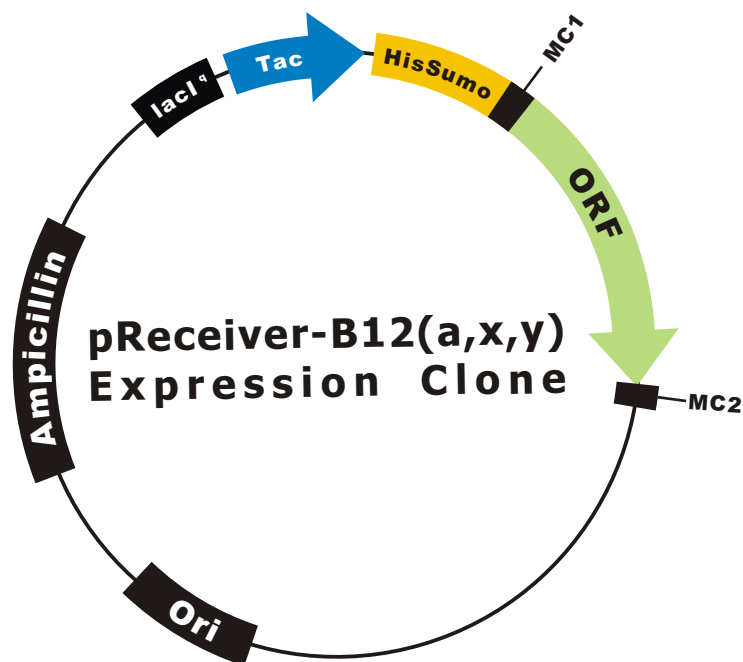


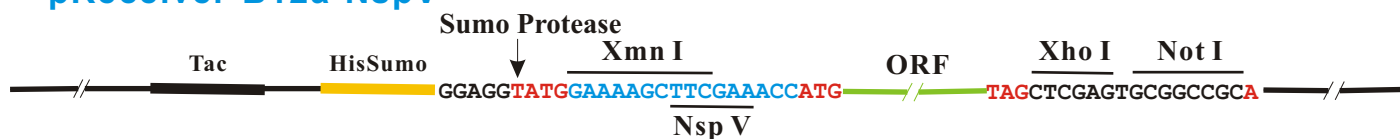
# OmicLink™ Expression Clone (Tac Promoter)



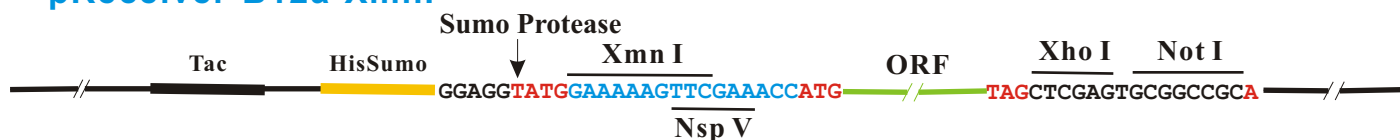
## Vector Features

Promoter	Tac
Host Cell	<i>E. Coli</i>
Bacterial selection antibiotic	Ampicillin
Tag	HisSumo

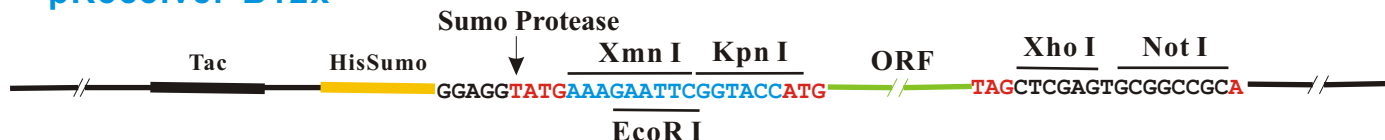
### pReceiver-B12a-NspV



### pReceiver-B12a-Xmnl



### pReceiver-B12x



### pReceiver-B12y

