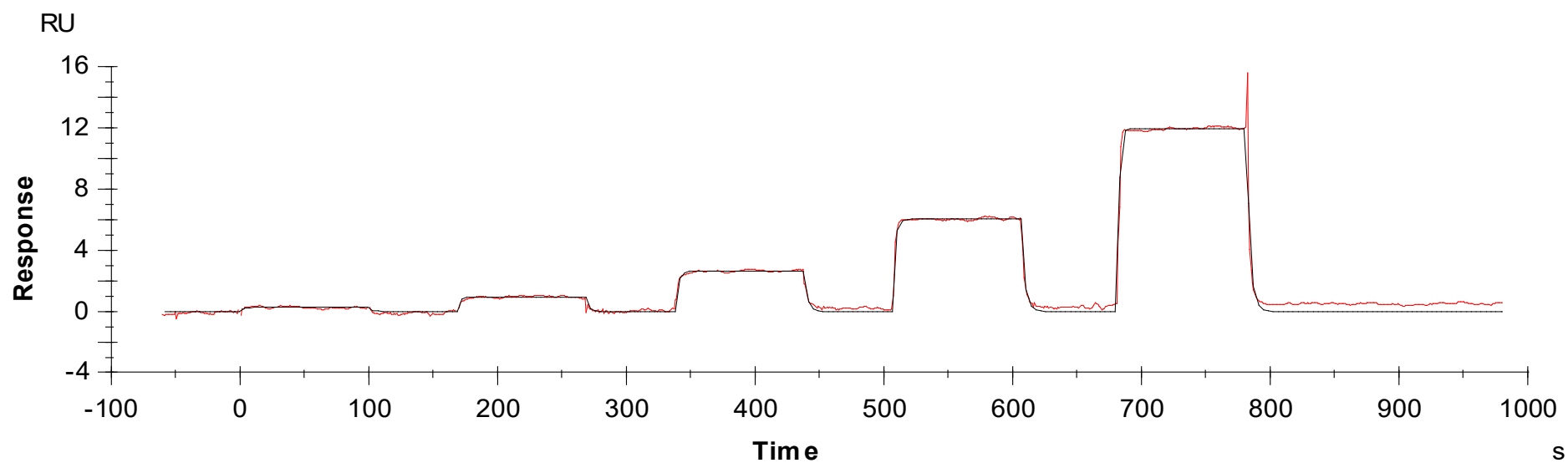




Binding Kinetics of Dinaciclib on CDK1(CDC2)



Data generated and provided by Carna Biosciences, Inc.