



**MCS1:** BstBI, MfeI, EcoRI, PmeI, AgeI, AfeI  
**MCS2:** PspXI, XhoI