

Binding kinetics of ponatinib on TRKB (unactivated)

