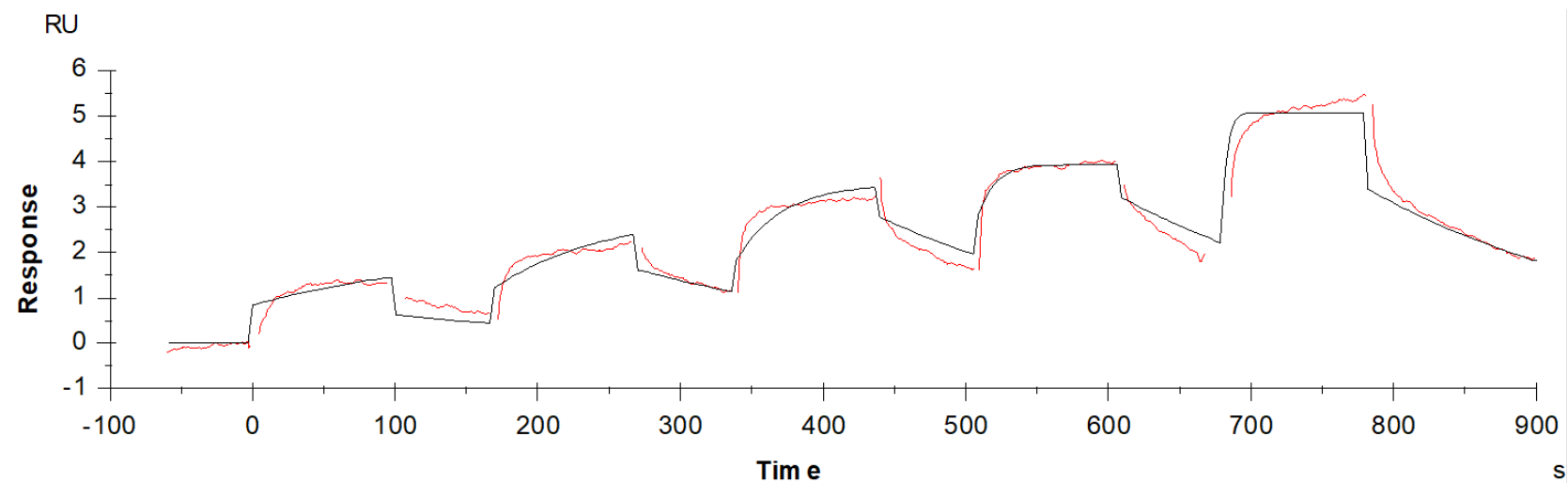


Binding Kinetics of Abemaciclib on CDK4



Data generated and provided by Carna Biosciences, Inc.