

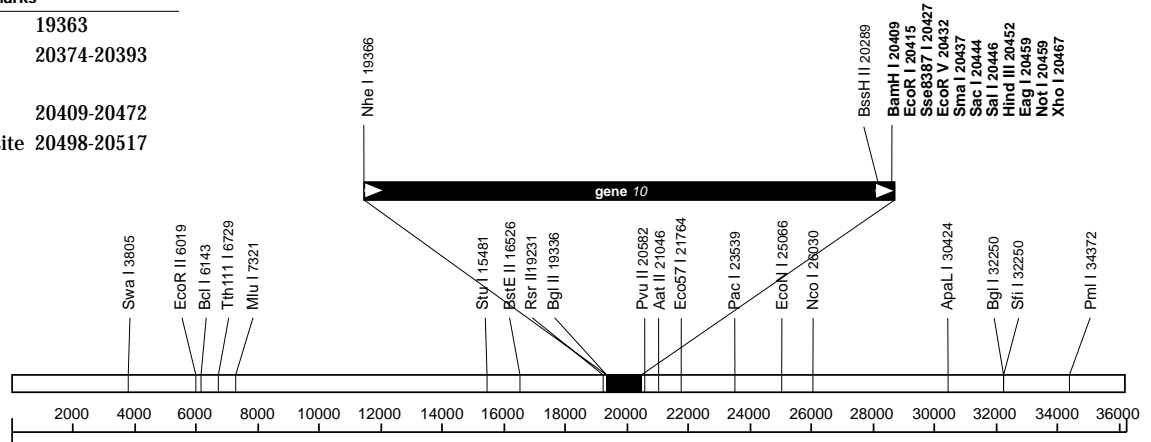
T7Select® 10-3b Vector

The T7Select® 10-3b vector is a protein display vector based on bacteriophage T7. The vector displays an average of 5–15 copies of peptides and proteins up to 1200 aa in size on the surface of the T7 capsid. Target sequences are fused to the C-terminus of the 10B capsid protein near amino acid 348. T7Select10 phage are grown on a complementing host (BLT5403) that supplies large amounts of the 10A capsid protein from a plasmid. Capsids thus contain mostly 10A protein, along with 5–15 copies of 10B fusion protein per virion.

The linear genome of T7Select10-3b is 36,249 bp in size. Unique restriction sites are shown on the maps below (cloning sites are shown in **bold** type on the line map).

T7Select10-3b sequence landmarks

10B translation start	19363
T7SelectUP priming site	20374-20393
Multiple cloning region (<i>Bam</i> H I - <i>Xho</i> I)	20409-20472
T7SelectDOWN priming site	20498-20517

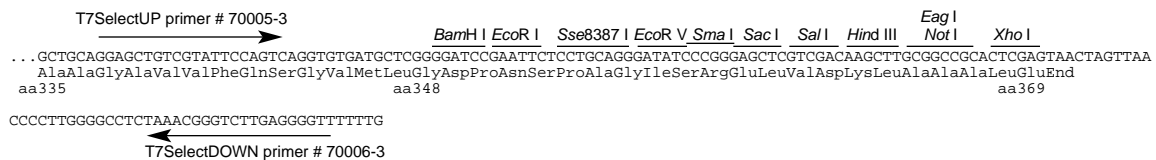


T7Select10-3b (36,249bp)

gene 10B @	insert	
left arm...GATCCG	AATTXXXX(N)XXXX	AGCTT...right arm
left arm...CTAGGCTTAA	XXXX(N)XXXXTCGA	A...right arm
...AspPro	AsnPhe...	
	AsnLeu...	
	AsnSer...	
	AsnTyr...	
	AsnCys...	
	AsnTrp...	

Reading frame of inserts cloned into *Eco*R I/*Hind* III T7Select10-3b vector arms.

The inserts require a 5'-AATT "sticky end" on the top strand (amino terminal side) and a 5'-AGCT "sticky end" on the bottom strand, either created with oligonucleotides or by *Eco*R I/*Hind* III digestion. The reading frame requires the AAT (Asn) initial codon, followed by a TXX codon (possible second amino acids are shown). Use of *Eco*R I cleavage products places C in the second position, resulting in a TCX codon (Ser).



