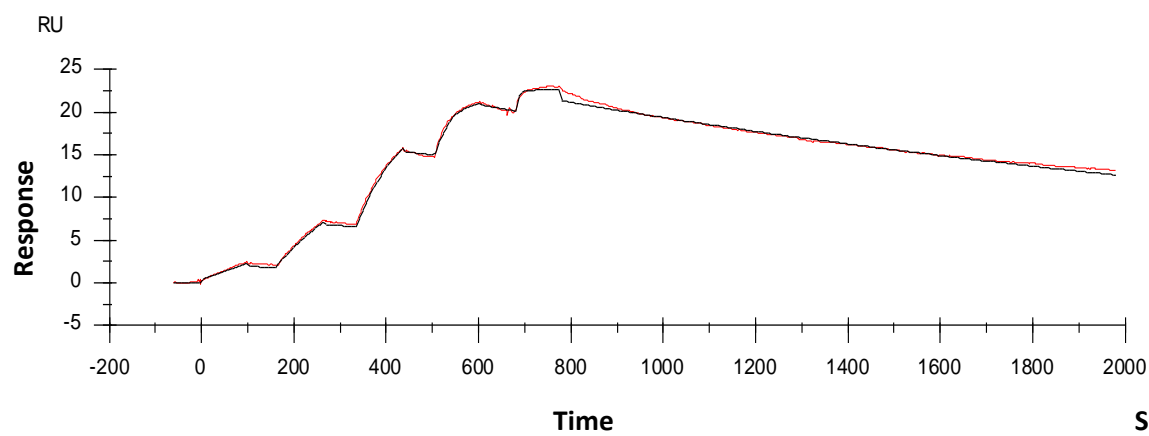


Binding kinetics of Ibrutinib on BTN-BRK(non-activated)

BTN-BRK Ibrutinib



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