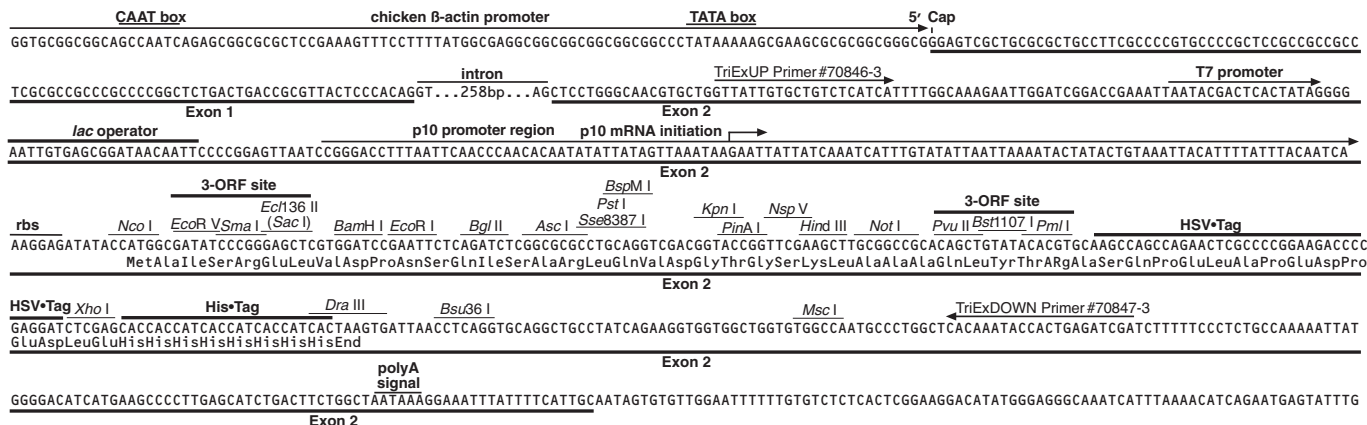
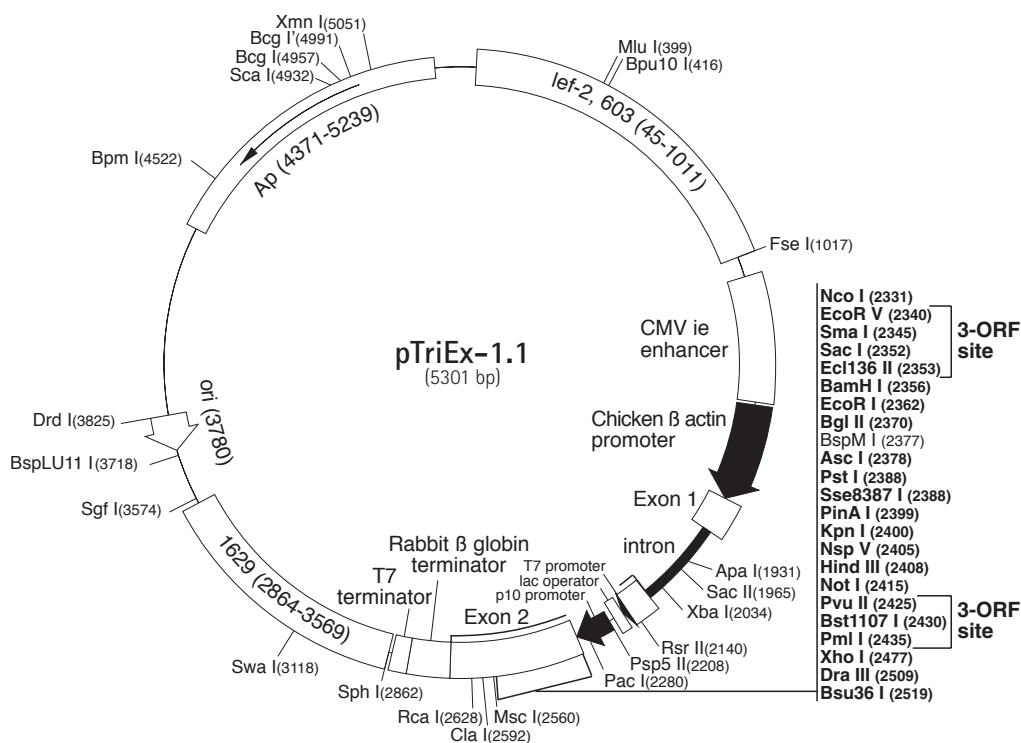


