

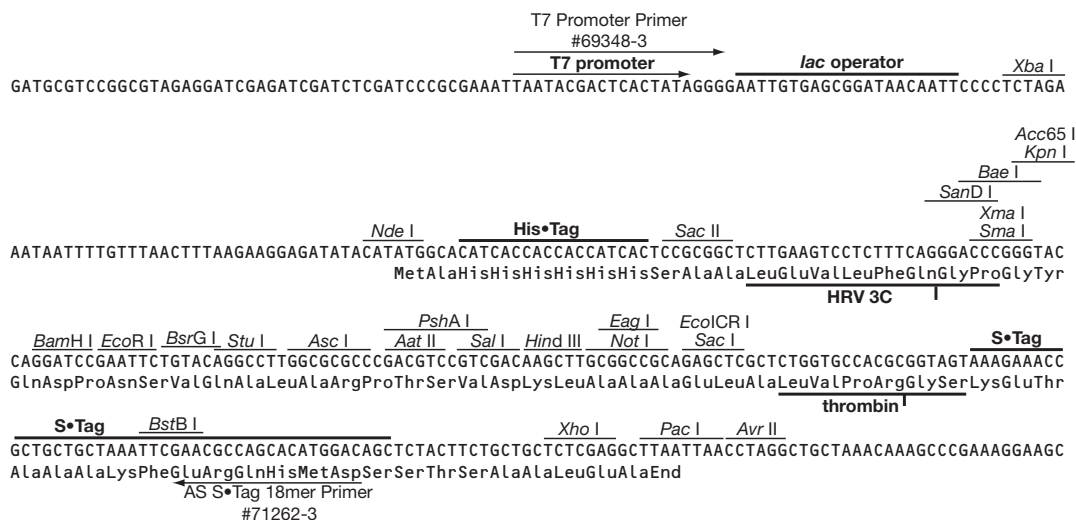
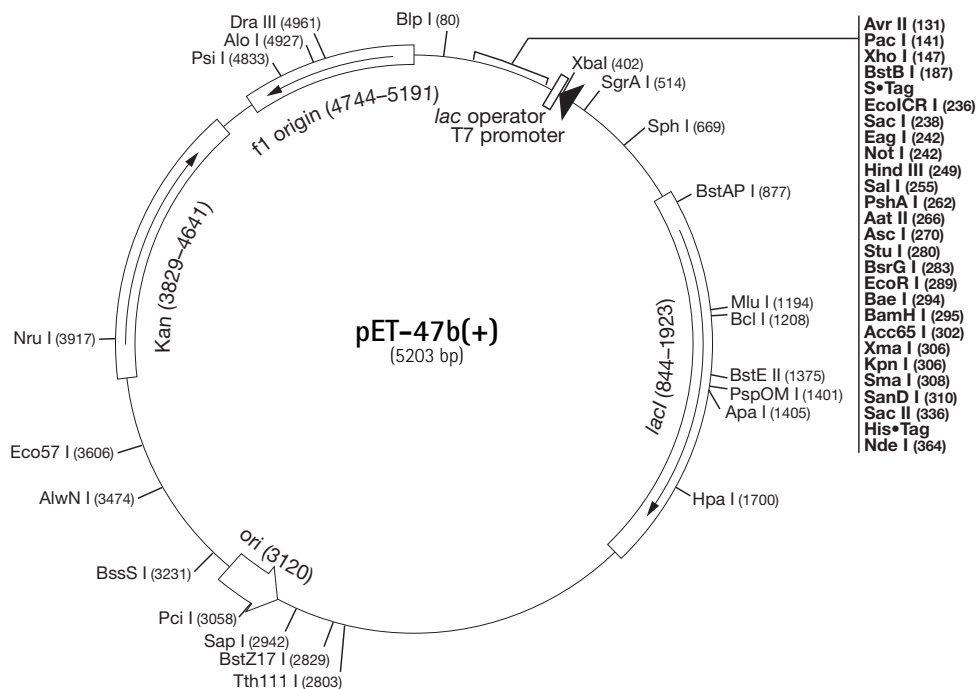
pET-47b(+)⁺ Vector

TB415 0804

	Cat. No.
pET-47b(+) ⁺ DNA	71461-3
pET-47b(+)⁺ sequence landmarks	
T7 promoter	437-453
T7 transcription start	436
His•Tag coding sequence	341-359
Multiple cloning sites (<i>SanD</i> I – <i>Avr</i> II)	131-315
S•Tag coding sequence	168-212
T7 terminator	26-73
<i>lacI</i> coding sequence	844-1923
pBR322 ori	3120
Kan coding sequence	3829-4641
f1 origin	4744-5191

The pET-47b(+)⁺ vector carries an N-terminal His•Tag[®] coding sequence followed by a recognition site for the human rhinovirus (HRV) 3C protease. This protease is highly specific for cleavage of the sequence LEVLFG↓GP (1), and is active at low temperatures (2). pET-47b(+)⁺ also contains an optional C-terminal thrombin recognition site followed by an S•Tag[™] coding sequence. Unique restriction 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 circle 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 the helper phage will produce virions containing single-stranded DNA that corresponds to the coding strand. Therefore, single-stranded sequencing should be performed using the AS S•Tag 18mer Primer (Cat. No. 71262-3).

