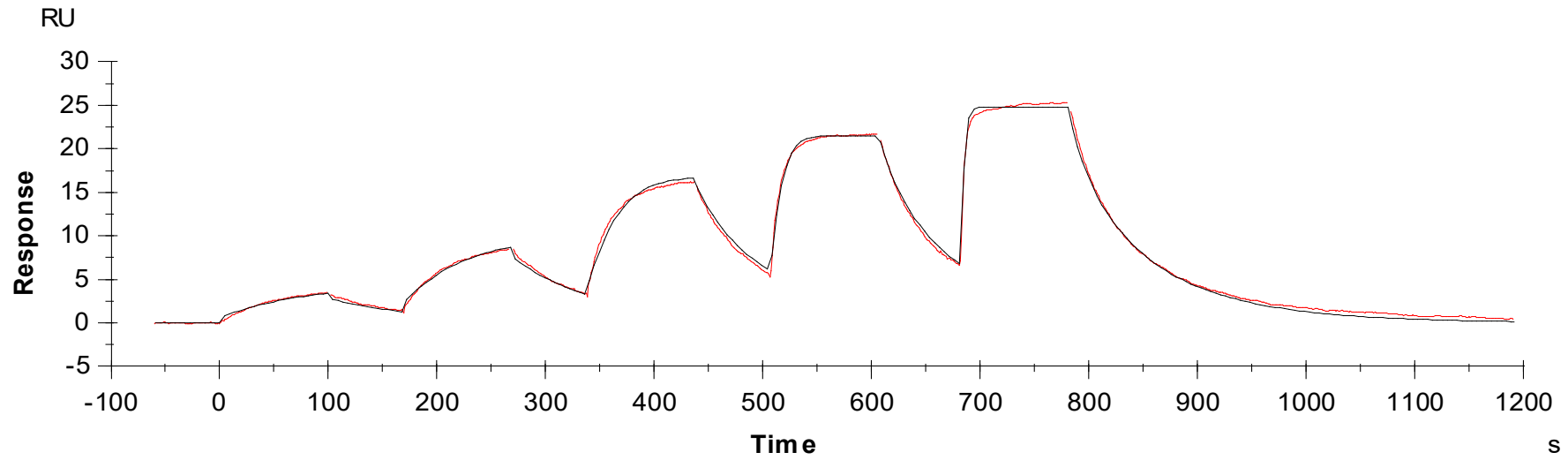




Binding Kinetics of Dinaciclib on CDK6



Data generated and provided by Carna Biosciences, Inc.