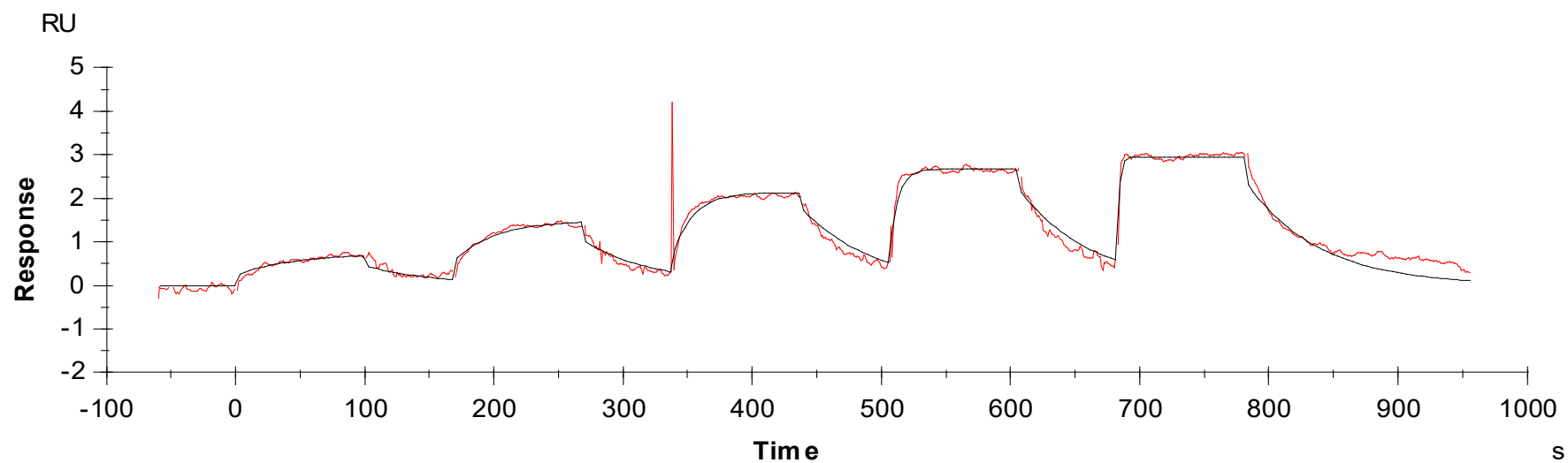


Binding Kinetics of Dinaciclib on CDK3



Data generated and provided by Carna Biosciences, Inc.