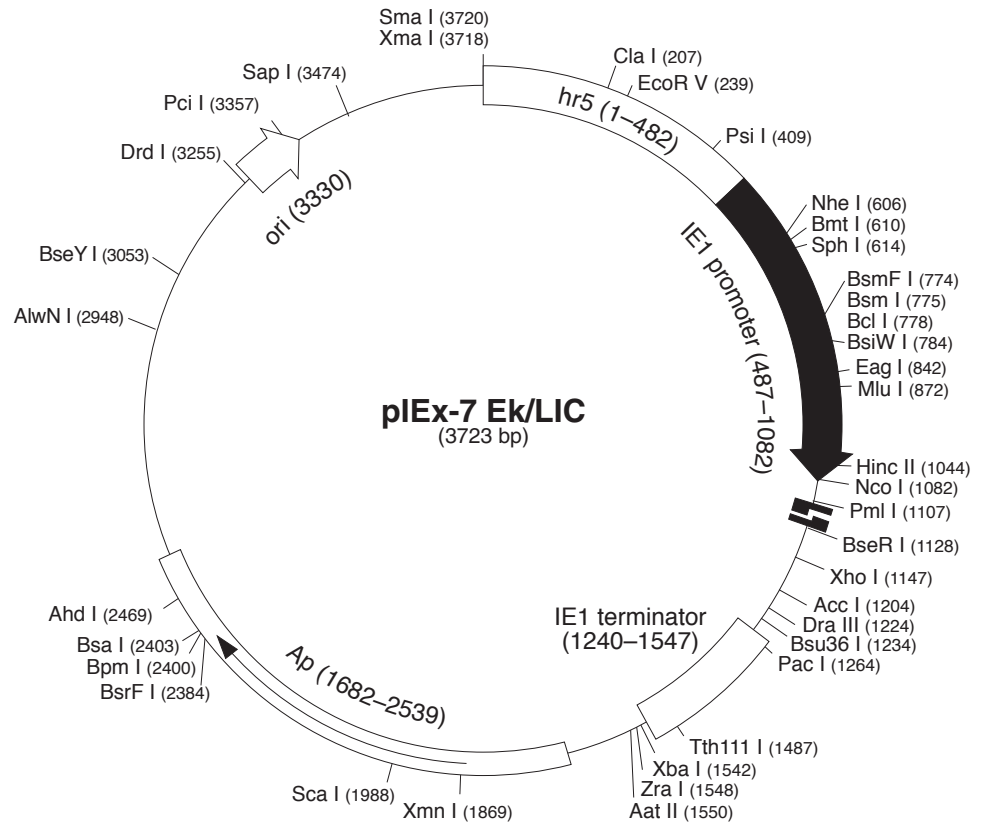


pIEx-7 Ek/LIC Vector

TB399 0903

	Cat. No.
pIEx-7 Ek/LIC Cloning Kit	71339-3
pIEx-7 Ek/LIC sequence landmarks	
hr5 enhancer	1-482
IE1 promoter	487-1082
IE1 transcription start	1034
His•Tag [®] coding sequence	1090-1107
Ek/LIC cloning site	1114-1142
S•Tag [™] coding sequence	1159-1203
IE1 terminator	1240-1547
<i>bla</i> (Ap ^R) coding sequence	1682-2539
pUC origin	3300

The pIEx[™]-7 Ek/LIC Vector is designed to allow rapid, high-level protein expression in insect cells by using transient transfection. This vector features the baculovirus hr5 enhancer and the IE1 (immediate early) promoter to direct expression in insect cells. This promoter/enhancer combination recruits endogenous insect cell transcription machinery, thereby avoiding baculovirus infection and the associated cytotoxic effects. Using specifically designed primers for amplification and the pIEx-7 Ek/LIC Cloning Kit (71339-3), inserts can be efficiently cloned without the need for restriction digestion or ligation. Fusion proteins contain a cleavable N-terminal His•Tag[®] sequence. Fusion to an optional C-terminal S•Tag[™] sequence can also be created for detection, purification and quantification of fusion proteins. Unique sites are shown on the circle map. Sequencing can be performed using the IE1 Promoter Primer (Cat. No. 69103-3) and AS S•Tag 18mer Primer (Cat. No. 71262-3).



pIEx-7 Ek/LIC cloning/expression region

