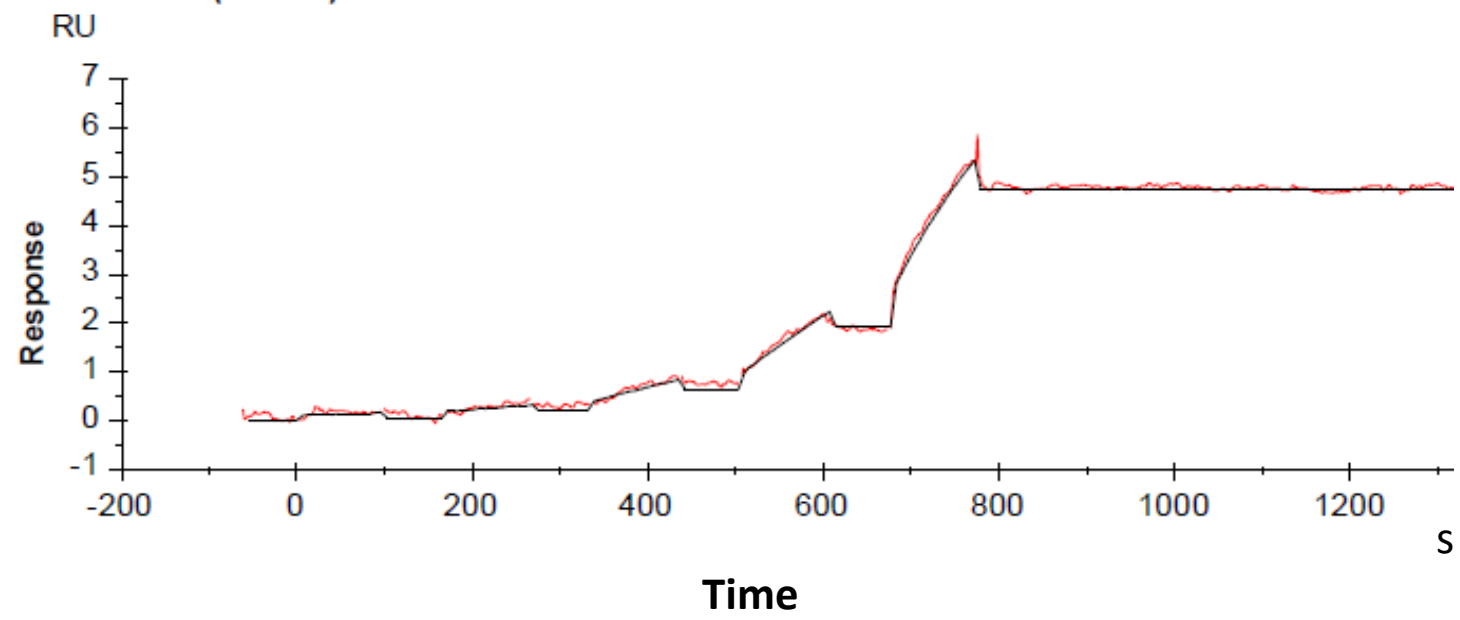


Binding kinetics of NTRC 0799-0 on IKK α (CHUK) inactive mutant

BTN-IKK α (CHUK) inactive mutant NTRC 0799-0



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