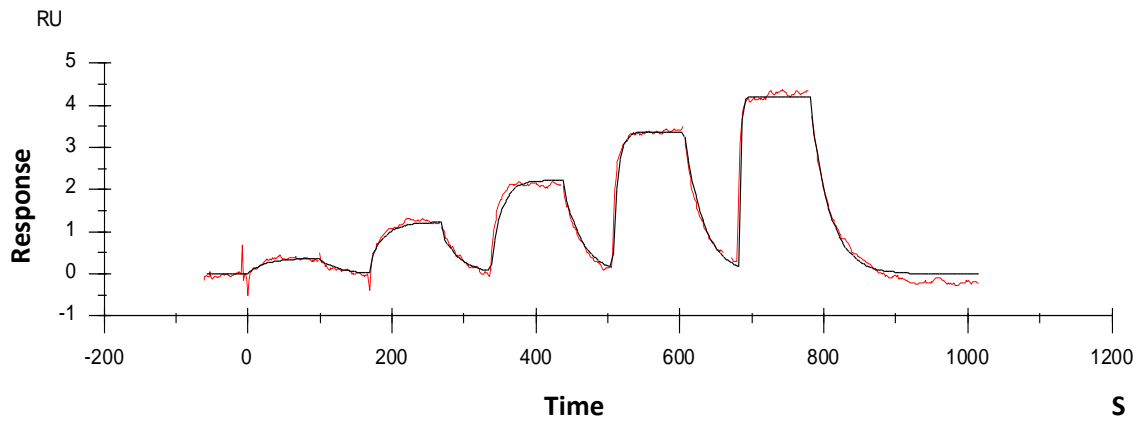


Binding kinetics of Dinaciclib on BTN-CDK4/CycD3

BTN-CDK4/CycD3 Dinaciclib



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