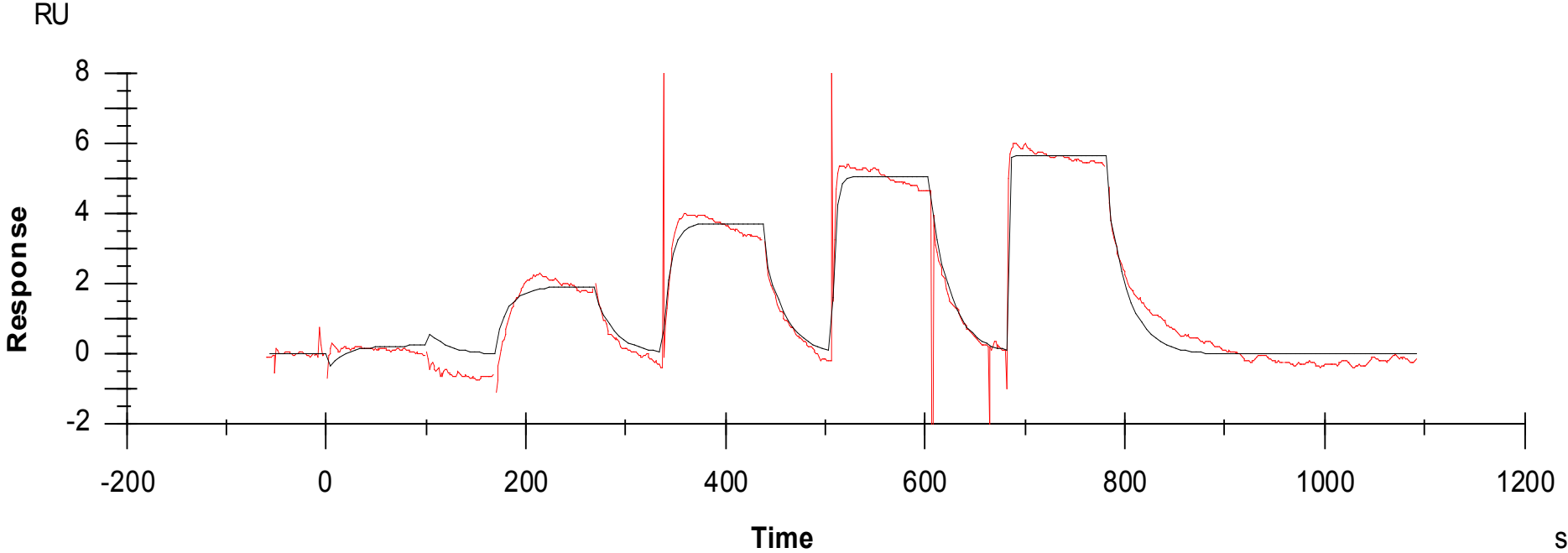


Binding kinetics of saracatinib on YES



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