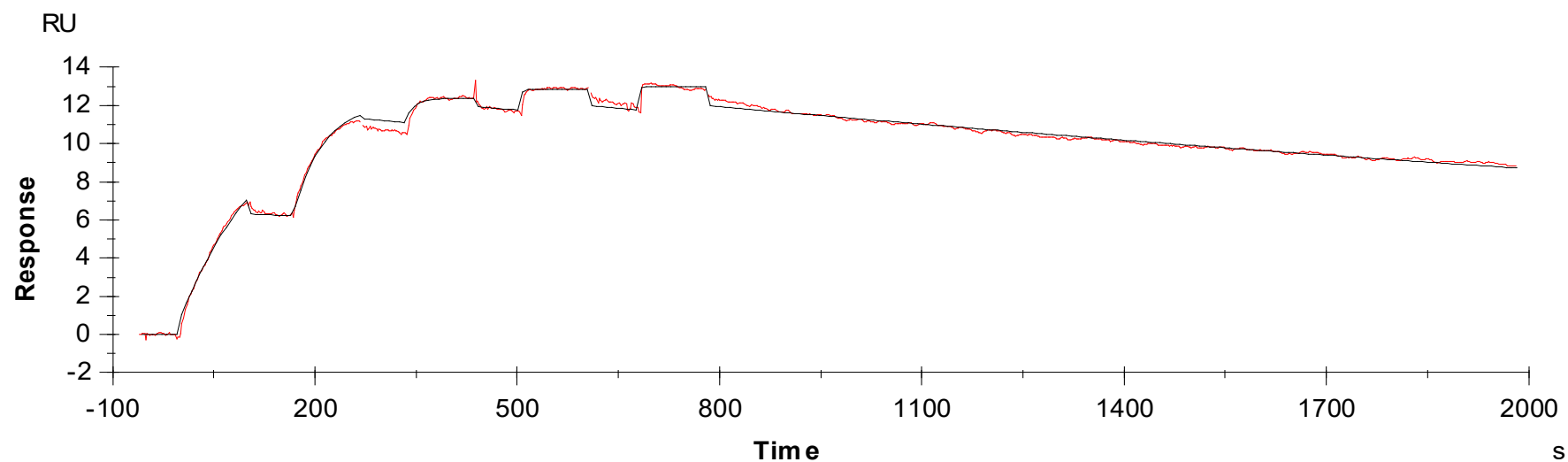




Binding Kinetics of Dinaciclib on CDK5/p35



Data generated and provided by Carna Biosciences, Inc.