1. Cordingley, M.G., Register, R.B., Callahan, P.L., Garsky, V.M., and Colonno, R.J. (1989) *J. Virol.* 63, 5037-5045.
2. Wang, Q.M., Johnson, R.B., Cox, G.A., Villarreal, E.C., and Loncharich, R.J. (1997) *Anal. Biochem.* 252, 238-245.



pET-47b(+)⁺ cloning/expression region

pET-47b(+) Restriction Sites

TB415 0804

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzymes that do not cut pET-47b(+):					
AatII	1	266	BtsI	4	1559 1927 4121 4208	AarI	AfII	AgeI	AhdI	AleI	AsiI
Acc65I	1	302	ClaI	2	471 3951	BbvCI	BglI	BglII	BmgBI	BmtI	BpII
AccI	2	256 2828	DraIII	1	4961	BsaI	BseRI	BsiWI	BspMI	Bsu36I	BtrI
AccII	3	856 2383 4746	DrdI	3	2751 3166 4916	DraI	EcoRV	FalI	FseI	FspAI	FspI
AfeI	2	599 2312	EaeI	4	242 502 634 1868	MfeI	MscI	NcoI	NheI	PinAI	PmeI
AflIII	2	1194 3058	EagI	1	242	PmlI	Psrl	PstI	PvuI	RsrII	SbfI
Alol	1	4927	EarI	3	812 2942 4073	Scal	SexAI	SfiI	SnaBI	SpeI	SrfI
AlwNI	1	3474	Ecil	3	985 3120 3266	Sse8387I	Swal				
ApaI	1	1405	Eco57I	1	3606						
ApaLI	3	1174 2872 3372	Eco57MI	4	1032 1521 2585 3606						
AscI	1	270	EcoCRI	1	236						
Asel	5	451 1879 1938 4459 4648	EcoNI	3	318 729 4172						
AvaI	2	147 306	EcoO109I	4	53 310 627 2064						
AvrII	1	131	EcoRI	1	289						
BaeI	1	294	HaeII	13							
BamHI	1	295	HincII	2	257 1700						
BanI	9	223 302 516 537 651	HindIII	1	249						
		1114 1833 1963 4998	HpaI	1	1700						
BanII	6	238 578 592 1405 3915	KasI	4	516 537 651 1833						
		5036	KpnI	1	306						
BbeI	4	520 541 655 1837	MluI	1	1194						
BbsI	3	1340 1679 2176	MslI	6	1246 1534 1564 2045 2240						
BceAI	6	713 1053 1680 3560 4579			2631						
		4986	NaeI	2	506 5064						
BcgI	3	236 1520 2635	NarI	4	517 538 652 1834						
BciVI	3	1651 3261 4655	NdeI	1	364						
BclI	1	1208	NgoMIV	2	504 5062						
BfrBI	2	4108 4374	NotI	1	242						
BlpI	1	80	NruI	1	3917						
Bme1580I	4	1178 1405 2876 3376	NsiI	2	4110 4376						
BmrI	5	723 1120 1357 1997 2797	NspI	4	669 2403 2695 3062						
BpmI	3	1032 1521 2585	NspV	1	187						
Bpu10I	2	2164 4277	Pacl	1	141						
BpuEI	5	21 2008 3149 3447 3688	PciI	1	3058						
BsaAI	2	2810 4961	PfiMI	3	179 776 4523						
BsaBI	3	467 477 2255	PfoI	2	761 2700						
BsaHI	6	263 517 538 652 1151	PpiI	2	3773 4927						
		1834	PpuMI	2	310 2064						
BsaWI	7	2 1513 2016 2247 3264	PshAI	1	262						
		3411 4395	PsiI	1	4833						
BsaXI	3	331 1869 4925	PspOMI	1	1401						
BseYI	3	1593 1728 3362	PvuII	3	1794 1887 2649						
BsgI	3	1045 1245 2218	SacI	1	238						
BsiEI	4	245 1979 2974 3398	SacII	1	338						
BsiHKAI	6	238 694 1178 2052 2876	Sall	1	255						
		3376	SanDI	1	310						
BsmAI	6	891 1296 1422 1809 2699	SapI	1	2942						
		4276	Sfcl	4	436 3323 3514 5180						
BsmBI	3	1809 2699 4276	SfoI	4	518 539 653 1835						
BsmFI	4	296 655 2329 5176	SgrAI	1	513						
BsmI	2	4144 4221	Smal	1	308						
Bsp1286I	11		SmII	6	36 147 1987 3164 3426						
BspCNI	9	93 114 1779 2156 2318			3703						
		2858 3346 3755 4290	SphI	1	669						
BspEI	2	2 2247	SspI	2	4185 4753						
BspHI	3	592 3778 4653	StuI	1	280						
BspLU11I	1	3058	StyI	3	57 131 275						
BsrBI	4	423 2991 4659 5105	TaqII	4	1993 2960 4514 4865						
BsrDI	2	1241 1607	TatI	2	283 2862						
BsrFI	5	504 513 880 4214 5062	TspGWI	5	274 2183 2501 4069 4081						
BsrGI	1	283	Tth111I	1	2803						
BssHII	2	270 1605	XbaI	1	402						
BssSI	1	3231	XcmI	3	1050 1566 1584						
Bst1107I	1	2829	XhoI	1	147						
BstAPI	1	877	XmaI	1	306						
BstBI	1	187	XmnI	2	2616 4649						
BstEII	1	1375	ZraI	1	264						
BstXI	3	996 1125 1248									
BstYI	7	295 758 1970 2250 3699									
		3710 4509									
BstZ17I	1	2829									
BtgI	2	335 631									