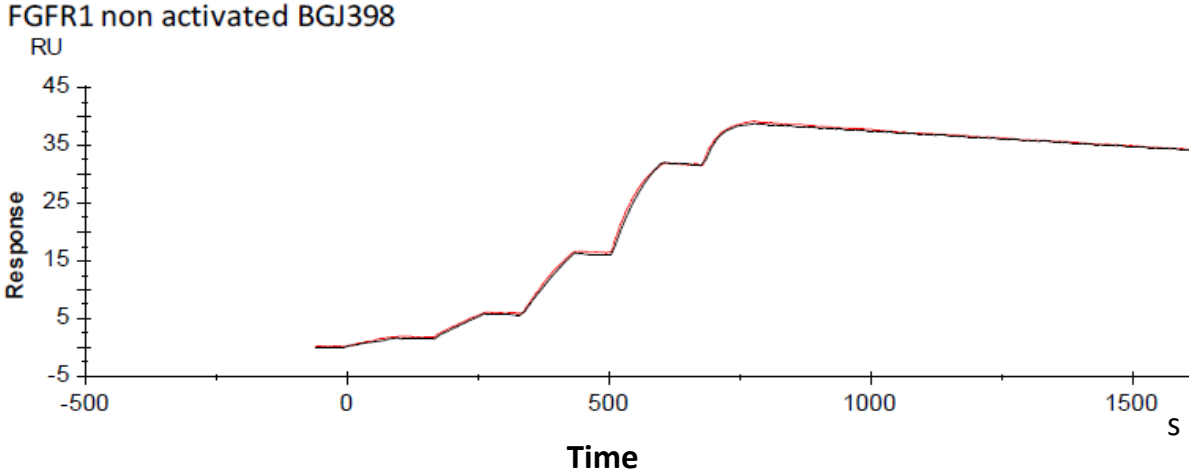


Binding kinetics of BGJ398 on FGFR1 non activated



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