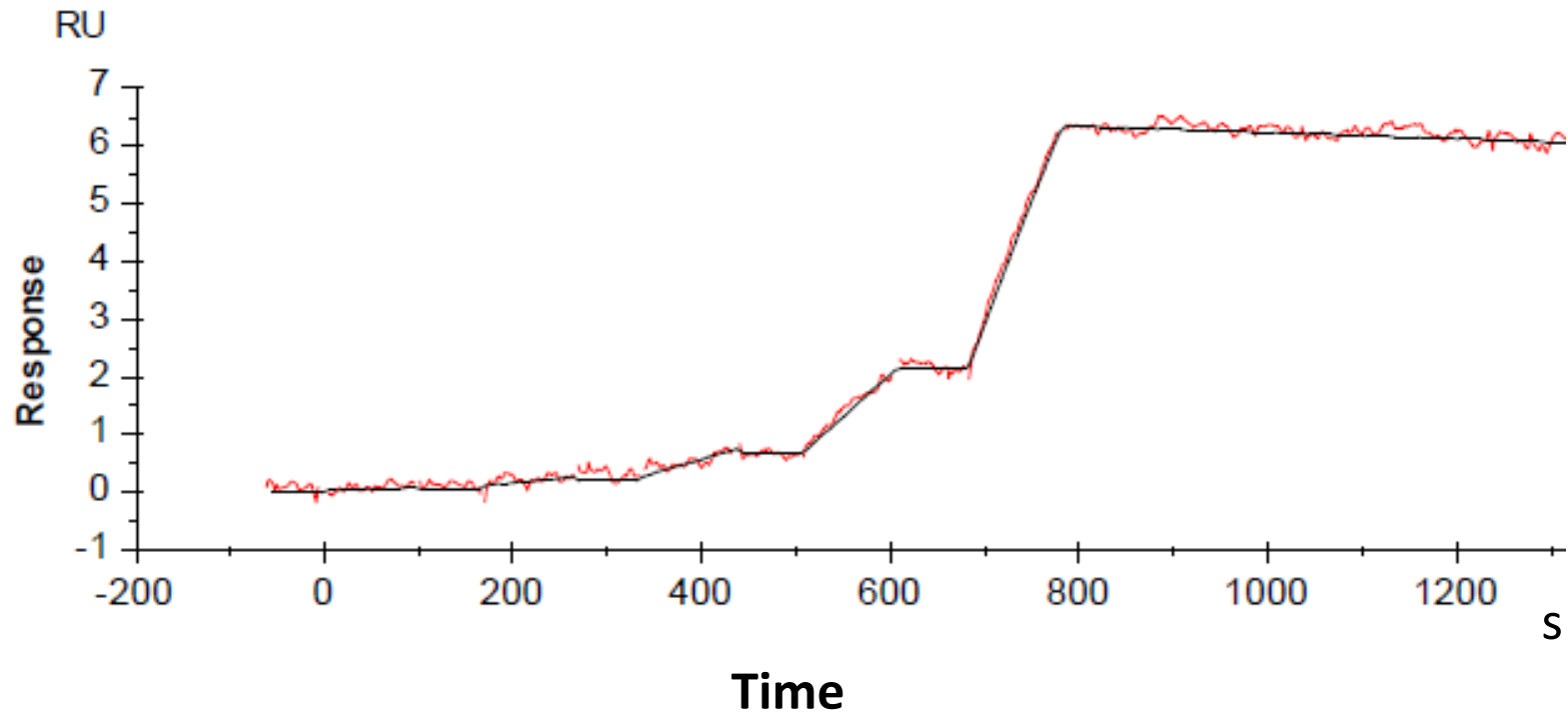


# Binding kinetics of GSK 1070916 on AurB /INCENP

GSK 1070916 on AURB /INCENP



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