pTriEx™ -1.1 Vector

	Cat No.
pTriEx-1.1 Vector DNA	70840-3
pTriEx-1.1 sequence landmarks	
CMV ie enhancer region	1079-1443
Chicken actin promoter region	1449-1726
Vertebrate transcription start	1727
T7 promoter	2150-2166
T7 transcription start	2167
lac operator	2171-2191
p10 promoter region	2205-2318
p10 transcription start	2249-2250
Multiple cloning sites (Nco I-Dra III)	2331-2512
HSV•Tag® coding sequence	2441-2476
His•Tag® coding sequence	2483-2506
Rabbit globin terminator region	2594-2800
T7 terminator	2804-2851
pUC origin	3780
bla coding sequence	4371-5239

The pTriEx™-1.1 vector¹ (cat. no. 70840-3) is uniquely 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 a hybrid promoter composed of the CMV immediate early enhancer fused to the chicken β-actin promoter. For expression in insect cells, pTriEx-1.1 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 λp_L and λp_R promoters. Alternatively, pTriEx recombinant plasmids can be transferred into a (DE3)pLacI host that allows IPTG based induction.

¹ US Patent 6,589,783



pTriEx-1.1 cloning/expression regions

pTriEx™-1.1 Restriction Sites

TB280 0607

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations
AatII	4	1204 1257 1340 3309	DsaI	2	1962 2331	ScrFI	19	
AccI	3	244 2391 2429	EaeI	5	1011 1015 2415 2558 4840	SfaNI	7	14 191 2653 3814 4708
AcII	64		EagI	2	1015 2415			4899 5148
AfIII	5	399 2432 3379 3529 3718	EarI	3	51 547 5247	Sfcl	6	2076 2162 2384 3982 4173
AhdI	2	499 4452	Ecl136II	1	2350			4693
AluI	18		Eco47III	2	1862 2986	Sgfl	1	3574
AlwI	11		Eco57I	2	4265 5119	SmaI	1	2345
Alw26I	7	285 949 1019 2116 2714	EcoO109I	4	1927 1928 2208 2820	SnaBI	2	1419 3028
		4513 5289	EcoRI	1	2362	SphI	1	2862
AlwNI	2	2084 4133	EcoRII	9	1167 1360 1576 2083 2567	Sse8387I	1	2388
Apal	1	1931			2887 3744 3864 3877	Sspl	4	425 3122 3319 5256
ApaLI	2	4031 5119	EcoRV	1	2340	StyI	2	2331 2815
ApoI	11		FauI	18		Swal	1	3118
AscI	1	2378	Fnu4HI	41		Tail	16	
AvaI	4	1951 2343 2468 2477	FokI	3	4418 4599 4886	TaqI	12	
Avall	5	2140 2208 2937 4590 4812	FseI	1	1017	TfiI	3	446 2991 3693
BamHI	1	2356	FspI	2	659 4674	Thal	20	
BanI	2	2396 4400	HaeII	3	1864 2988 3965	TseI	19	
BanII	4	1460 1920 1931 2352	HaeIII	18		Tsp45I	3	2021 4708 4919
BbsI	3	498 2469 3487	HgaI	9	146 503 966 1036 2928	Tsp509I	40	
