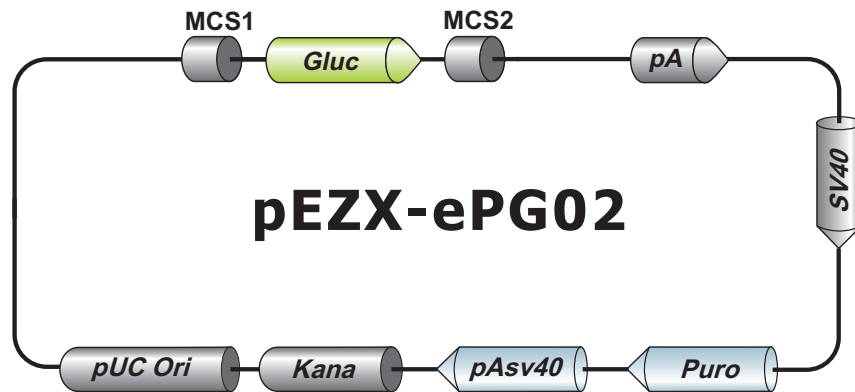


GLuc-ON™ Promoter Reporter Clones



MCS1: BstBI, BglII, EcoRI, EcoRV, HindIII
MCS2: NotI, XbaI