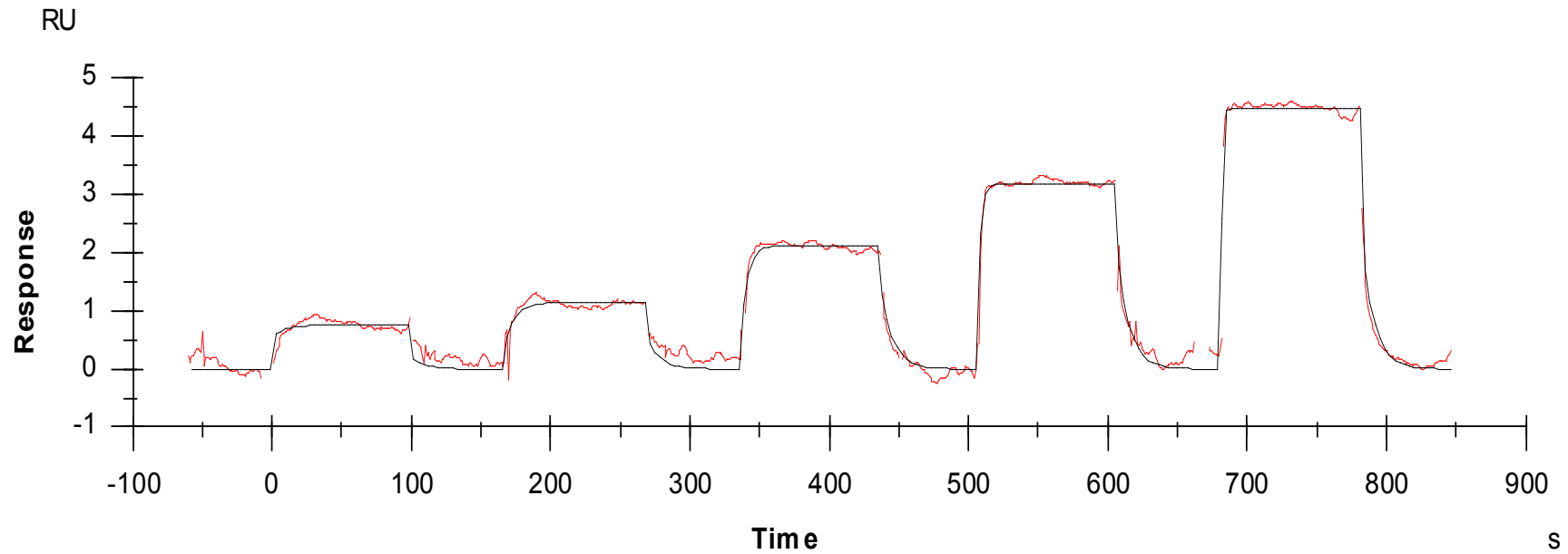


Binding Kinetics of PIK3CA[P539R]/PIK3R1

Idelalisib



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