

**Binding kinetics of BMS-777607 on BTN-TYRO3[non-activated]**

