

# Binding kinetics of acalabrutinib on BTK[T316A]

BTK[T316A] Acalabrutinib  
RU

