

Binding kinetics of AZ23 on TRKA

