

# Binding kinetics of SNS-032 on CDK5/p25

