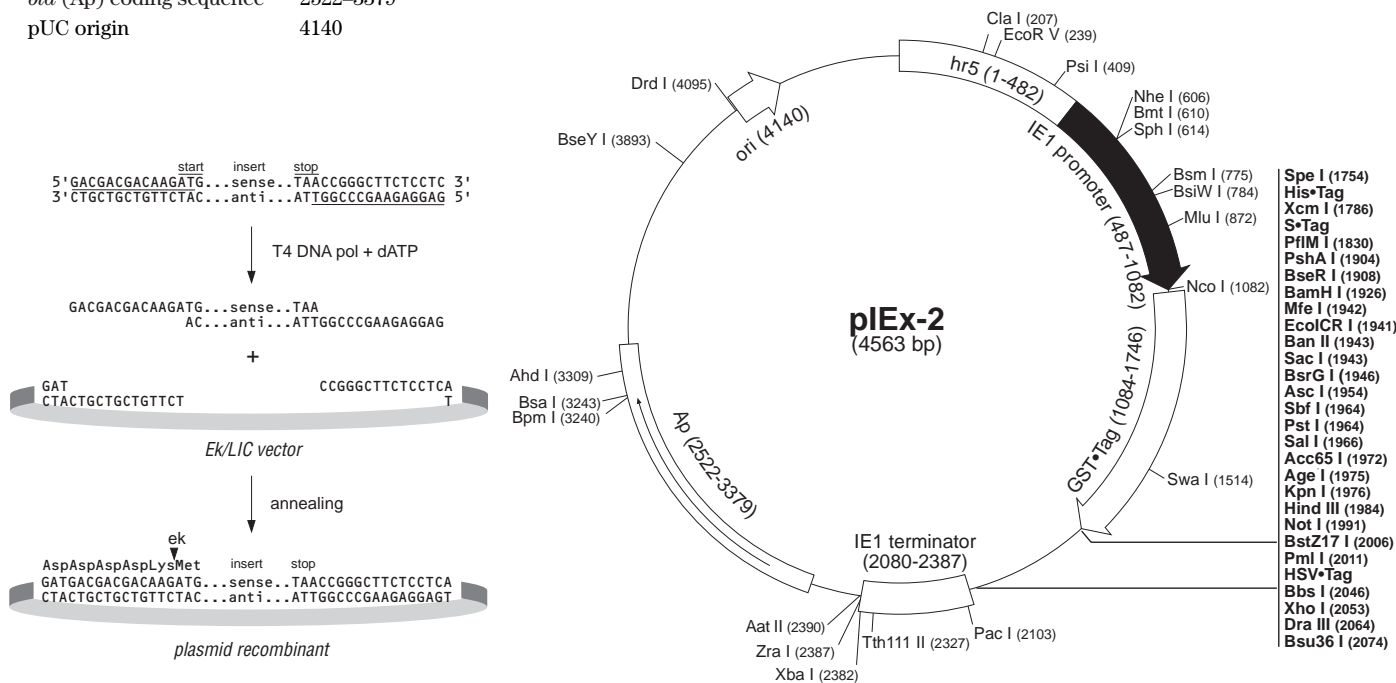


pIEx-2 Vector

	Cat. No.
pIEx-2 DNA	71238-3
pIEx-2 Ek/LIC Vector	71239-3
pIEx-2 sequence landmarks	
hr5 enhancer	1-482
IE1 promoter	487-1082
IE1 transcription start	1034
GST•Tag coding sequence	1084-1746
His•Tag® coding sequence	1768-1785
S•Tag™ coding sequence	1795-1839
Multiple cloning sites (<i>PshA I-Dra III</i>)	1900-2079
HSV•Tag® coding sequence	2017-2052
IE1 terminator	2080-2387
<i>bla</i> (Ap) coding sequence	2522-3379
pUC origin	4140

The pIEx™ vectors are designed for cloning and high-level expression of proteins in transiently transfected *Spodoptera*-derived insect cells. Transcription is driven by the AcNPV-derived hr5 enhancer and immediate early promoter, IE1. pIEx-2 contains N-terminal GST•Tag, His•Tag, and S•Tag coding sequences, and a C-terminal HSV•Tag coding sequence. Unique sites are shown on the circle map. N-terminal vector encoded sequences can be completely removed by cloning into the *PshA I* site and cleaving the fusion protein with enterokinase.

pIEx-2 is also available as an Ek/LIC vector (Cat. No. 71239-3). The Ek/LIC vector is prepared for rapid directional cloning of PCR-amplified DNA. Using specifically designed primers for amplification, inserts can be efficiently cloned without the need for restriction enzyme digestion or ligation (below left).



