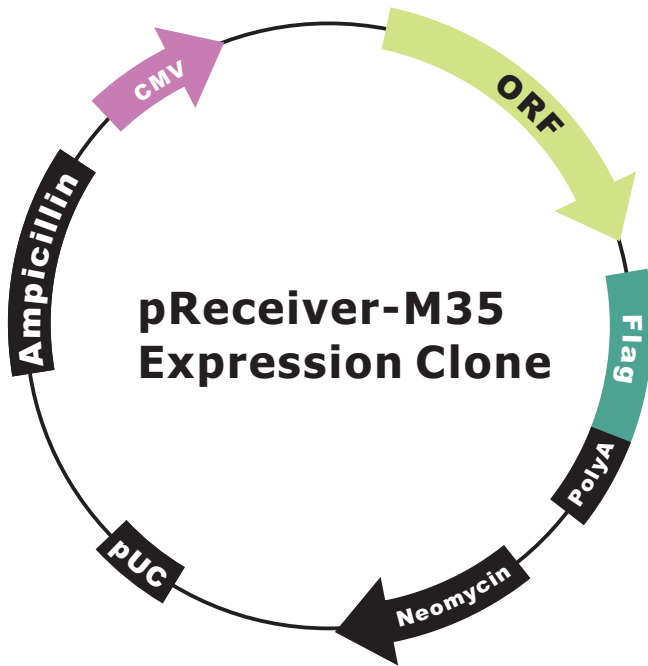


# OmicLink™ Expression Clone (CMV Promoter)



## Vector Features

Promoter	CMV
Host Cell	Mammalian
Bacterial selection antibiotic	Ampicillin
Mammalian selection marker	Yes
Tag	Flag