

Binding kinetics of AZ23 on TRKA (NTRK1) non-activated

