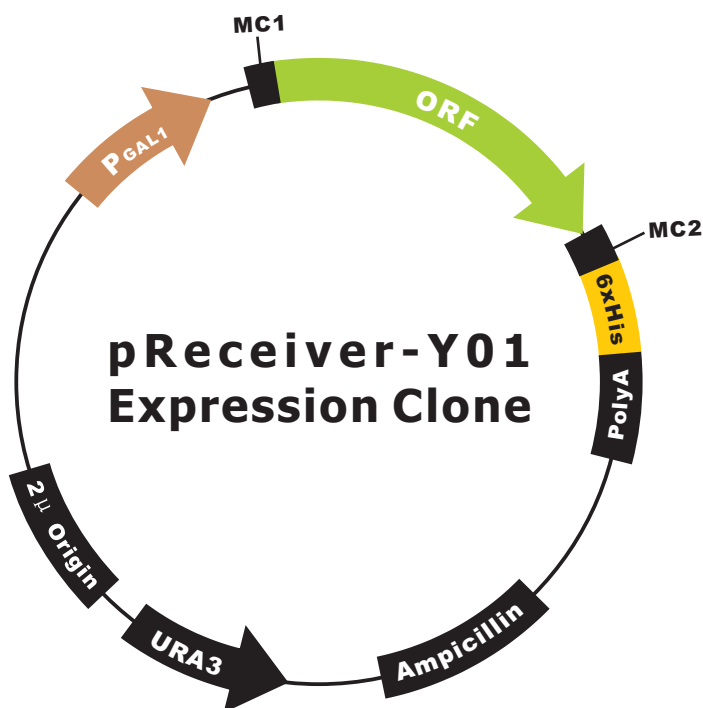


OmicLink™ Expression Clone (GAL1 Promoter)



Vector Features

Promoter	<i>GAL1</i>
Host Cell	<i>S. cerevisiae</i>
Bacterial selection antibiotic	Ampicillin
Tag	C-His