

## Binding kinetics of ARQ-531 on BTK [T474S]

