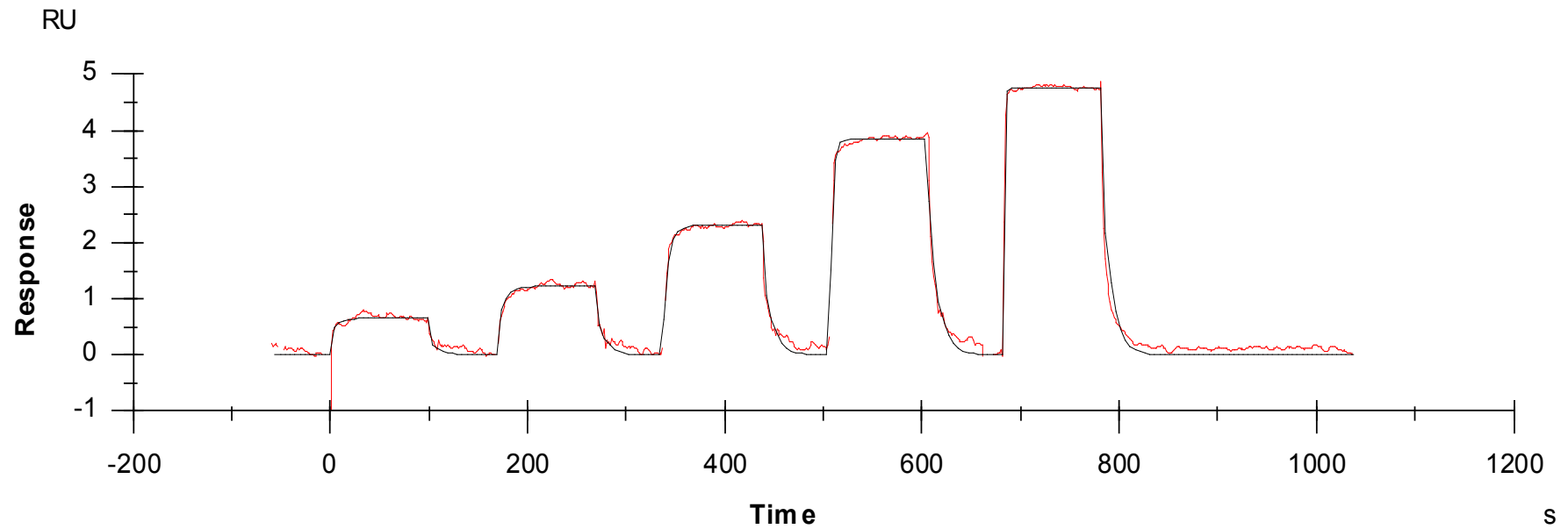


Binding Kinetics of PIK3CA[R88Q]/PIK3R1

Idelalisib



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