

Binding kinetics of AZ23 on TRKC (NTRK3)

