

# Binding kinetics of ARQ-531 on BTK [T474I]

