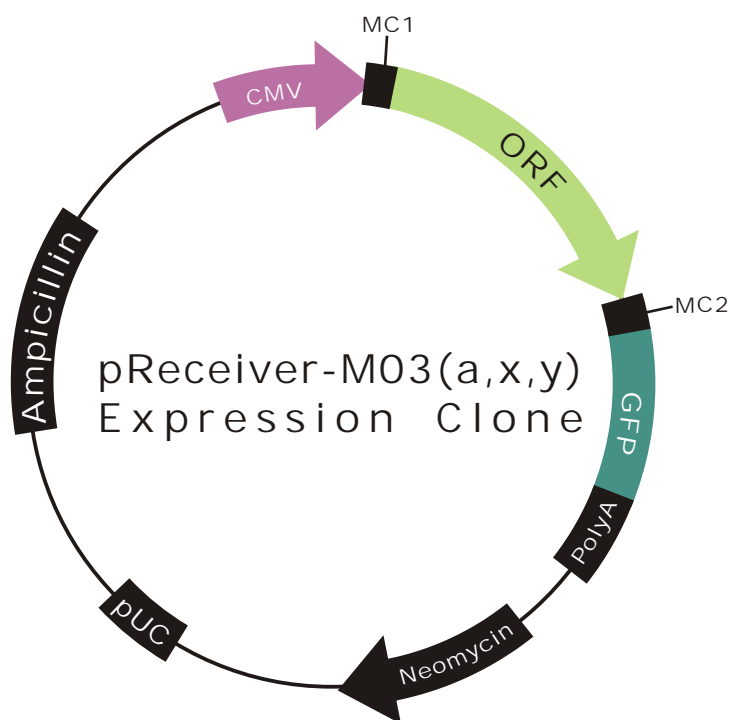


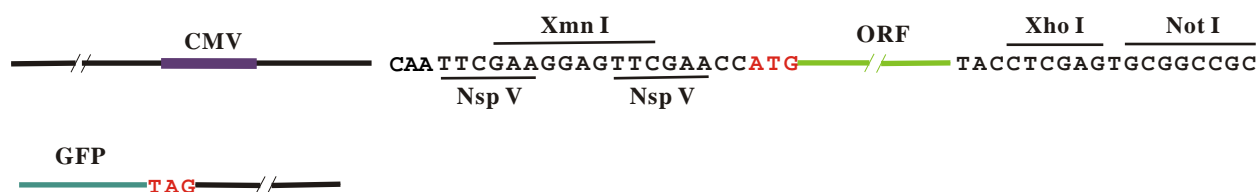
OmicLink™ Expression Clone (CMV Promoter)



Vector Features

Promoter	CMV
Host Cell	Mammalian
Bacterial selection antibiotic	Ampicillin
Mammalian selection marker	Yes
Tag	C-GFP

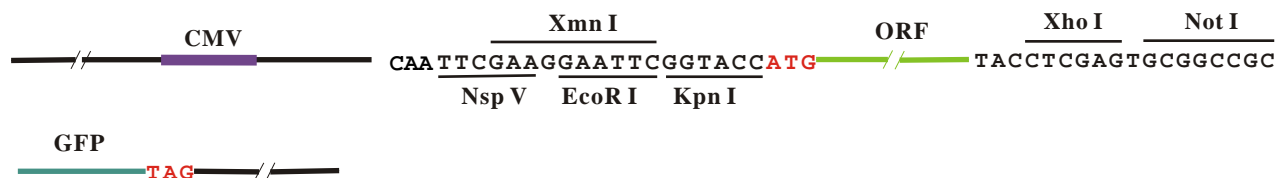
pReceiver-M03a-Xmnl



pReceiver-M03a-NspV



pReceiver-M03x



pReceiver-M03y

