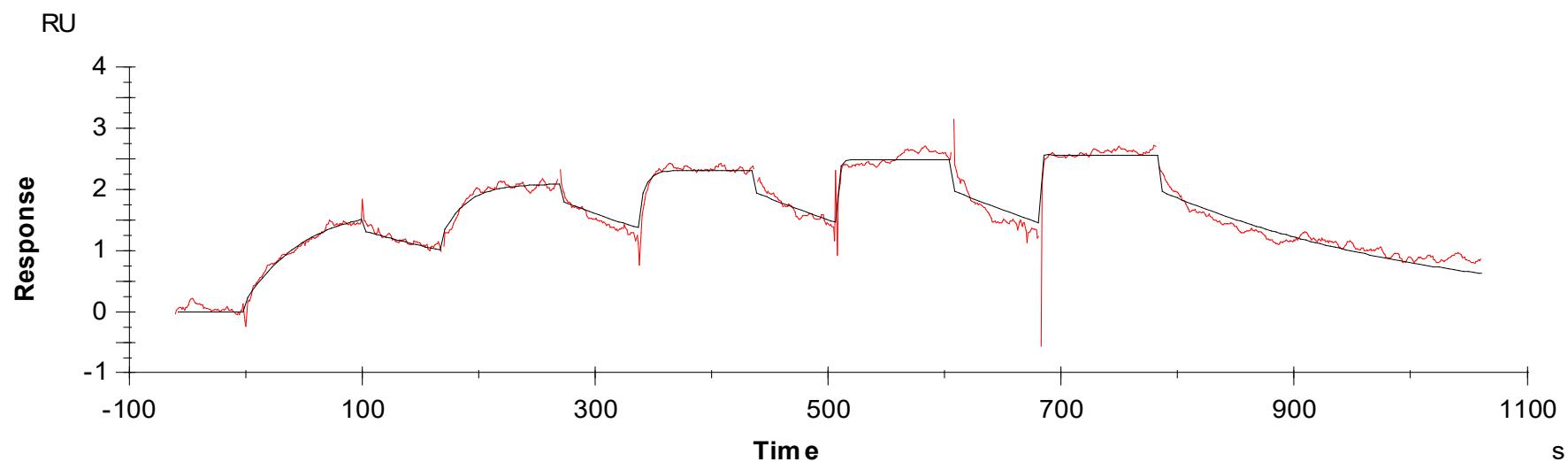




Binding Kinetics of Dinaciclib on CDK5



Data generated and provided by Carna Biosciences, Inc.