

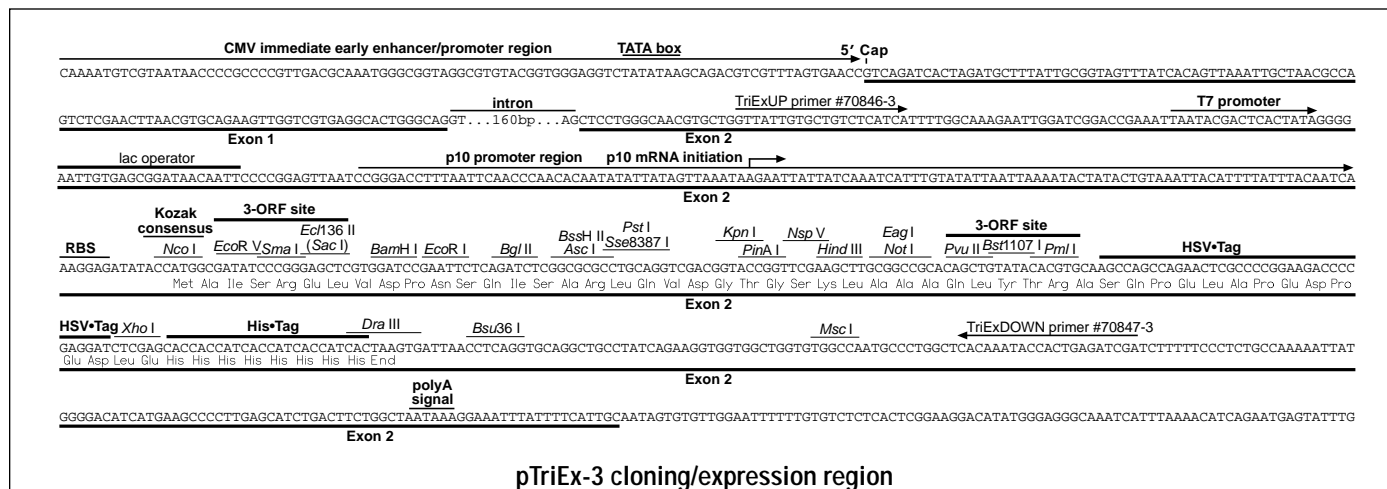
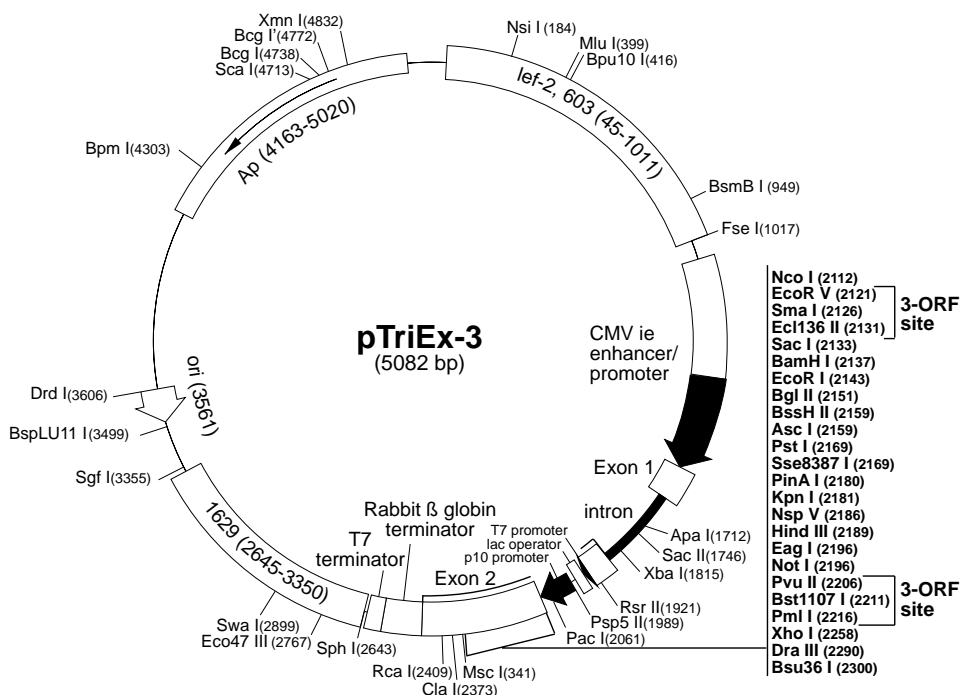
pTriEx-3 Vector

