

# pT7Blue-2 Vector

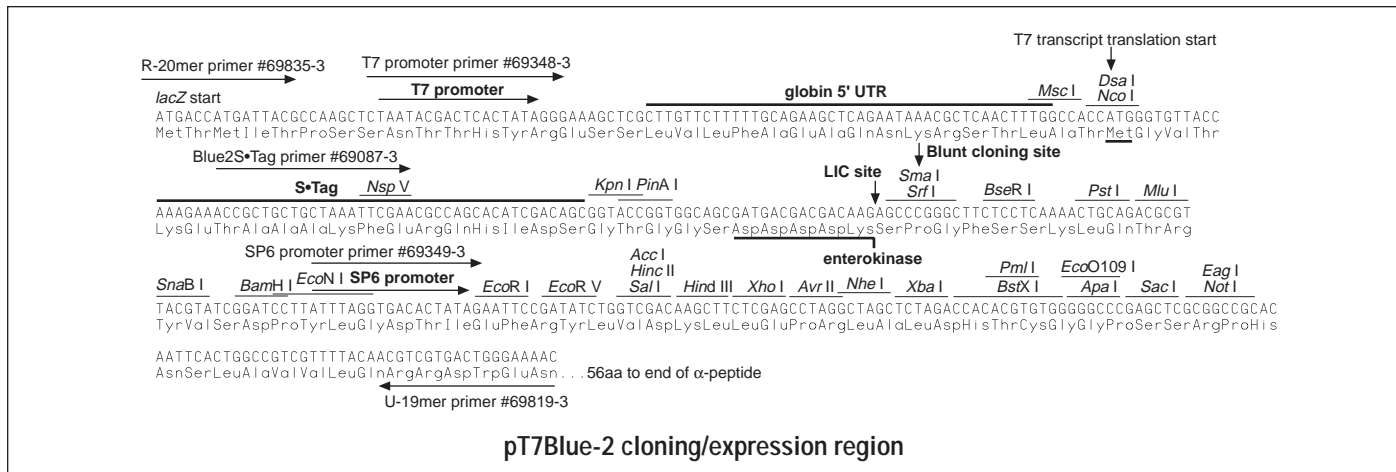
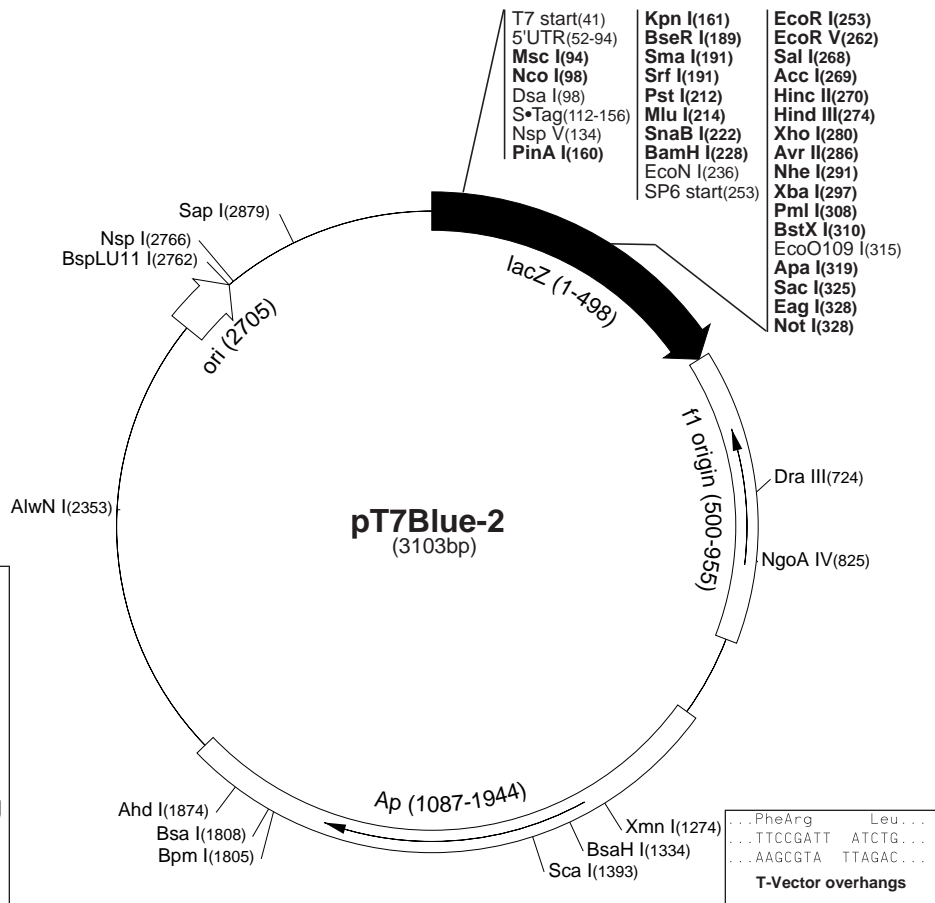
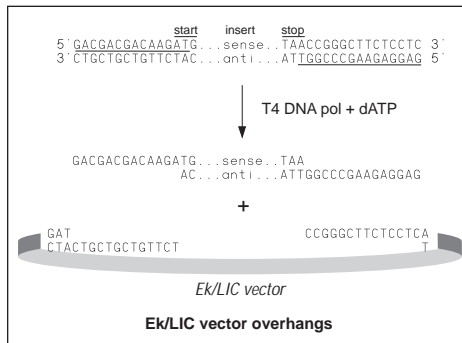
pT7Blue-2 contains the pUC19 backbone (including high-copy number origin of replication and *lacZ* sequences), f1 origin of replication and modified multiple cloning region designed for optimal expression of cloned inserts by *in vitro* translation. The multiple cloning region contains a T7 promoter upstream of the *Xenopus*  $\beta$ -globin 5' UTR translation enhancer and S•Tag™ sequences, followed by an enterokinase cleavage site and extensive region of cloning sites. An SP6 promoter is placed within the multiple cloning sites to allow the synthesis of essentially unfused RNA from DNA inserted downstream. Unique sites are shown on the circle map. The coding strand for transcription from the T7 promoter is shown below. The f1 origin in pT7Blue-2 is oriented so that infection with helper phage will produce virions containing single stranded DNA that is complementary to the strand shown below. Therefore, single stranded sequencing should be performed using the T7 promoter primer (Cat. No. 69348-3), Blue2S•Tag primer (Cat. No. 69087-3), or SP6 promoter primer (Cat. No. 69349-3), depending on the location of the insert.

