

# Binding kinetics of AZ23 on TRKB

TRKB AZ23