BbvI	19				3088 3514 3828 4978	TspRI	10	690 2590 2939 3352 4125
BcgI	1	4957	HhaI	27				4138 4400 4505 4852 4879
BcgI'	1	4991	HincII	2	245 2392	VspI	5	1086 2149 2276 3513 4624
Bfal	9	491 804 1079 2035 2804	HindIII	1	2408	XbaI	1	2034
		2988 3258 4212 4642	Hinfl	9	156 446 1728 2156 2991	XhoI	1	2477
BglI	4	1169 1291 1362 4572			3367 3693 4088 4447	XmnI	1	5051
BglIII	1	2370	HphI	10	183 879 1465 1831 2483			
Bpml	1	4522			2489 4522 4938 5144 5179			
Bpu10I	1	416	KpnI	1	2400			
BsaI	2	285 4513	MaellI	14				
BsaAI	3	1419 2435 3028	MbolI	14				
BsaHI	7	495 1201 1254 1337 3306	MluI	1	399			
		3506 4989	MnlI	30				
BsaJI	12		MscI	1	2560			
BsaWI	5	833 2399 3923 4070 4743	MseI	37				
BsgI	2	1560 2545	MsiI	6	950 2056 2757 4704 4863			
BsiEI	7	1018 2418 3574 3634 4057			5222			
		4822 4971	MspI	22				
BsiHKA1	5	2352 2484 4035 5038 5123	MspA1I	6	655 1964 2425 4059 4304			
BsII	10	264 1128 1937 1985 2829			5087			
		3195 3740 3758 3923 4202	MunI	2	3017 3362			
BsmBI	2	949 1019	MwoI	27				
BsmFI	8	1254 1405 1844 1983 2002	NciI	10	1789 1922 2194 2206 2344			
		2221 2636 3434			2345 2459 4097 4635 4986			
Bsp1286I	9	1460 1758 1920 1931 2352	NcoI	1	2331			
		2484 4035 5038 5123	NdeI	4	1313 2728 2788 2796			
BspLU11I	1	3718	NgoAIV	2	781 1013			
BspMI	1	2377	NlaIII	15				
BsrI	12		NlaIV	12				
BsrBI	6	1647 1761 1824 1947 2179	NotI	1	2415			
		3651	Nsil	2	184 1445			
BsrDI	4	79 2679 4513 4687	NspI	2	2862 3722			
BsrFI	6	42 781 1013 2025 2399	NspV	1	2405			
		4532	Pacl	1	2280			
BsrGI	3	49 768 3358	PinAI	1	2399			
BssHII	6	1570 1572 1650 1713 1736	PleI	6	150 1736 2150 3361 4096			
		2378			4441			
BssSI	3	2351 3890 5116	PmlI	1	2435			
Bst1107I	1	2430	Psp1406I	2	4678 5051			
BstXI	2	167 2568	Psp5II	1	2208			
BstYI	7	2356 2370 2473 4358 4369	PstI	1	2388			
		5077 5094	PvuI	2	3574 4822			
Bsu36I	1	2519	PvuII	1	2425			
Cac8I	26		RcaI	1	2628			
Clal	1	2592	RsaI	14				
CviJI	76		RsrII	1	2140			
Ddel	8	416 2367 2506 2519 2586	SacI	1	2352			
		3992 4409 4949	SacII	1	1965			
Dpnl	19		Sall	2	243 2390			
DraI	4	429 2748 3118 5029	Sau3AI	19				
DraIII	1	2509	Sau96I	14				
DrdI	1	3825	Scal	1	4932			

Enzymes that do not cut pTriEx-1.1:

AfIII	AvrII	BclI	Bpu1102I	BsaBI	BseRI
		BsmI	BspEI	BstEII	EcoNI
		HpaI	NarI	NheI	NruI
		PmeI	PshAI	SanDI	SapI
		SfiI	SgrAI	SpeI	SrfI
		SunI	Tth111I	UbaEI	XcmI