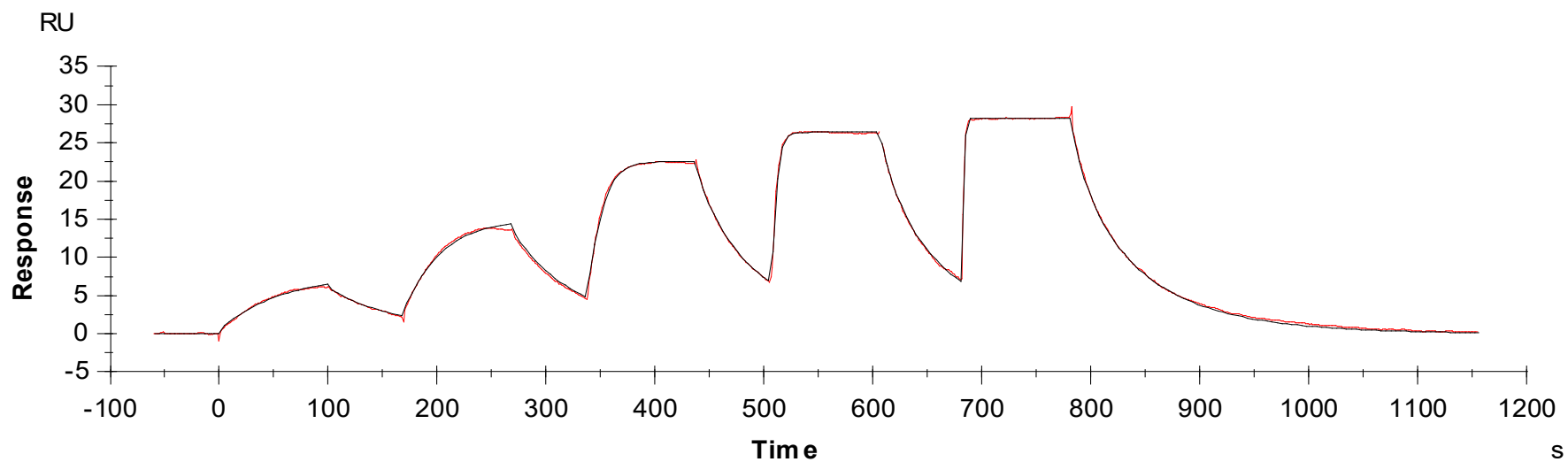




Binding Kinetics of Dinaciclib on CDK2



Data generated and provided by Carna Biosciences, Inc.