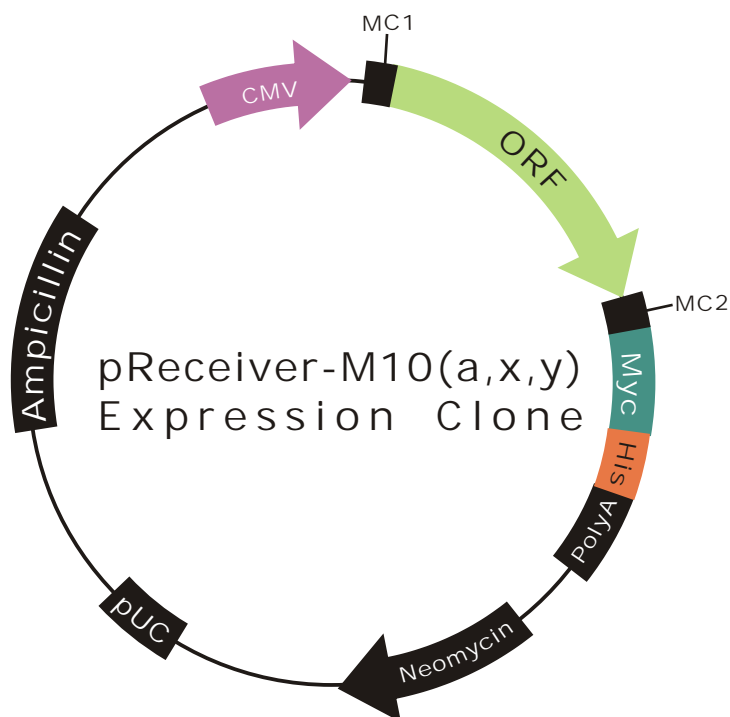


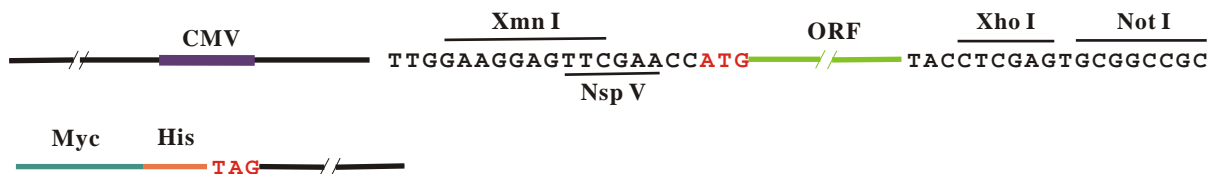
OmicLink™ Expression Clone (CMV Promoter)



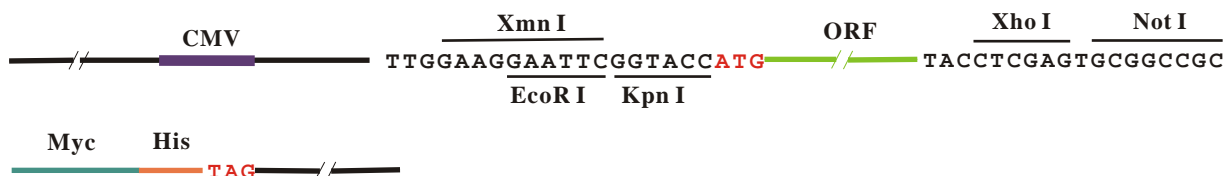
Vector Features

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

pReceiver-M10a



pReceiver-M10x



pReceiver-M10y