## pT7Blue-2 sequence landmarks

<i>lacZ</i> start codon	1
<i>lacZ</i> $\alpha$ -peptide	1-498
T7 promoter	24-40
T7 transcription start	41
<i>Xenopus</i> globin 5'UTR	52-94
S•Tag coding sequence	112-156
multiple cloning region	
( <i>MscI</i> - <i>NotI</i> )	92-334
SP6 promoter	236-252
SP6 transcription start	253
f1 origin	500-955
<i>bla</i> coding sequence	1087-1944
pUC origin	2705

Notes: the *SrfI* and *SmaI* sites are destroyed during Ligation Independent Cloning.

Primer sequence extensions required for LIC compatibility are underlined in the diagram below.



# pT7Blue-2 Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations
AccI	1	269	EaeI	5	92 328 344 1481 2923	ScrFI	10	190 191 380 1338 1689
AccIII	3	564 1570 2810	EagI	1	328			2385 2603 2616 2737 3025
Acil	36		Eam1105I	1	1874	SfaNI	4	1173 1422 1613 2665
AfiIII	4	214 305 307 2762	EarI	3	437 1075 2879	Sfcl	7	36 208 248 943 1628
AgeI	1	160	Ecil	3	1718 2546 2692			2306 2497
AluI	20		Eco57I	2	1208 2220	SmaI	1	191
AlwI	10	223 236 1235 1239 1556	EcoNI	1	236	SnaBI	1	222
		2019 2020 2116 2118 2204	EcoO109I	1	315	SrfI	1	191
Alw21I	4	325 1206 1291 2452	EcoRI	1	253	Sspl	2	516 1069
Alw44I	2	1202 2448	EcoRII	5	378 2601 2614 2735 3023	StyI	2	98 286
AlwNI	1	2353	EcoRV	1	262	TaqI	7	134 149 269 281 757
Apal	1	319	FauI	5	457 861 930 2922 2964			1220 2664
Apol	4	130 253 526 537	FokI	4	398 1435 1722 1903	TaqII	7	91 628 1172 1189 1342
AvaI	3	189 280 318	FspI	2	478 1651			1527 2866
Avall	2	1510 1732	GdIII	4	328 344 1481 2923	TfiI	2	2788 2928
AvrII	1	286	HaeI	4	94 2288 2740 2751	ThaI	13	
BaeI	1	156	HaeII	4	875 883 2522 2892	TseI	15	
BamHI	1	228	HaeIII	15		Tsp45I	5	241 364 897 1401 1612
BanI	4	157 761 1921 3018	HgaI	5	222 941 1342 2072 2650	Tsp509I	12	
BanII	4	190 319 325 799	HgiEII	1	2180	Tth111III	3	2139 2171 2178
BbvI	15		HhaI	20		UbaII	10	328 331 926 940 1130
BccI	5	714 731 1425 1712 1836	Hin4I	3	264 1799 1873			1359 1481 2565 2720 2838
Bce83I	4	1270 2138 2379 2677	HincII	1	270	VspI	3	1699 2934 2993
BceII	3	332 750 2263	HindIII	1	274	XbaI	1	297
Bcgl	2	1336 1370	Hinfl	8	30 651 673 1875 2392	XhoI	1	280
