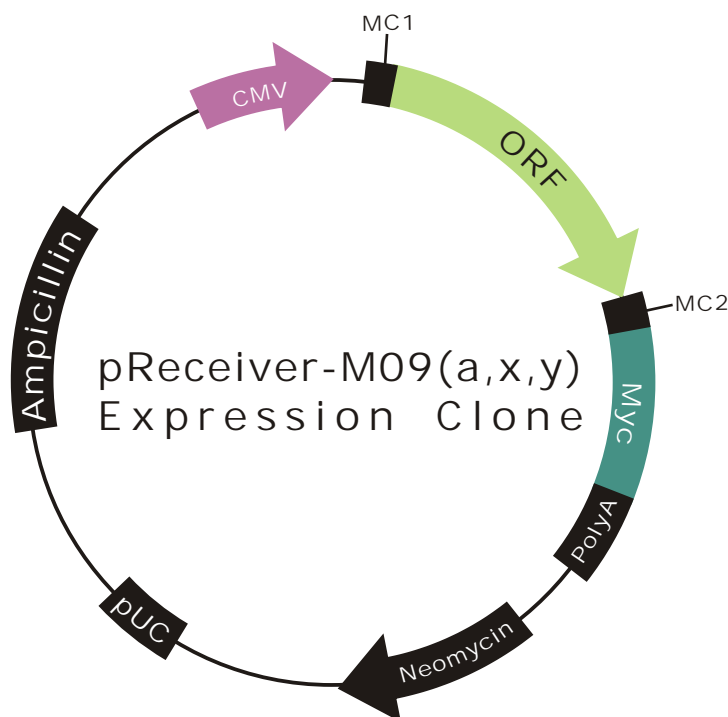


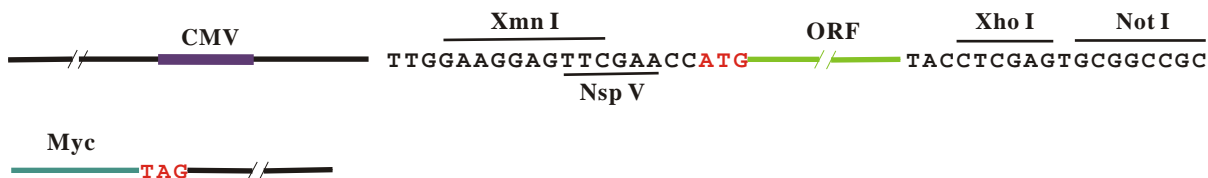
# OmicLink™ Expression Clone (CMV Promoter)



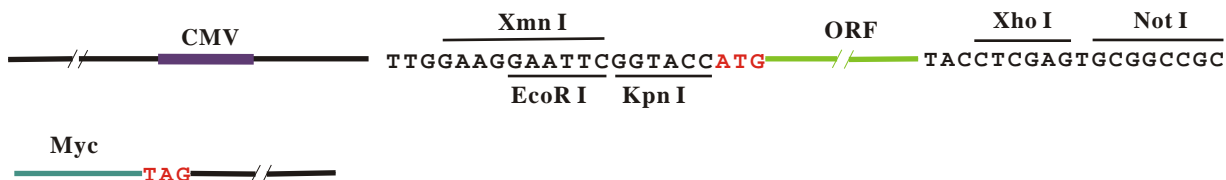
## Vector Features

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

## pReceiver-M09a



## pReceiver-M09x



## pReceiver-M09y

