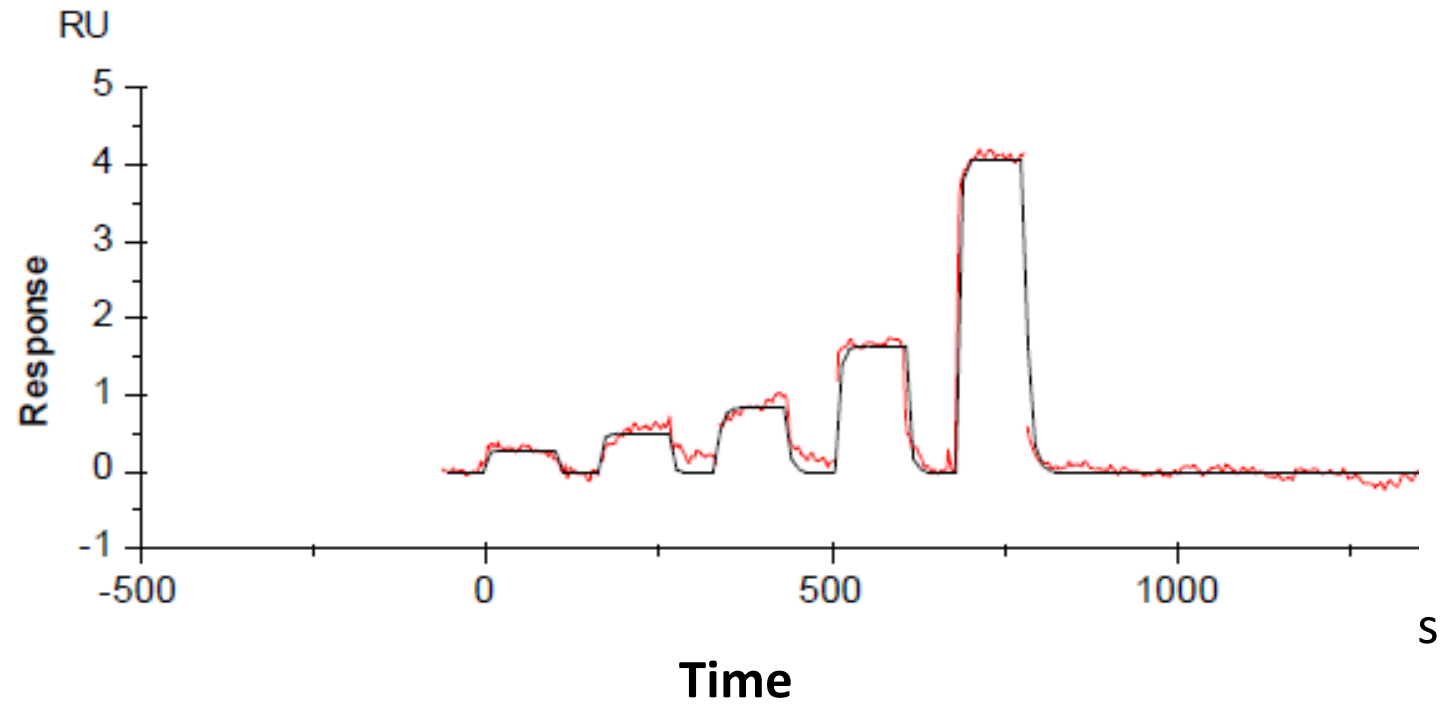


# Binding kinetics of dasatinib on TGFBR1

TGFBR1 Dasatinib



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