Baculovirus Locus	polh
Promoters	CMV immediate early p10 T7lac
C-terminal fusion options	HSV•Tag His•Tag
Cloning options	polylinker

The pTriEx™-3 vector¹ (cat. no. 70823-3) is designed to allow rapid characterization of target genes in multiple expression systems. With this vector a single recombinant plasmid can be used to test expression in *E. coli*, insect and vertebrate cells. Transient vertebrate expression is mediated by the CMV immediate early enhancer and promoter². For expression in insect cells, pTriEx-3 contains flanking baculovirus sequences to permit the generation of recombinant baculoviruses using the BacVector™ System. In baculovirus-infected insect cells, expression is driven by the very late p10 promoter. Expression in *E. coli* is regulated by the tightly controlled T7lac promoter. Expression can be induced in hosts such as NovaBlue by infecting with λCE6, a phage that constitutively expresses T7 RNA polymerase from the λ_{p_l} and λ_{p_l} promoters. Alternatively, pTriEx recombinant plasmids can be transferred into a (DE3)pLacI host that allows IPTG based induction.

pTriEx-3 sequence landmarks

CMV ie enhancer/promoter	1021–1597
Vertebrate transcription start	1598
T7 promoter	1931–1947
T7 transcription start	1948
lac operator	1952–1972
p10 promoter region	1986–2099
p10 transcription start	2030–2031
Multiple cloning sites (Nco I–Dra III)	2112–2290
HSV•Tag® coding sequence	2222–2257
His•Tag® coding sequence	2264–2287
Rabbit globin terminator region	2375–2581
T7 terminator	2585–2632
pUC origin	3561
bla coding sequence	4163–5020



