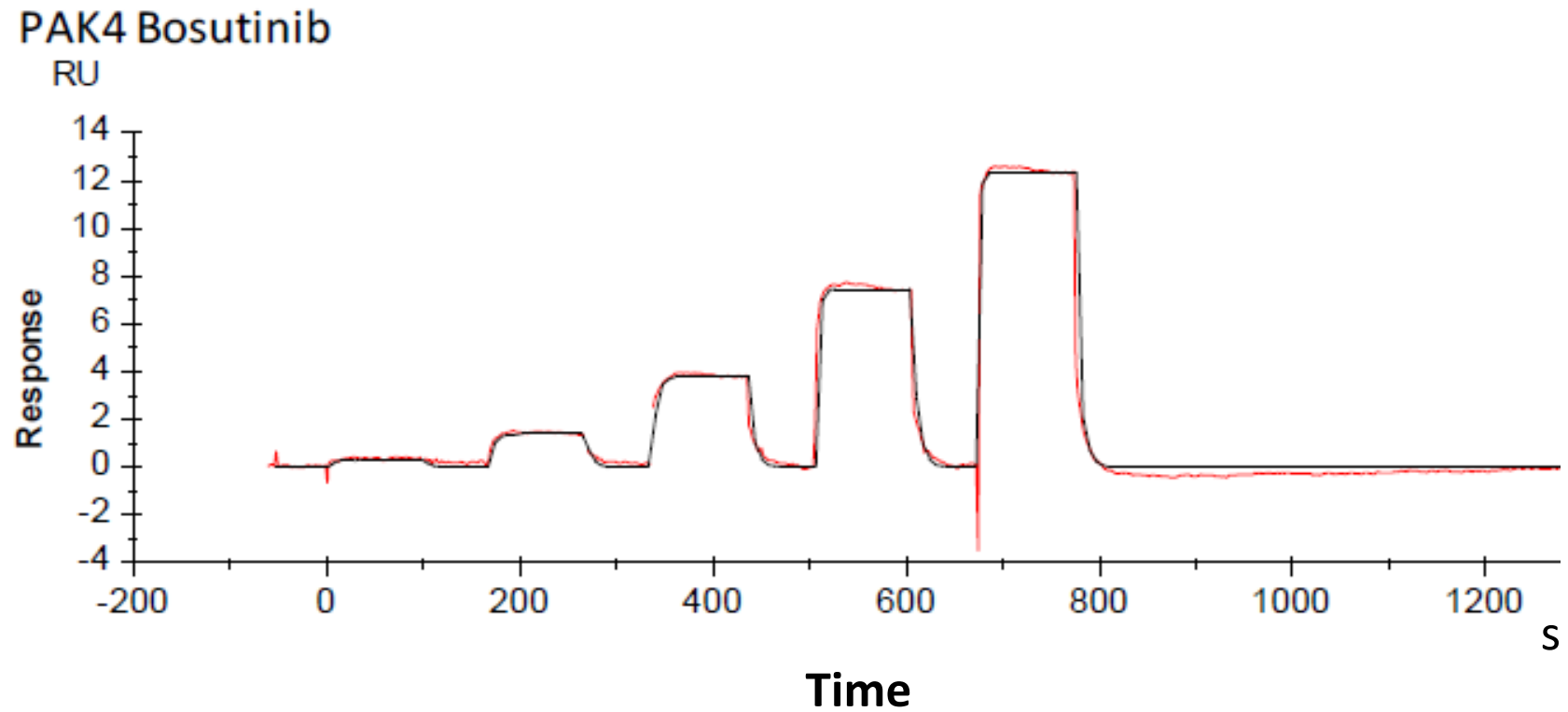


# Binding kinetics of bosutinib on PAK4



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