

Structure of Rous Sarcoma Virus Genome

