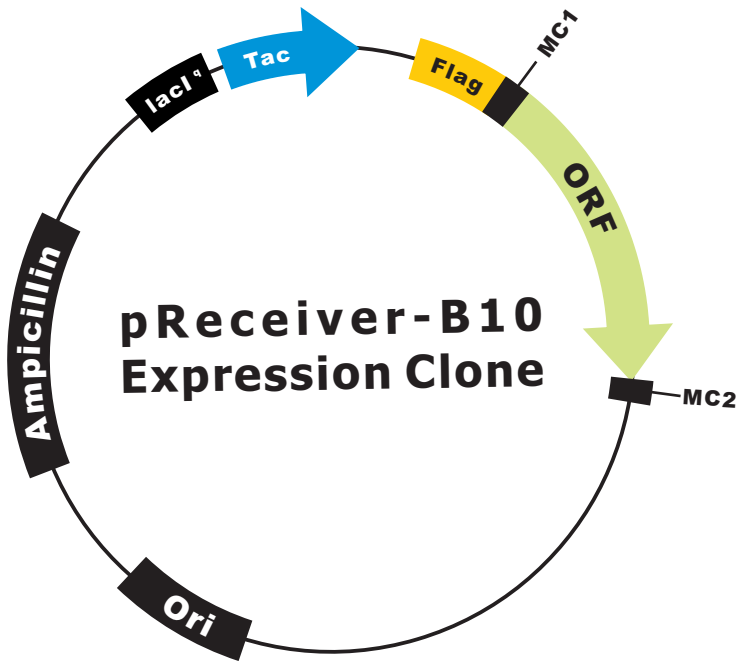


OmicsLink™ Expression Clone (Tac Promoter)



Vector Features

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