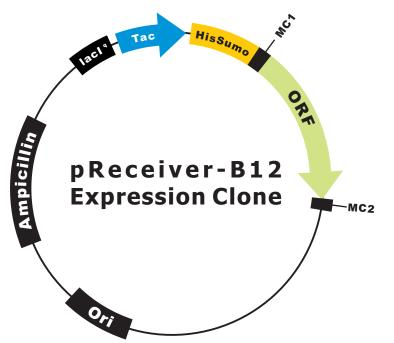


OmicsLink[™] Expression Clone (Tac Promoter)



Vector Features

Promoter	Tac
Host Cell	E. Coli
Bacterial selection antibiotic	Ampicillin
Tag	HisSumo