Bfal	7	287 292 298 875 1681			2788 2863 2928	XmnI	1	1274
		2016 2269	HphI	7	253 725 1147 1182 1388			
BglI	2	488 1756			1804 2031			
BmgI	1	317	KpnI	1	161			
BpmI	1	1805	Maell	11				
BsaI	1	1808	MaellI	15				
BsaAI	3	222 308 724	MbolI	8	454 864 1092 1201 1279			
BsaHI	1	1334			2034 2105 2896			
BsaJI	6	98 189 286 378 2602	MluI	1	214			
		3023	MmeI	3	701 2369 2553			
BsaWI	4	160 1578 2409 2556	MnlI	12				
BsaXI	2	672 2908	MscI	1	94			
Bsbl	1	631	MseI	18				
BscGI	6	818 1326 1848 1872 2093	MslI	4	308 1103 1462 1621			
		2439	MspI	13				
BseRI	1	189	MspA1I	7	121 155 428 1238 2179			
Bsil	2	1205 2589			2424 2942			
BsiEI	6	331 459 1356 1505 2428	MwoI	15				
		2852	NciI	5	190 191 1338 1689 2385			
BsII	9	238 264 620 946 2284	NcoI	1	98			
		2563 2729 2747 2921	NgoAIV	1	825			
BsmAI	2	1032 1808	NheI	1	291			
BsoFI	25		NlaIII	9	9 102 1038 1431 1467			
Bsp24I	4	2068 2100 2246 2278			1545 1555 2046 2766			
Bsp1286I	7	190 319 325 799 1206	NlaIV	15				
		1291 2452	Nott	1	328			
BspLU11I	1	2762	NspI	1	2766			
BsrI	12		NspV	1	134			
BsrBI	4	868 1032 2833 3074	Pfi1108I	1	1855			
BsrDI	2	1640 1814	PleI	6	24 659 667 1883 2386			
BsrFI	3	160 825 1789			2871			
BstXI	1	310	PmlI	1	308			
BstYI	7	228 1227 1244 2012 2024	Psp1406I	3	509 1272 1645			
		2110 2121	PstI	1	212			
Cac8I	17		PvuI	2	459 1505			
CjeI	14		PvuII	2	428 2942			
CjePI	10	1470 1503 1942 1975 2062	RcaI	2	1034 2042			
		2095 2208 2240 2241 2273	RleAI	1	300			
CviJI	56		RsaI	2	159 1393			
CviRI	11		SacI	1	325			
DdeI	5	71 1373 1913 2079 2488	Sall	1	268			
DpnI	15		SapI	1	2879			
DraI	3	1296 1988 2007	Sau96I	8	315 316 447 715 1510			
DrallI	1	724			1732 1749 1828			
DrdI	2	679 2660	Sau3AI	15				
DrdII	1	729	Scal	1	1393			
Dsal	1	98						

Enzymes that do not cut pT7Blue-2:

AatII	AflII	ApaBI	AscI	BbsI
BclI	BglII	Bpu10I	Bpu1102I	BsaBI
BsgI	BsmI	BsmBI	BsmFI	BspEI
BspGI	BspMI	BsrGI	BssHII	Bst1107I
BstEII	Bsu36I	Clal	Eco47III	FseI
HpaI	MunI	NarI	NdeI	NruI
NsiI	Pacl	PfiMI	PmeI	PshAI
Psp5II	RsrII	SacII	SexAI	SfiI
SgfI	SgrAI	SpeI	SphI	Sse8387I
StuI	SunI	Swal	Tth111I	XcmI