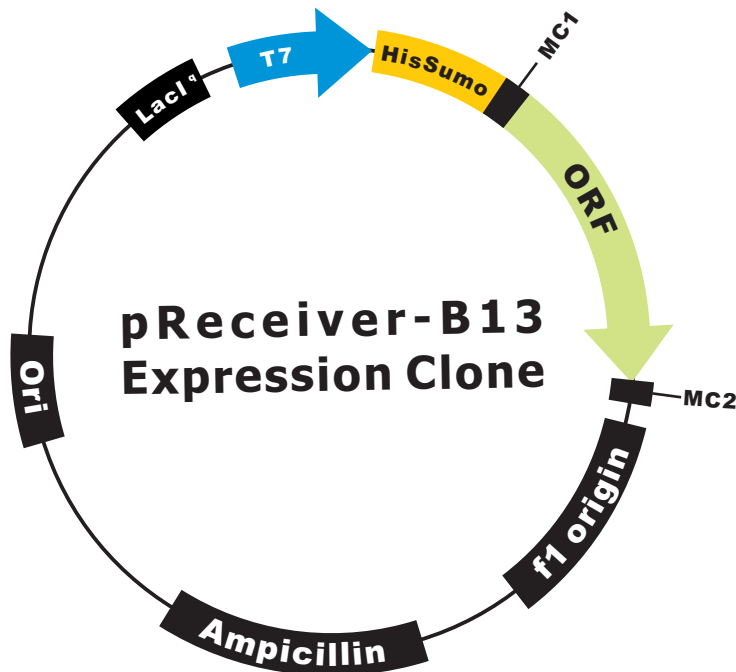


OmicLinkTM Expression Clone (T7 Promoter)



Vector Features

Promoter	T7
Host Cell	<i>E. Coli</i>
Bacterial selection antibiotic	Ampicillin
Tag	His Sumo