

Binding kinetics of ponatinib on TRKA (unactivated)

