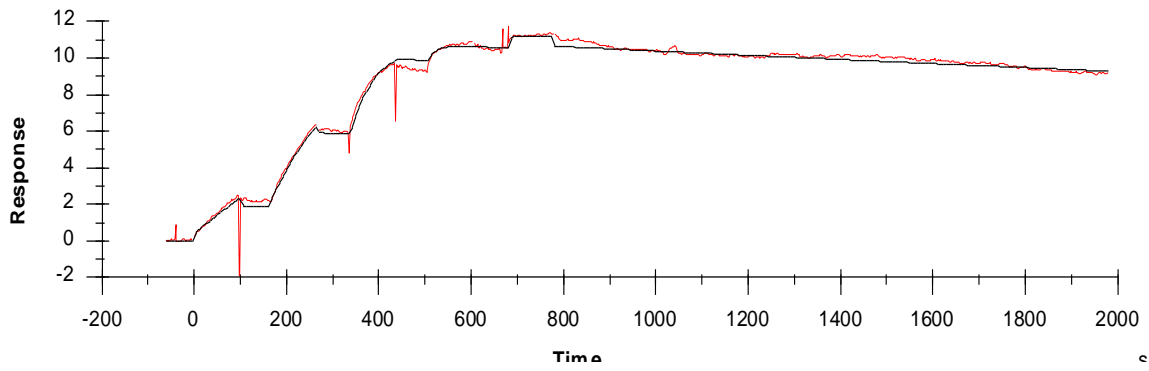


Binding kinetics of dinaciclib on CDK9/CycT1

CDK9/CycT1

dinaciclib

RU



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