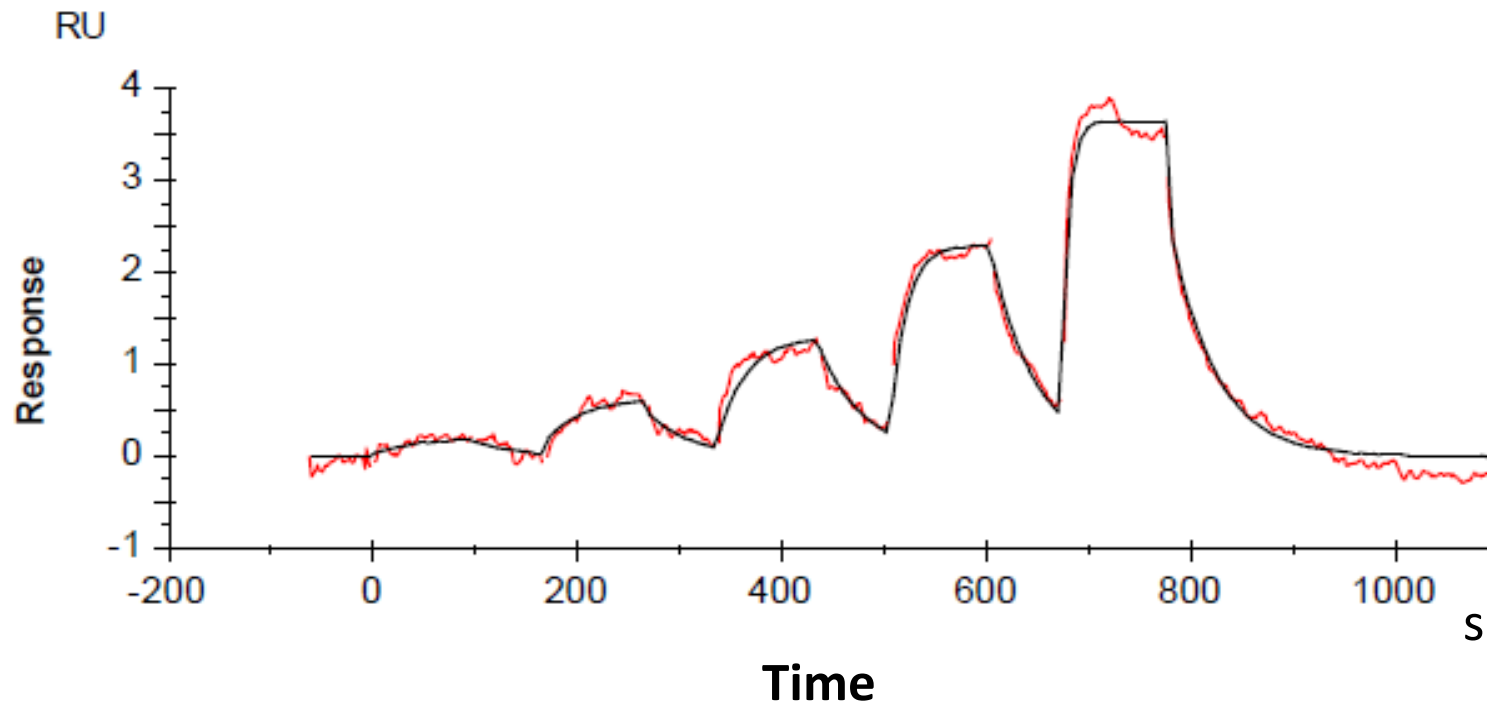


Binding kinetics of sorafenib on TRKC

TRKC Sorafenib



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