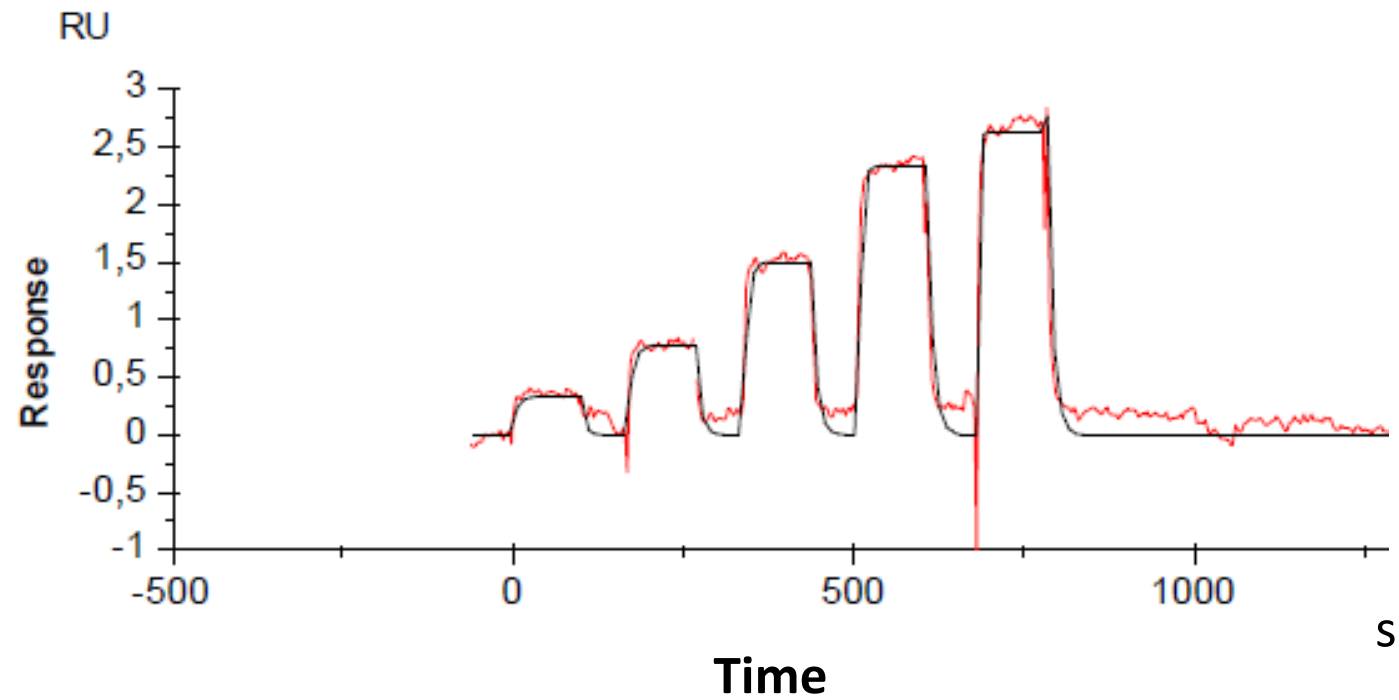


Binding kinetics of sunitinib on FGFR3

FGFR3 - Sunitinib



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