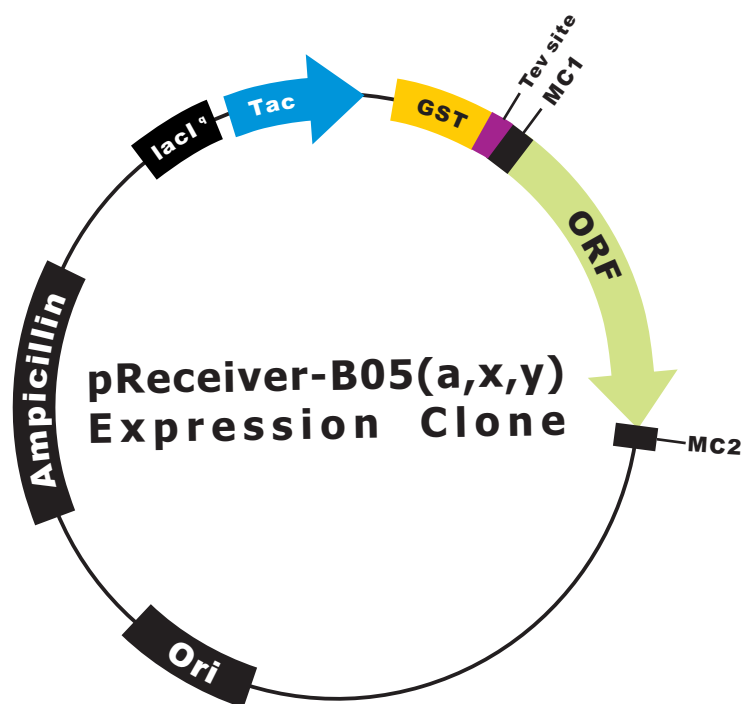


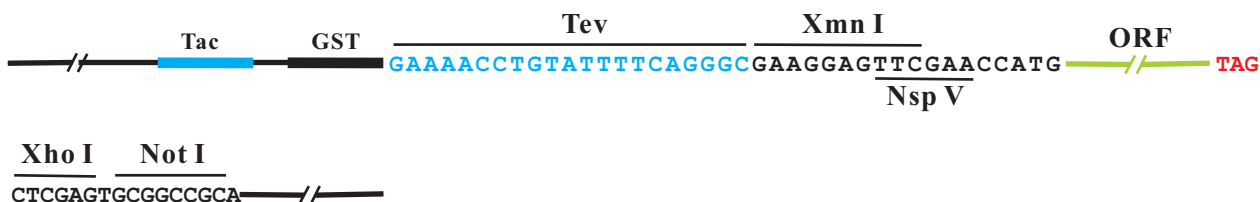
OmicLinkTM Expression Clone (Tac Promoter)



Vector Features

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

pReceiver-B05a



pReceiver-B05x



pReceiver-B05y