pTriEx-3 cloning/expression region

pTriEx-3 Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AatII	6	1140 1193 1276 1462 1583 3090	EaeI	4	1011 2196 2339 4621	ScrFI	16			
AccI	3	244 2172 2210	EagI	1	2196	SfaNI	9	14 191 1372 1599 2434 3595 4489 4680 4929		
AcII	43		EarI	3	51 547 5028	SfiI	6	1857 1943 2165 3763 3954 4474		
AflIII	5	399 2213 3160 3310 3499	Ecl136II	1	2131	Sgfi	1	3355		
AhdI	2	499 4233	Eco47III	1	2767	Smal	1	2126		
AluI	18		Eco57I	2	4046 4900	SnaBI	2	1355 2809		
AlwI	11		EcoO109I	3	1709 1989 2601	SphI	1	2643		
Alw26I	8	285 949 1449 1659 1897 2495 4294 5070	EcoRI	1	2143	Sse8387I	1	2169		
AlwNI	2	1865 3914	EcoRII	8	1103 1296 1864 2348 2668 3525 3645 3658	Sspl	4	425 2903 3100 5037		
ApaI	1	1712	EcoRV	1	2121	StyI	2	2112 2596		
ApaLI	2	3812 4900	FauI	7	1108 1134 1301 1529 1712 1738 1775	Swal	1	2899		
ApoI	11		Fnu4HI	24		TaiI	18			
AscI	1	2159	FokI	3	4199 4380 4667	TaqI	12			
AvaI	4	1732 2124 2249 2258	FseI	1	1017	TfiI	3	446 2772 3474		
AvaII	5	1921 1989 2718 4371 4593	FspI	2	659 4455	ThaI	12			
BamHI	1	2137	HaeII	2	2769 3746	TseI	14			
BanI	3	1480 2177 4181	HaeIII	16		Tsp45I	3	1802 4489 4700		
BanII	2	1712 2133	HgaI	9	146 503 966 1541 2709 2869 3295 3609 4759	Tsp509I	38			
BbsI	3	498 2250 3268	HhaI	16		TspRI	11			
BbvI	14		HincII	3	245 1531 2173	VspI	5	1022 1930 2057 3294 4405		
BcgI	1	4738	HindIII	1	2189	XbaI	1	1815		
BcgI'	1	4772	Hinfl	9	156 446 1425 1937 2772 3148 3474 3869 4228	XhoI	1	2258		
Bfal	9	491 804 1606 1816 2585 2769 3039 3993 4423	HphI	8	183 879 2264 2270 4303 4719 4925 4960	XmnI	1	4832		
BglI	4	1105 1227 1298 4353	KpnI	1	2181					
BglIII	1	2151	MaeIII	13		Enzymes that do not cut pTriEx-3:				
Bpml	1	4303	MboII	15		AflIII	AvrII	BclI	Bpu1102I	BsaBI
Bpu10I	1	416	MluI	1	399	BseRI	BsmI	BspEI	BstEII	EcoNI
Bsal	2	285 4294	MnII	23		EheI	HpaI	NarI	NheI	NruI
BsaAI	3	1355 2216 2809	MscI	1	2341	PfiI	PmeI	PshAI	SanDI	SapI
BsaHI	9	495 1137 1190 1273 1459 1580 3087 3287 4770	MseI	37		SexAI	SfiI	SgrAI	SpeI	SrfI
BsaJI	11		MslI	7	950 1380 1837 2538 4485 4644 5003	StuI	SunI	Tth111I	UbaEI	XcmI
BsaWI	5	833 2180 3704 3851 4524	MspI	20						
BsgI	2	1689 2326	MspA1I	6	655 1745 2206 3840 4085 4868					
BsiEI	6	2199 3355 3415 3838 4603 4752	MunI	2	2798 3143					
BsiHKAI	5	2133 2265 3816 4819 4904	MwoI	18						
BsII	10	264 1064 1718 1766 2610 2976 3521 3539 3704 3983	NciI	8	1975 1987 2125 2126 2240 3878 4416 4767					
BsmBI	1	949	NcoI	1	2112					
BsmFI	8	1190 1341 1509 1764 1783 2002 2417 3215	NdeI	4	1249 2509 2569 2577					
Bsp1286I	6	1712 2133 2265 3816 4819 4904	NgoAIV	2	781 1013					
BspLU11I	1	3499	NlaIII	14						
BspMI	2	1686 2158	NlaIV	12						
BsrI	13		NotI	1	2196					
BsrBI	3	1728 1960 3432	Nsil	1	184					
BsrDI	4	79 2460 4294 4468	NspI	2	2643 3503					
BsrFI	6	42 781 1013 1806 2180 4313	NspV	1	2186					
BsrGI	3	49 768 3139	Pacl	1	2061					
BssHII	1	2159	PinAI	1	2180					
BssSI	3	2132 3671 4897	PleI	6	150 1419 1931 3142 3877 4222					
Bst1107I	1	2211	PmlI	1	2216					
BstXI	2	167 2349	Psp1406I	2	4459 4832					
BstYI	7	2137 2151 2254 4139 4150 4858 4875	Psp5II	1	1989					
Bsu36I	1	2300	PstI	1	2169					
Cac8I	18		PvuI	2	3355 4603					
Clal	1	2373	PvuII	1	2206					
CviJI	66		RcaI	1	2409					
Ddel	8	416 2148 2287 2300 2367 3773 4190 4730	Rsal	15						
Dpnl	20		RsrII	1	1921					
Dral	4	429 2529 2899 4810	SacI	1	2133					
DrallI	1	2290	SacII	1	1746					
DrdI	1	3606	Sall	2	243 2171					
Dsal	2	1743 2112	Sau3AI	20						
			Sau96I	12						
			Scal	1	4713					