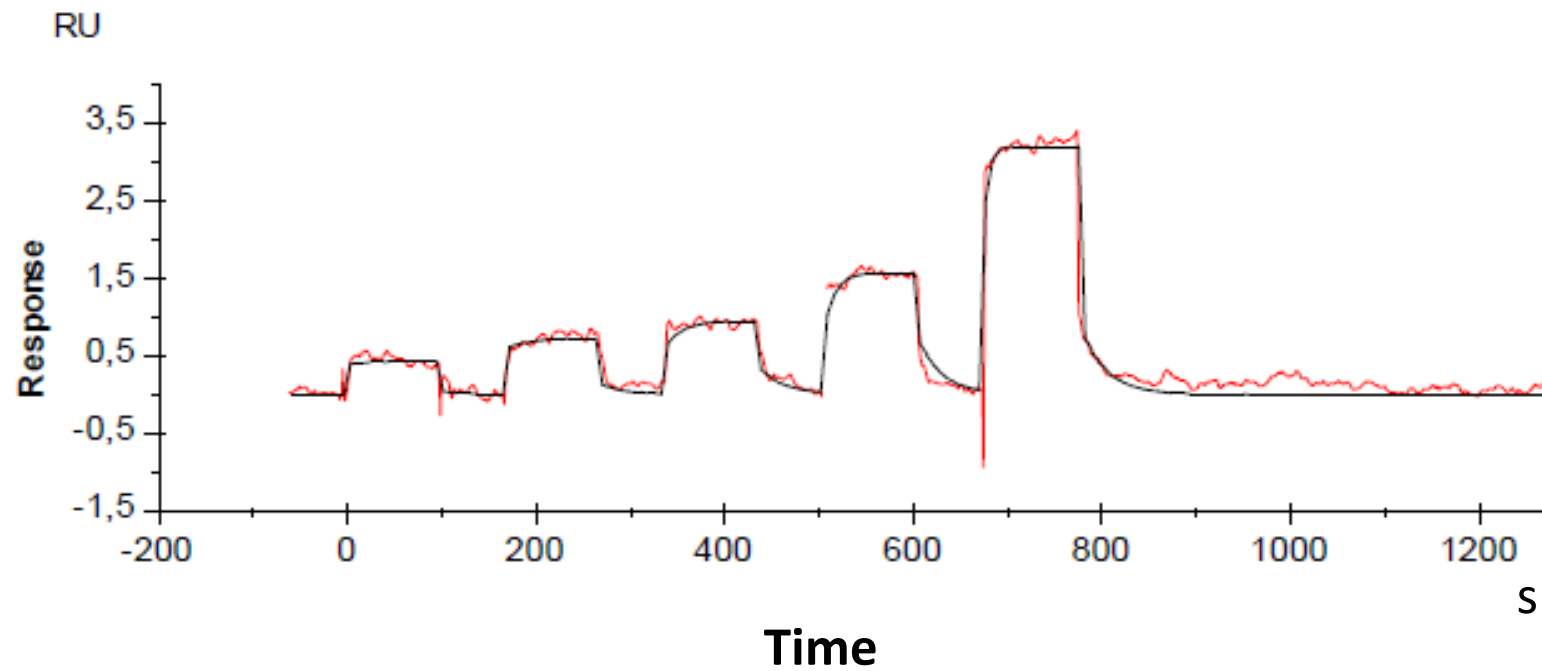


Binding kinetics of vandetanib on KDR

KDR-Vandetanib



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