

Binding kinetics of SB 431542 on TGFBR1

