

Binding kinetics of BMS-777607 on BTN-MER

