	16 334 1449 3675 3749	Enzymes that do not cut pET-14b:				
Bsbl	2	2499 4219	HincII	2	961 4217	AflII	Agel	Apal	AscI	AvrII
BscGI	13		HindIII	1	29	BaeI	BclI	BmgI	BsaXI	BseRI
BsgI	1	1943	Hinfl	11		BsrGI	BssHII	BstEII	BstXI	Bsu36I
Bsil	3	2956 4340 4647	HphI	12		DrallI	DrdII	FseI	HpaI	KpnI
BstEI	7	289 964 1250 2699 3123 4046 4195	MaeII	10	1209 1265 1854 1878 2108 2534 3486 3902 4275 4595	MluI	MunI	NotI	NsiI	NspV
BsII	21		MaeIII	17		Pacl	PmeI	PmlI	RleAI	RsrII
Bsml	1	1667	MbolI	11		Sacl	SacII	SexAI	SfiI	Sgfl
BsmAI	4	644 2424 3737 4513	Mmel	4	222 309 2998 3182	Smal	SnaBI	SpeI	SrfI	Sse8387I
BsmBI	1	2424	MnII	31		Stul	SunI	Swal	XcmI	
BsmFI	4	860 1181 1406 2054	MscI	1	1754					
BsoFI	52		MseI	18						
Bsp24I	8	689 721 3276 3308 3454 3486 4580 4612	Msil	7	1339 1770 1965 2356 3928 4087 4446					
Bsp1286I	10	280 783 797 899 1486 1777 2601 3101 4262 4347	MspI	28						
BspEI	2	380 1972	MspAII	8	462 548 1449 2374 2493 3125 3370 4311					
BspGI	3	1367 1444 2309	MwoI	36						
BspLU11I	1	2783	NarI	4	722 743 857 1514					
BspMI	1	1362	NciI	10	171 843 1567 1793 2121 2427 2462 3163 3859 4210					
BsrI	19		NcoI	1	580					
BsrBI	2	2716 4517	NdeI	1	522					
BsrDI	2	3737 3911	NgoAIV	4	709 1077 1237 1591					
BsrFI	7	160 709 718 1077 1237 1591 3756	NheI	1	229					
Bst1107I	1	2554	NlaIII	27						
			NlaIV	26						
			NruI	1	1282					
			NspI	4	874 2128 2420 2787					
			Pfi1108I	2	1066 3694					
			PfiMI	2	1629 1678					
			PleI	5	660 948 2677 3162 3665					