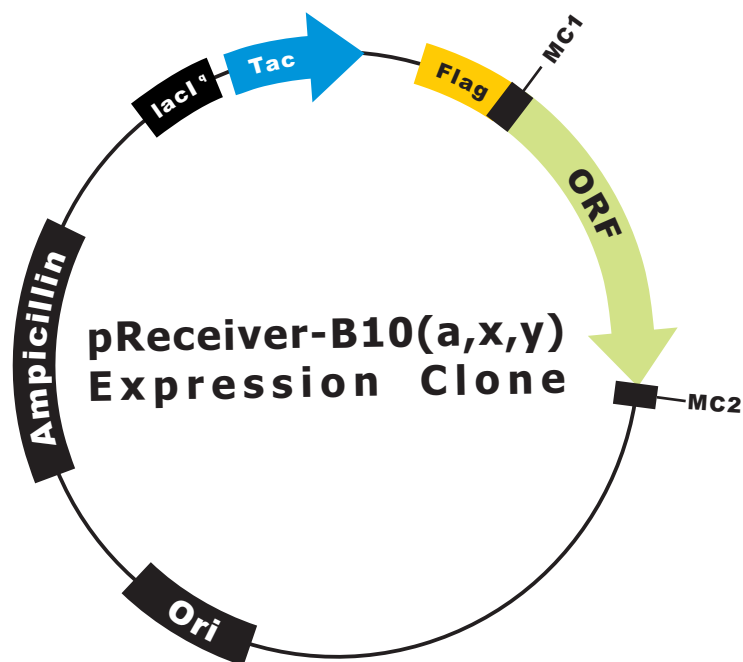


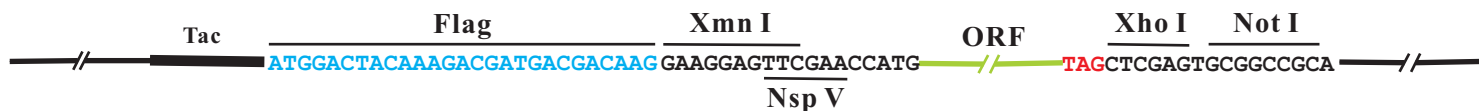
OmicLink™ Expression Clone (Tac Promoter)



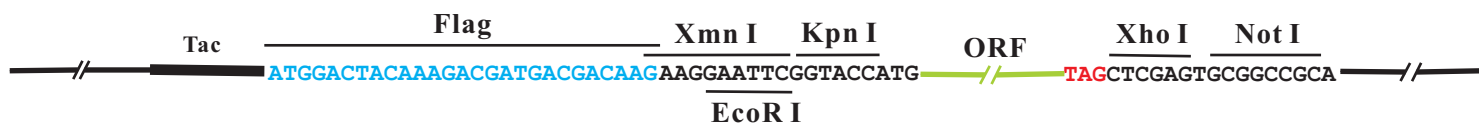
Vector Features

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

pReceiver-B10a



pReceiver-B10x



pReceiver-B10y

