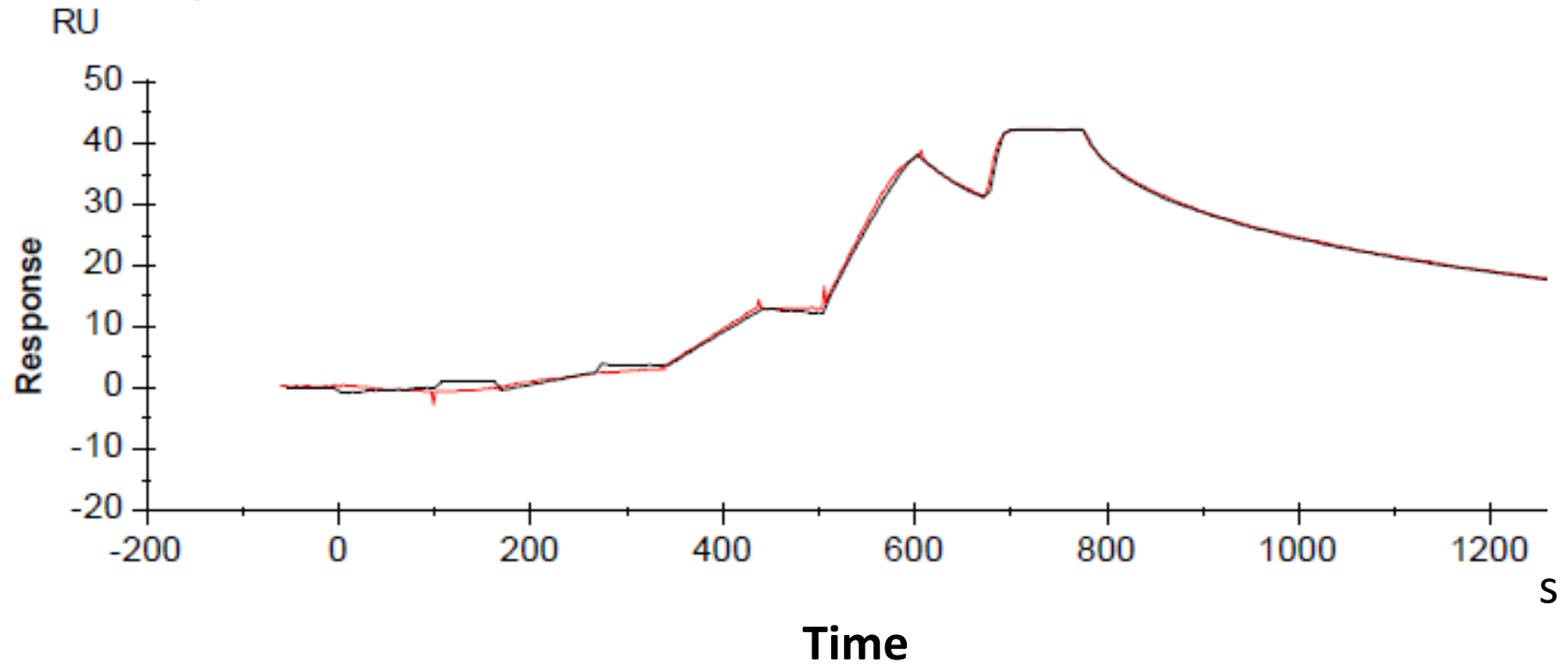


Binding kinetics of PH797804 on p38 (inactive)

inactive p38 PH797804



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