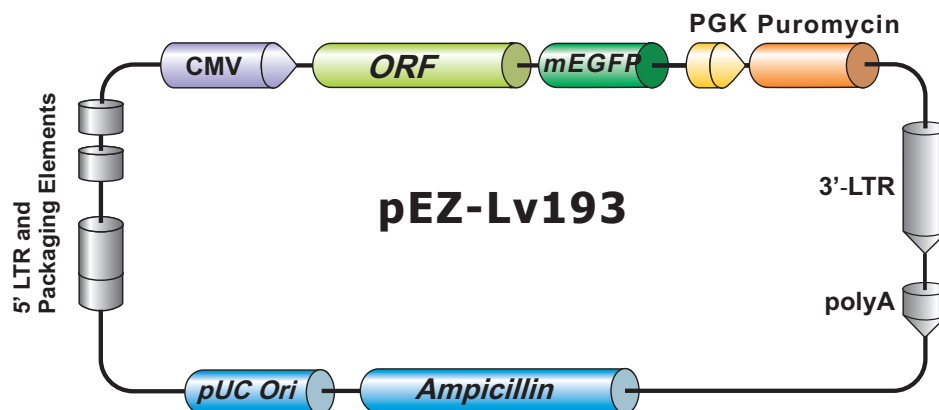
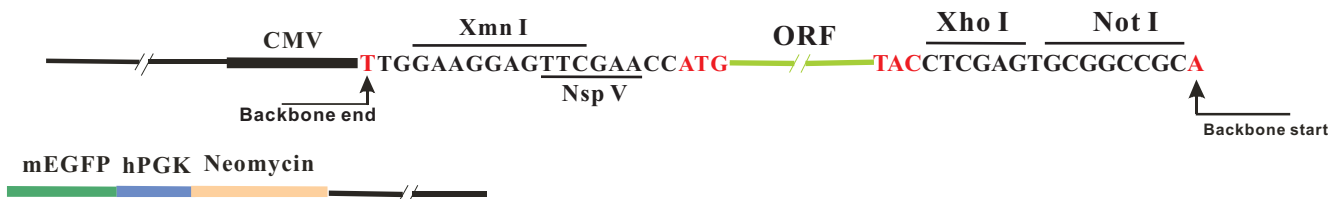


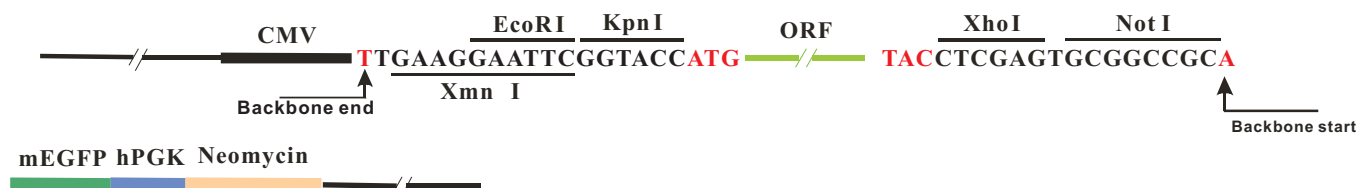
OmicLink™ Expression Clone (CMV Promoter)



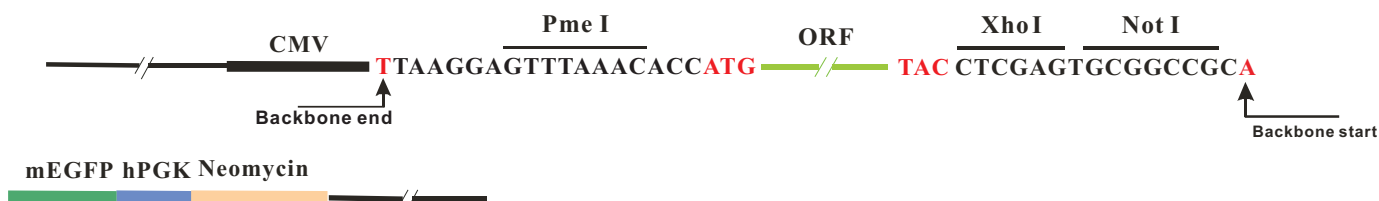
pEZ-Lv193a



pEZ-Lv193x

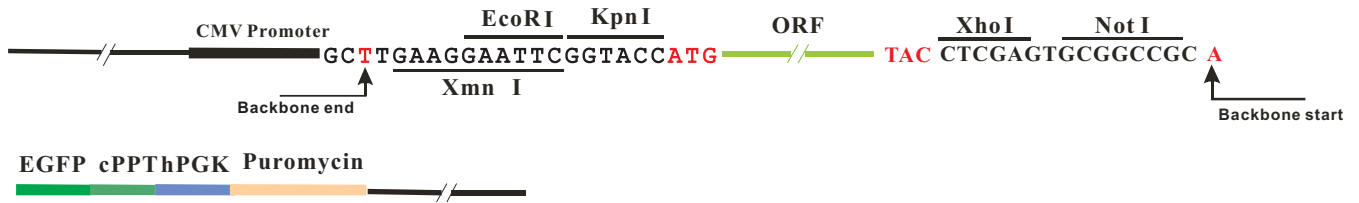


pEZ-Lv193y

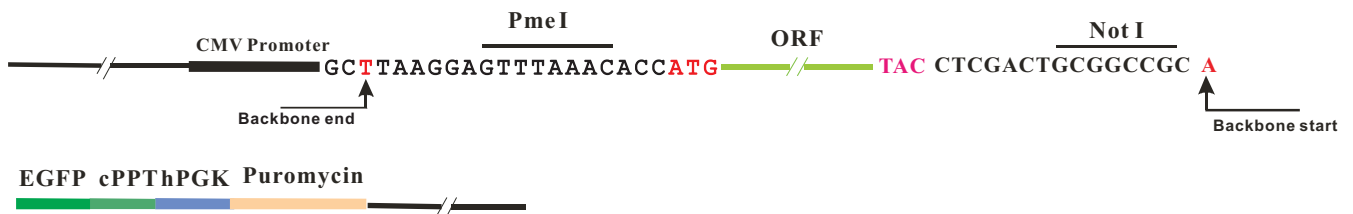


OmicLink™ Expression Clone (CMV Promoter)

pReceiver-Lv122x-NotI



pReceiver-Lv122y-XhoI



pReceiver-Lv122y-NotI

