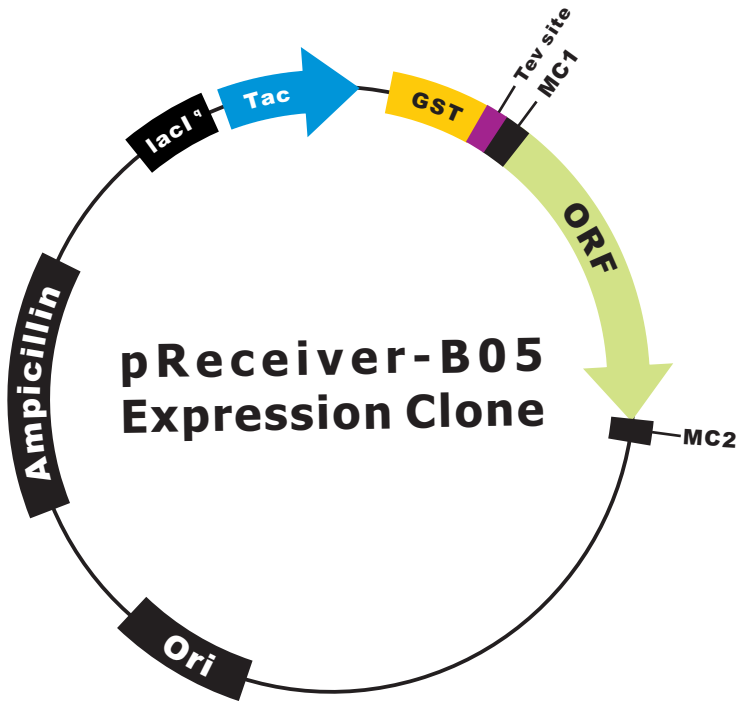


# OmicLink™ Expression Clone (Tac Promoter)



## Vector Features

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