

Binding kinetics of BMS-777607 on EPHA2

