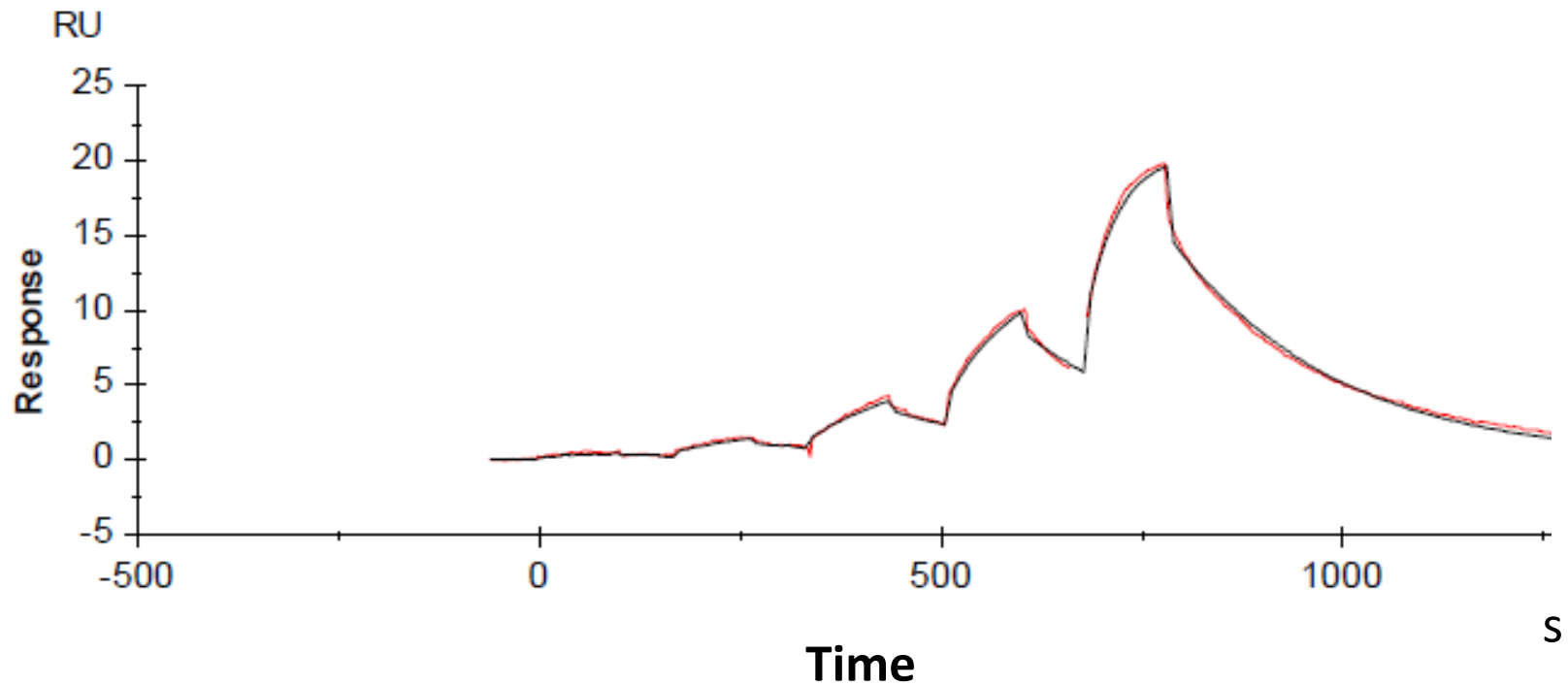


Binding kinetics of dasatinib on FLT3 (non-activated)

FLT3 Dasatinib



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