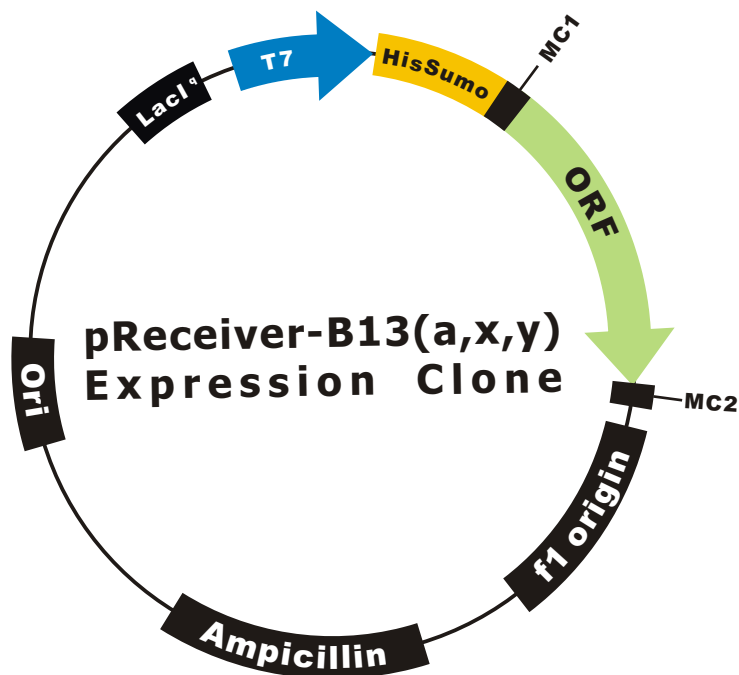


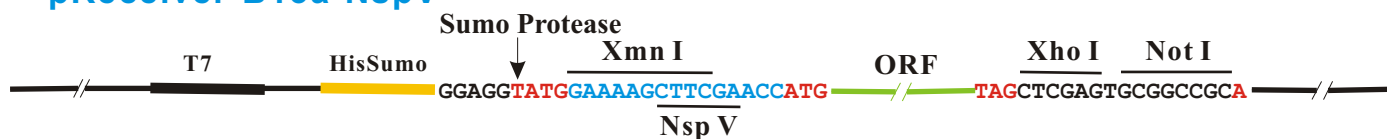
OmicLink™ Expression Clone (T7 Promoter)



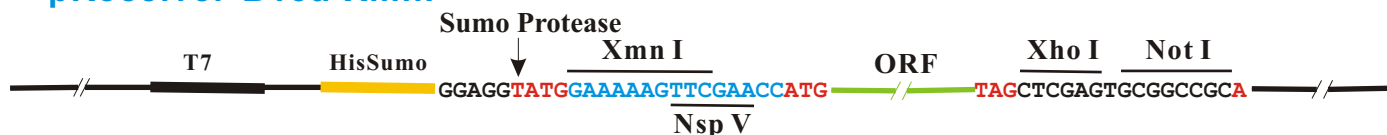
Vector Features

Promoter	T7
Host Cell	<i>E. Coli</i>
Bacterial selection antibiotic	Ampicillin
Tag	His Sumo

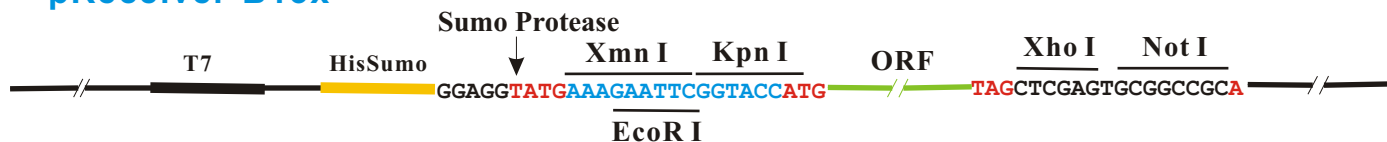
pReceiver-B13a-NspV



pReceiver-B13a-XmnI



pReceiver-B13x



pReceiver-B13y

