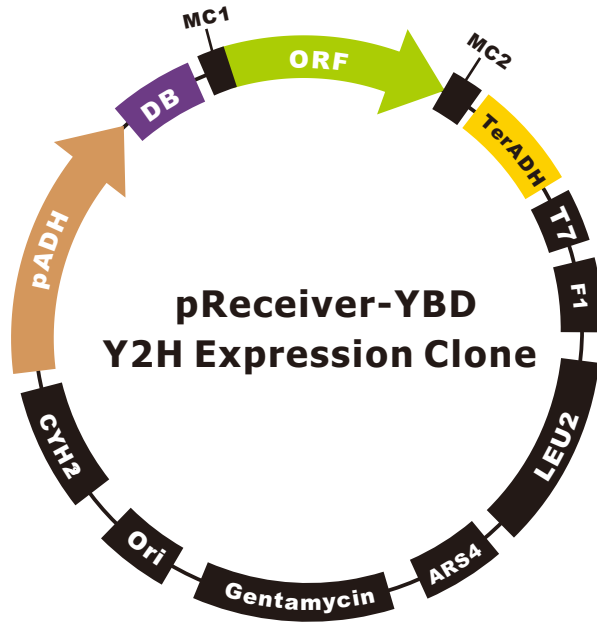


OmicLink™ Y2H Expression Clone



Vector Features

Promoter	<i>pADH</i>
Host Cell	<i>Yeast</i>
Bacterial selection antibiotic	Gentamycin
Tag	DB