



TR002

MCS: SpeI, SwaI