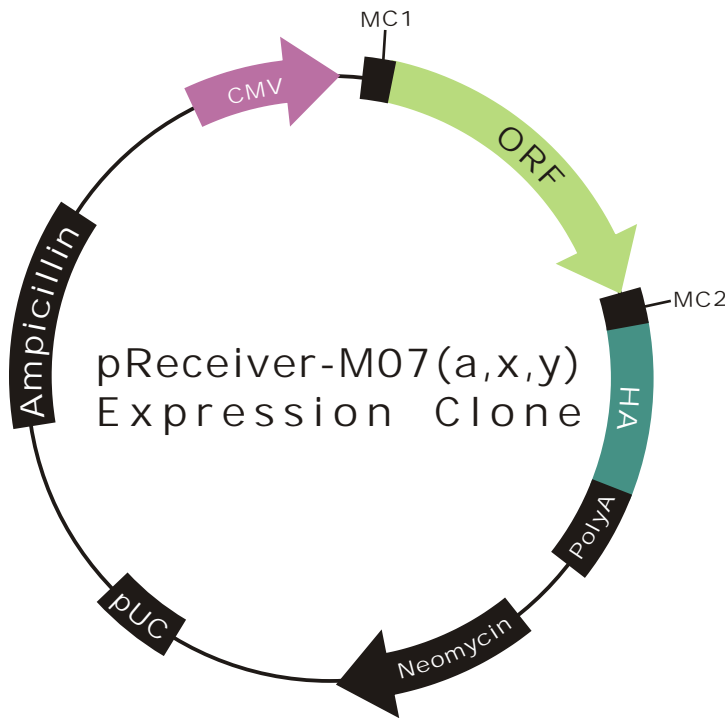


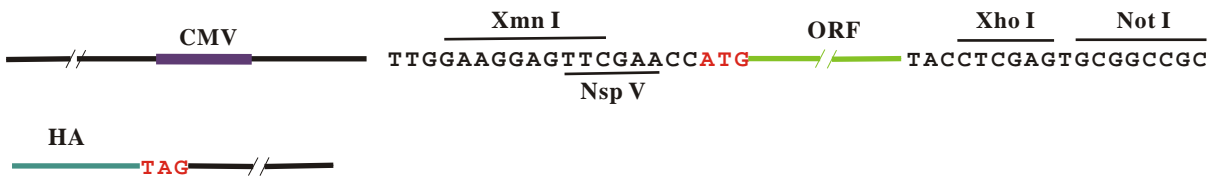
OmicLink™ Expression Clone (CMV Promoter)



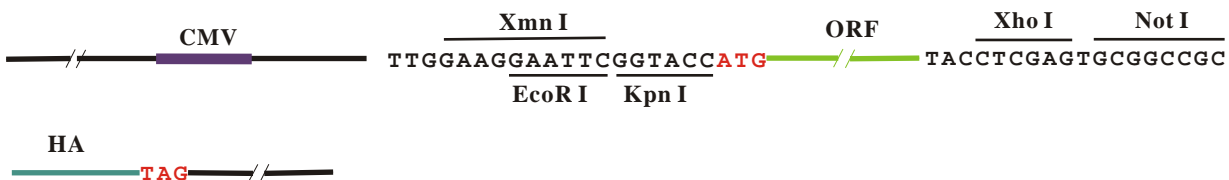
Vector Features

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

pReceiver-M07a



pReceiver-M07x



pReceiver-M07y

