

Binding kinetics of BMS-777607 on TYRO3 non-activated

