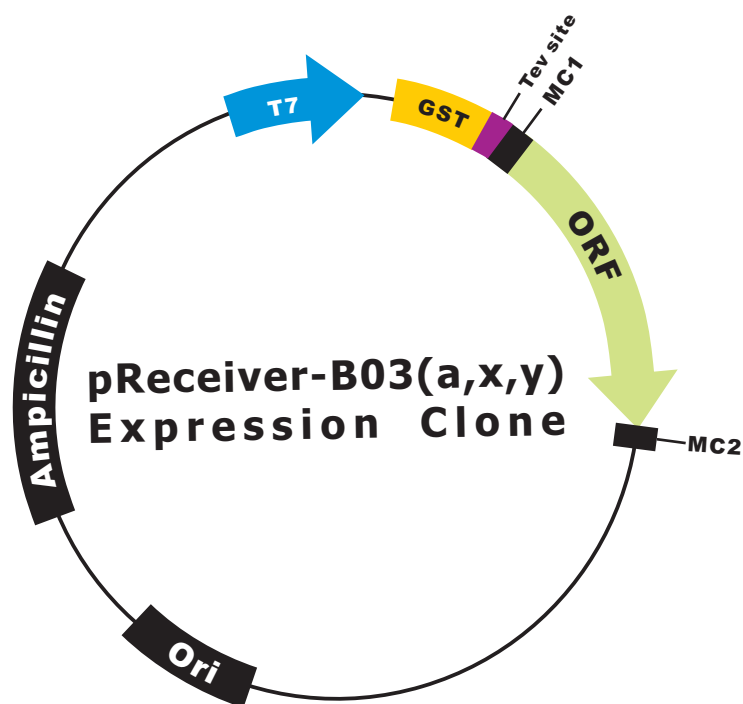


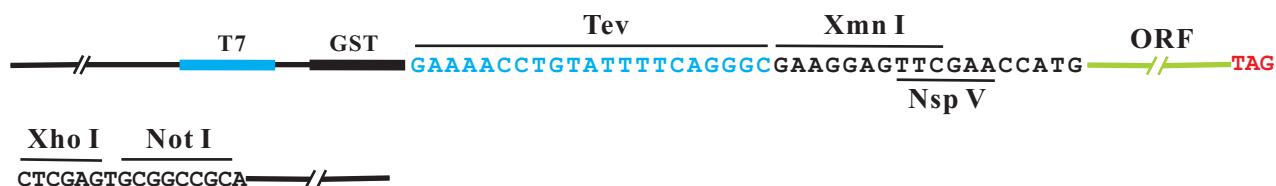
# OmicLink™ Expression Clone (T7 Promoter)



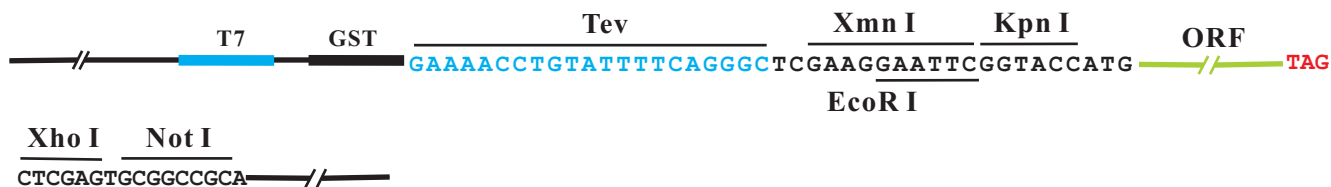
## Vector Features

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

## pReceiver-B03a



## pReceiver-B03x



## pReceiver-B03y

