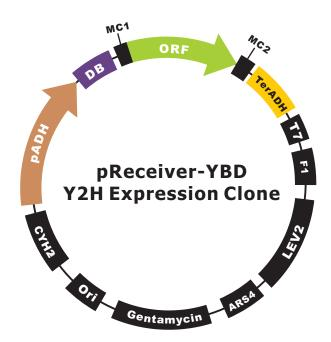


OmicsLink[™] Y2H Expression Clone



Vector Features

Promoter	pADH
Host Cell	Yeast
Bacterial selection antibiotic	Gentamycin
Tag	DB