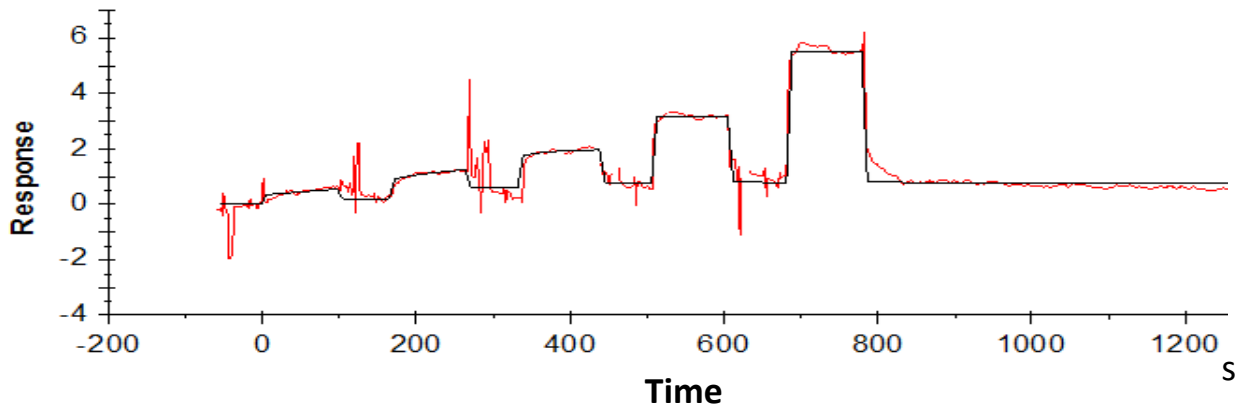


# Binding kinetics of ibrutinib on BTK[T474I] non activated

BTK[T474I] non activated Ibrutinib  
RU



The data was generated and provided by  
**Netherlands Translational Research Center B.V. (NTRC)**