

Binding kinetics of ponatinib on TRKB (NTRK2) non-activated

