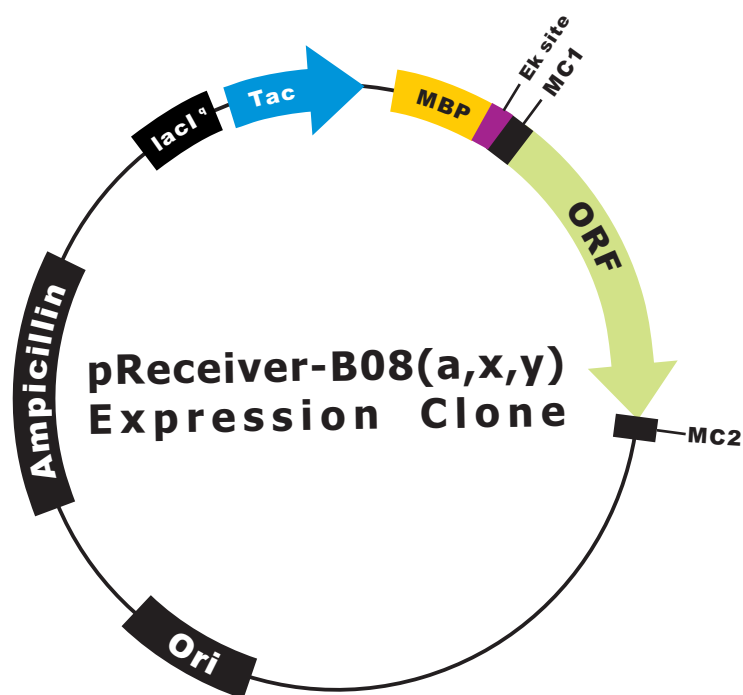


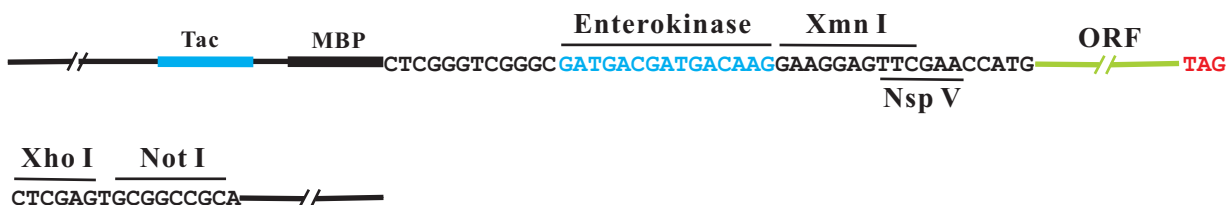
OmicLink™ Expression Clone (Tac Promoter)



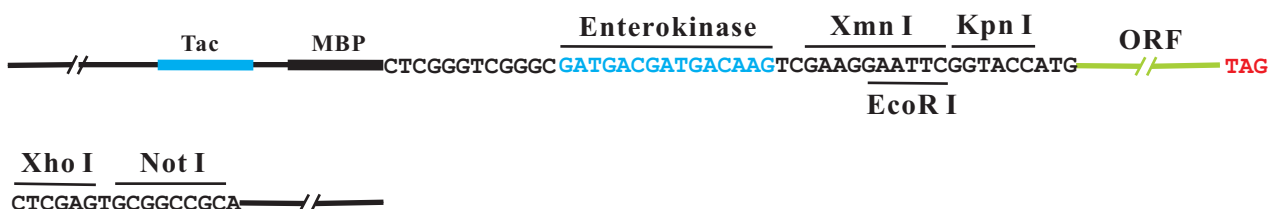
Vector Features

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

pReceiver-B08a



pReceiver-B08x



pReceiver-B08y