T7Select10-3b cloning region

T7Select® 10-3b Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AatII	1	21046	DrarI	9	276 441 3805 4333 8555 13331 16758 27683 36060	RsrII	1	19231		
AccI	31		DrarII	15		SacI	1	20444		
Acil	176		DrdI	10	538 4041 4796 13663 14048 14515 28472 29322 31059 31233	Sall	1	20446		
AfIII	16		Dsal	25		SanDI	3	11865 22707 24905		
AfIII	21		Eael	3	20459 31850 32251	SapI	4	7063 11249 16607 33367		
AhdI	11		Eagl	1	20459	Sau3AI	8	6143 6246 8994 10814 19336 20409 31996 32399		
AluI	125		EarI	44		Sau96I	78			
AlwI	3	20404 20417 32407	Eco57I	1	21764	Scal	4	11977 30414 35164 35306		
Alw26I	89		EcoNI	1	25066	ScrFI	10	537 6021 11159 12134 18300 19589 20436 20437 24674 34034		
AlwNI	16		EcoO109I	22		SfaNI	83			
ApaLI	1	30424	EcoRI	1	20415	Sfcl	45			
ApoI	13		EcoRII	1	6019	Sfil	1	32250		
AvaI	7	8344 12362 20404 20435 20467 22431 26394	EcoRV	1	20432	SmaI	1	20437		
Avall	54		FauI	21		SnaBI	13			
AvrII	3	16464 27875 29534	Fnu4HI	140		SpeI	3	5526 13739 20475		
BamHI	1	20409	FokI	87		Sse8387I	1	20427		
BanI	32		Fspl	6	1970 12125 21766 25178 26116 31627	Sspl	6	7446 23544 28843 31802 32267 32546		
BanII	2	16066 20444	HaeII	25		SitI	1	15481		
BbsI	35		aeIII	62		StyI	35			
BbvI	103		HgaI	60		Swal	1	3805		
Bcgl	19		HhaI	92		TalI	156			
Bcgl'	19		HincII	57		TaqI	97			
BclI	1	6143	HindIII	1	20452	TfiI	97			
Bfal	51		Hinfl	200		Thal	60			
BglI	1	32250	HpaI	18		Tsel	103			
BglII	1	19336	HphI	93		Tsp45I	99			
Bpml	21		KpnI	5	41 3449 7024 20141 36130	Tsp509I	71			
Bpu10I	38		MaeIII	196		TspRI	82			
Bpu1102I	19		MbolI	131		Tth111I	1	6729		
BsaI	27		MluI	1	7321	UbaEI	5	1780 6602 7218 7787 14195		
BsaAI	35		MnlI	309		VspI	10	387 435 8326 14987 15849 21133 23568 27154 31072 35523		
BsaBI	7	870 1015 4982 8993 15364 16449 33934	MscI	2	31852 32253	XbaI	2	10309 30606		
BsaHI	7	339 8773 8916 20603 21043 21968 31817	MseI	187		XcmI	7	13850 17255 17417 25607 25682 26028 26649		
BsaJI	82		MslI	38		XhoI	1	20467		
BsaWI	32		MspI	56		XmnI	12			
BseRI	10	10085 10097 13966 16422 18878 27318 27747 35396 35560 35706	MspA1I	28		Enzymes that do not cut T7Select10-3b:				
BsgI	18		MunI	8	3571 11038 14199 17888 17994 30474 31759 32699	Apal	Ascl	BspEI	Eco47III	FseI
BsiEI	17		MwoI	147		NgoAIV	PvuI	SacII	SexAI	Sgfl
BsiHKA1	26		NarI	2	8916 31817	SgrAI	SphI	SrfI	SunI	
BsII	86		NciI	9	537 11159 12134 18300 19589 20436 20437 24674 34034					
BsmI	13		NcoI	1	26030					
BsmBI	14		NdeI	6	3507 4195 10767 16415 24038 30043					
BsmFI	43		NheI	1	19366					
Bsp1286I	40		NlaIII	127						
BspLU11I	6	799 11802 15602 15723 15769 33883	NlaIV	95						
BspMI	17		NotI	1	20459					
BsrI	111		NruI	3	1103 23878 25031					
BsrBI	14		Nsil	7	6585 16450 18439 26313 34637 35450 35609					
BsrDI	15		NspI	21						
BsrFI	3	19789 32022 33777	NspV	5	3849 9639 15675 19322 27749					
BsrGI	10	1949 3348 8546 17337 21651 22727 23153 28624 32174 34814	Pacl	1	23539					
BssHII	1	20289	PfIMI	7	7148 15226 22169 31137 32341 32704 34346					
BssSI	30		PinAI	2	32022 33777					
Bst1107I	8	5511 15973 19727 20776 23451 24803 25083 35941	PleI	103						
BstEII	1	16526	PmeI	2	276 8555					
BstXI	9	1702 7465 7937 11857 12196 24271 25755 30975 35277	PmlI	1	34372					
BstYI	3	19336 20409 32399	PshAI	4	12529 25316 26045 34930					
Bsu36I	28		Psp1406I	17						
Cac8I	97		Psp5II	12						
Clal	3	19317 23412 33392	PstI	2	20373 20427					
CviJI	515		PvuII	1	20585					
Ddel	264		RcaI	10	2903 4800 8906 12941 14716 17049 24161 25220 30745 35217					
Dpnl	8	6145 6248 8996 10816 19338 20411 31998 32401	RsaI	157						