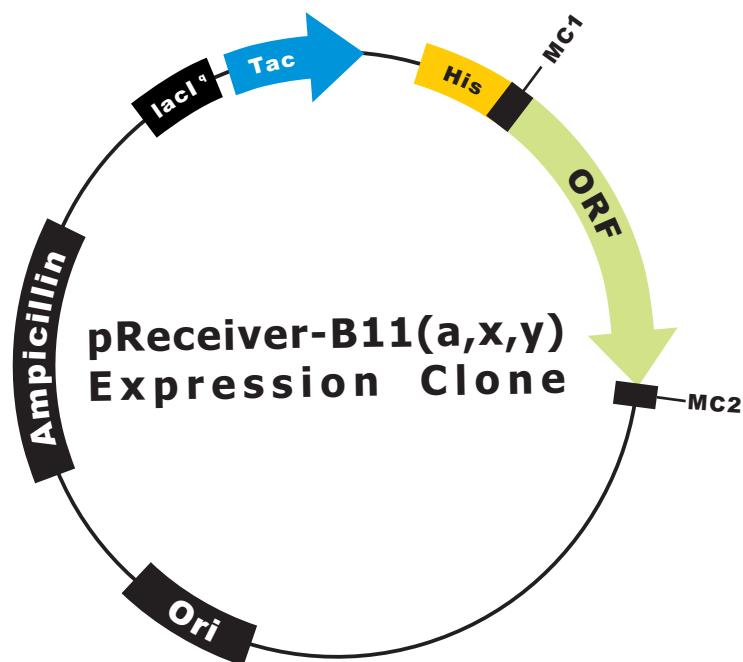


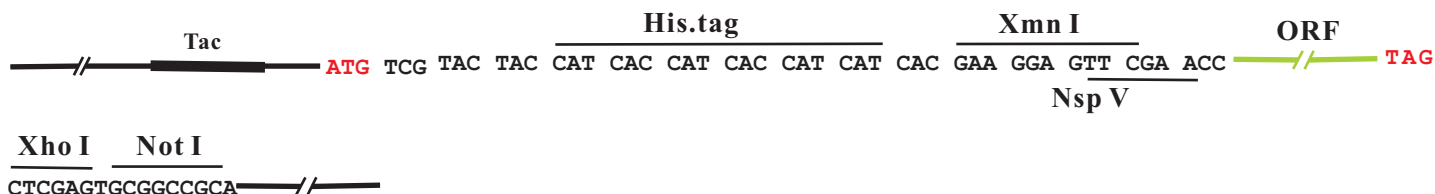
# OmicLink™ Expression Clone (Tac Promoter)



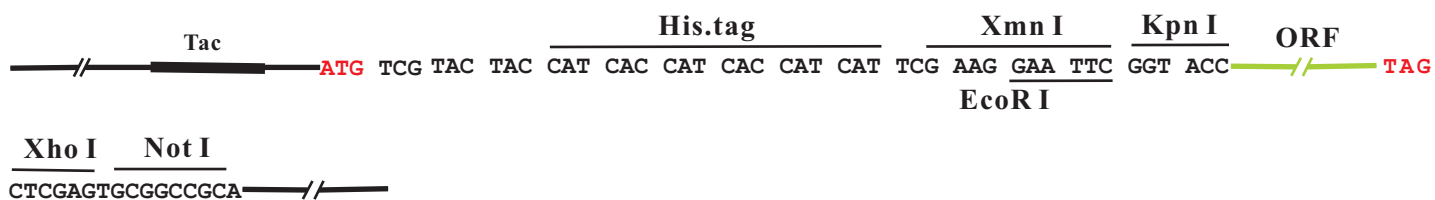
## Vector Features

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

## pReceiver-B11a



## pReceiver-B11x



## pReceiver-B11y

