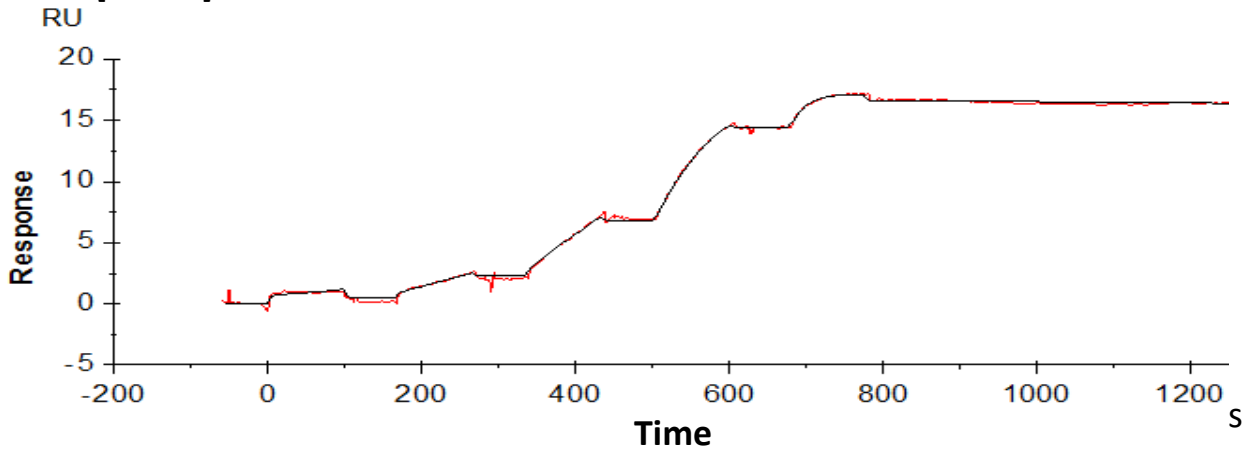


Binding kinetics of GDC 0853-0 on BTK[C481S] non activated

BTK[C481S] non activated GDC 0853-0



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