



MCS1: AfeI, AgeI, PmeI, EcoRI, MfeI, BstBI

MCS2: XhoI, PspXI