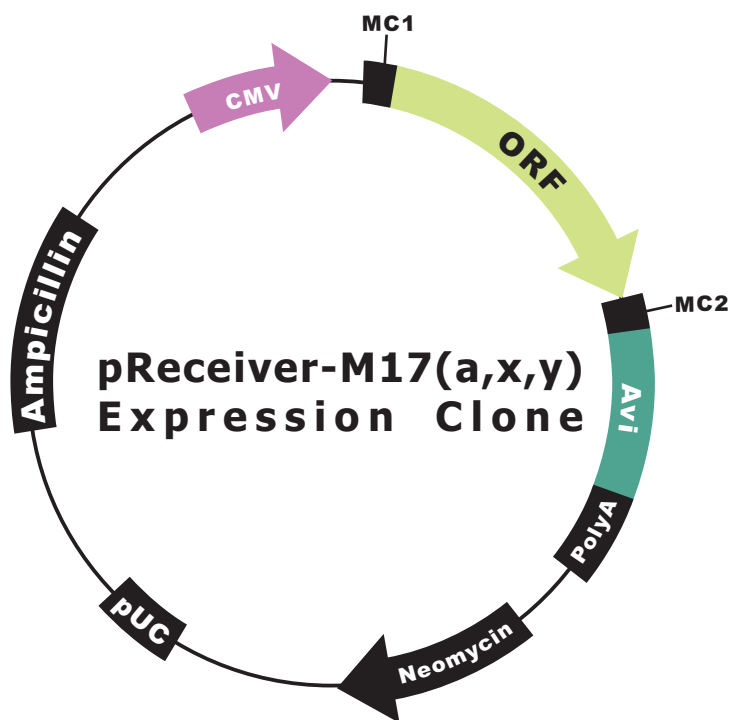


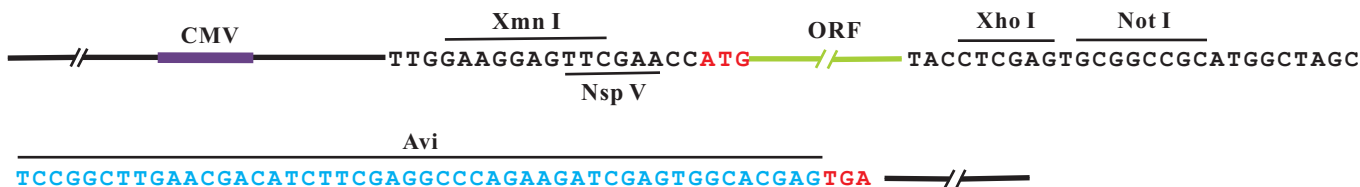
OmicLink™ Expression Clone (CMV Promoter)



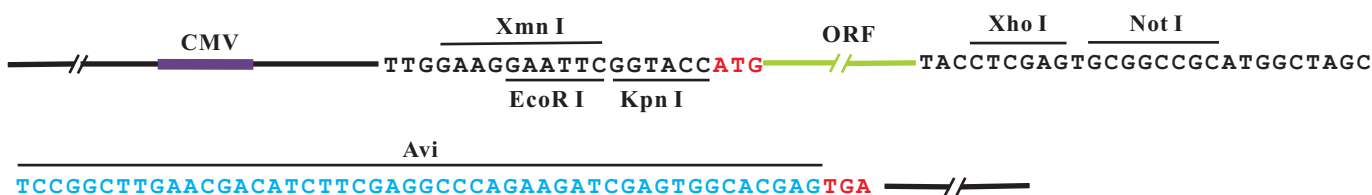
Vector Features

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

pReceiver-M17a



pReceiver-M17x



pReceiver-M17y