pIEx-2 cloning/expression region

pIEx-2 Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AatII	1	2390	Eco57MI	3	2643 3240 3655		
Acc65I	1	1972	EcoICRI	1	1941		
AccI	2	1967 2005	EcoNI	1	1097		
AccI	3	2337 2707 3080	EcoO109I	1	1119		
AflIII	6	192 372 872 1325 2008	EcoRI	6	83 185 292 365 472		
		4197			4553		
AgeI	1	1975	EcoRV	1	239		
AhdI	1	3309	FspI	2	654 3086		
AlwNI	2	1879 3788	HaeII	2	3957 4327		
ApaLI	2	2637 3883	HincII	2	1044 1968		
AscI	1	1954	HindIII	1	1984		
Asel	5	831 2103 3134 4369 4428	KpnI	1	1976		
AvaI	4	1862 2044 2053 4558	MfeI	1	1932		
BamHI	1	1926	MluI	1	872		
BanI	3	1972 3356 4453	MscI	1	1294		
BanII	1	1943	MspI	4	1181 2538 2897 3056		
BbsI	1	2045	NcoI	1	1082		
BceAI	6	68 174 731 1967 2136	NheI	1	606		
		3697	NotI	1	1991		
BcgI	3	1117 2221 2771	NspI	4	614 691 1329 4201		
BciVI	4	533 868 2472 3999	NspV	2	1484 1981		
BclI	2	778 1521	PacI	1	2103		
BglI	2	995 3191	PciI	2	1325 4197		
Bme1580I	2	2641 3887	PfiMI	1	1830		
BmrI	2	2365 3269	PinAI	1	1975		
BmtI	1	610	PmlI	1	2011		
Bpml	1	3240	Ppil	2	2681 3491		
BpuEI	4	2705 3573 3814 4112	PshAI	1	1904		
BsaAI	4	193 373 784 2011	PsiI	1	409		
BsaHI	2	2387 2769	Psri	3	89 298 371		
Bsal	1	3243	PstI	1	1964		
BsaWI	4	1975 3013 3844 3991	PvuI	2	207 2940		
BsaXI	3	1930 2234 4359	PvuII	3	776 2001 4377		
BseRI	1	1908	SacI	1	1943		
BseYI	1	3893	Sall	1	1966		
BsgI	1	1373	SapI	2	1166 4314		
BsiEI	7	207 845 1994 2791 2940	SbfI	1	1964		
		3863 4287	Scal	2	1660 2828		
BsiHKAI	5	1943 2060 2641 2726 3887	Sfcl	5	1960 2247 3063 3741 3932		
BsiWI	1	784	SmaI	2	1864 4560		
BsmAI	2	2467 3243	SmlI	5	2053 2684 3552 3829 4091		
BsmFI	2	774 1843	SpeI	1	1753		
BsmI	1	775	SphI	1	614		
Bsp1286I	5	1943 2060 2641 2726 3887	Sse8387I	1	1964		
BspCNI	6	838 2087 2821 3340 3506	SspI	3	973 2264 2504		
		3915	StyI	2	568 1082		
BspHI	2	2469 3477	Swal	1	1514		
BspMI	2	1864 1953	TaqII	3	2777 2962 4301		
BsrBI	5	863 1378 2467 4268 4509	TatI	4	724 1658 1946 2826		
BsrDI	2	3075 3249	TspGWI	7	552 587 636 790 1939		
BsrFI	2	1975 3224			2528 2870		
BsrGI	1	1946	Tth111I	1	2327		
BssHII	3	502 763 1954	XbaI	1	2382		
BssSI	4	793 1942 2640 4024	XcmI	1	1786		
Bst1107I	1	2006	XhoI	1	2053		
BstBI	2	1484 1981	XmaI	2	1862 4558		
BstYI	8	1926 2049 2662 2679 3447	XmnI	2	1480 2709		
		3459 3545 3556	ZraI	1	2388		
BstZ17I	1	2006					
Bsu36I	1	2074	Enzymes that do not cut pIEx-2:				
BtgI	1	1082	AarI	AfeI	AfIII	AleI	Alol
BtsI	4	727 2890 2910 4417	Apal	AsiSI	AvrII	BaeI	BbeI
ClaI	1	207	BbvCI	BfrBI	BglII	BclI	BmgBI
DraI	5	1514 1623 2731 3423 3442	BpII	Bpu10I	BsaBI	BsmBI	BspEI
DraIII	1	2064	BstAPI	BstEII	BstXI	BtrI	FalI
DrdI	1	4095	FseI	FspAI	HpaI	KasI	NaeI
EaeI	5	842 1292 1991 2916 4358	NarI	NdeI	NgoMIV	NruI	NsiI
EagI	2	842 1991	PfoI	PmeI	PpuMI	PspOMI	RsrII
EarI	4	1166 1791 2510 4314	SacII	SanDI	SexAI	SfiI	SfoI
Ecil	4	916 3167 3995 4141	SgrAI	SnaBI	SrfI	StuI	
Ecl136II	1	1941					
Eco57I	2	2643 3655					