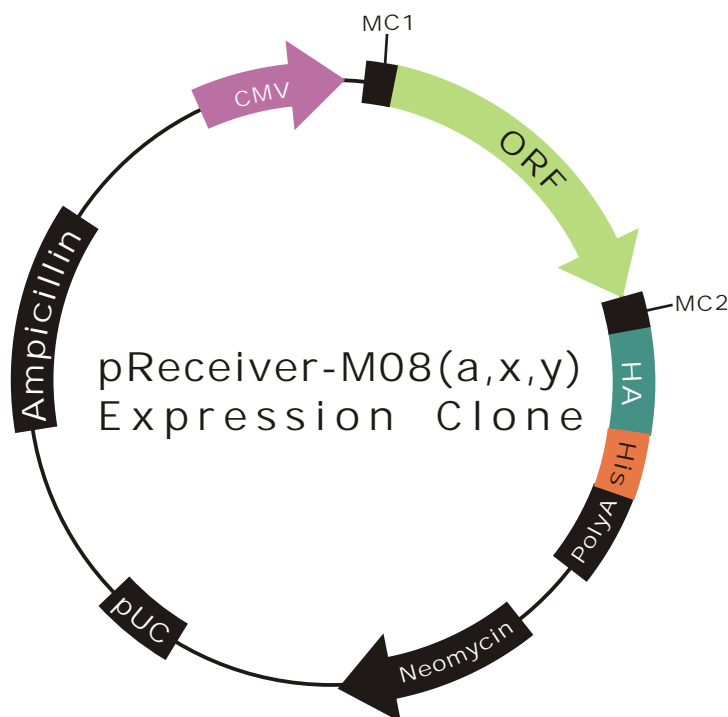


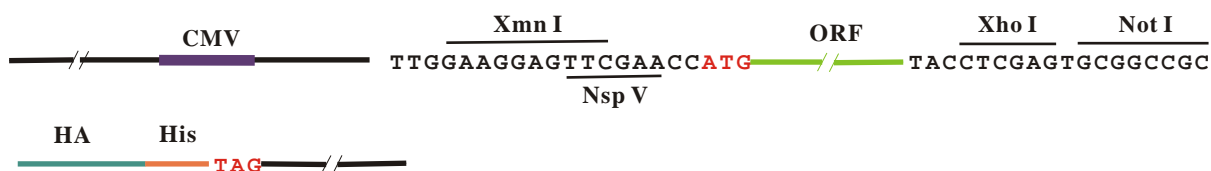
# OmicLink™ Expression Clone (CMV Promoter)



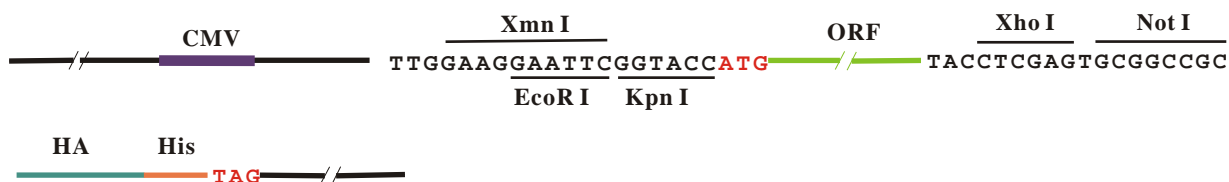
## Vector Features

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

## pReceiver-M08a



## pReceiver-M08x



## pReceiver-M08y

