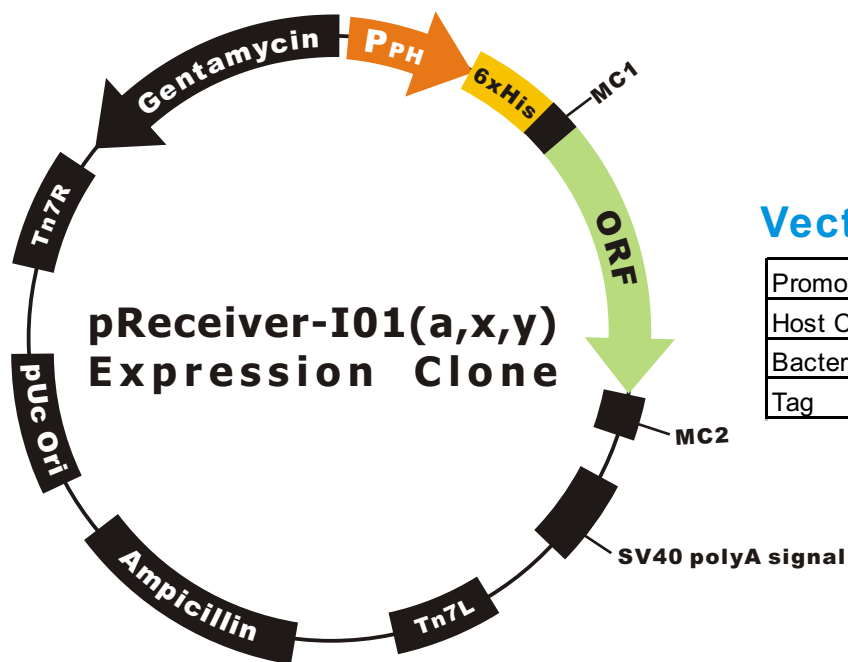


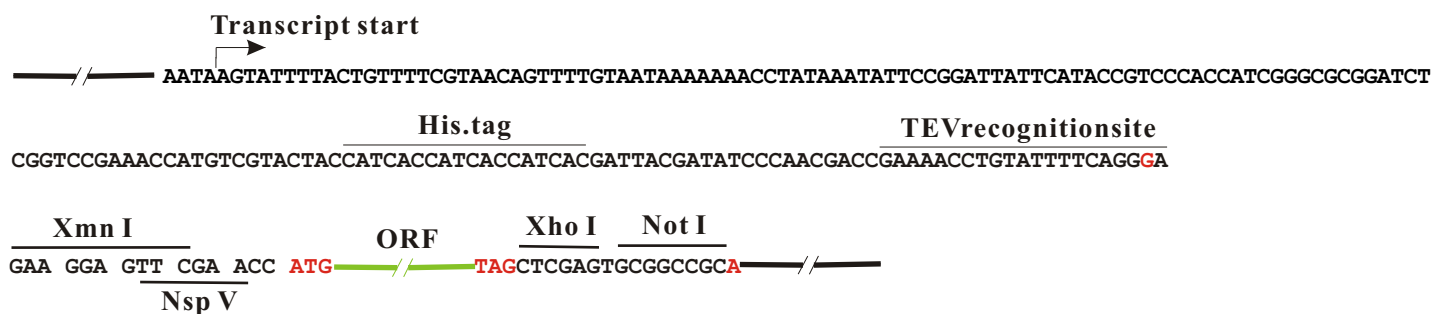
OmicsLink™ Expression Clone (AcMNPV polyhedrin promoter)



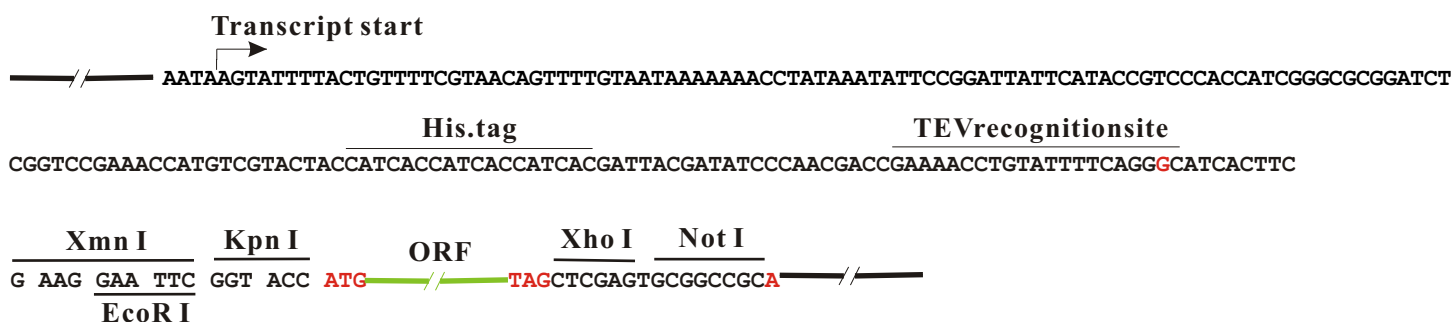
Vector Features

Promoter	AcMNPV polyhedrin
Host Cell	Insect cell
Bacterial selection antibiotic	Ampicillin
Tag	N-His

pReceiver-I01a



pReceiver-I01x



pReceiver-I01y

