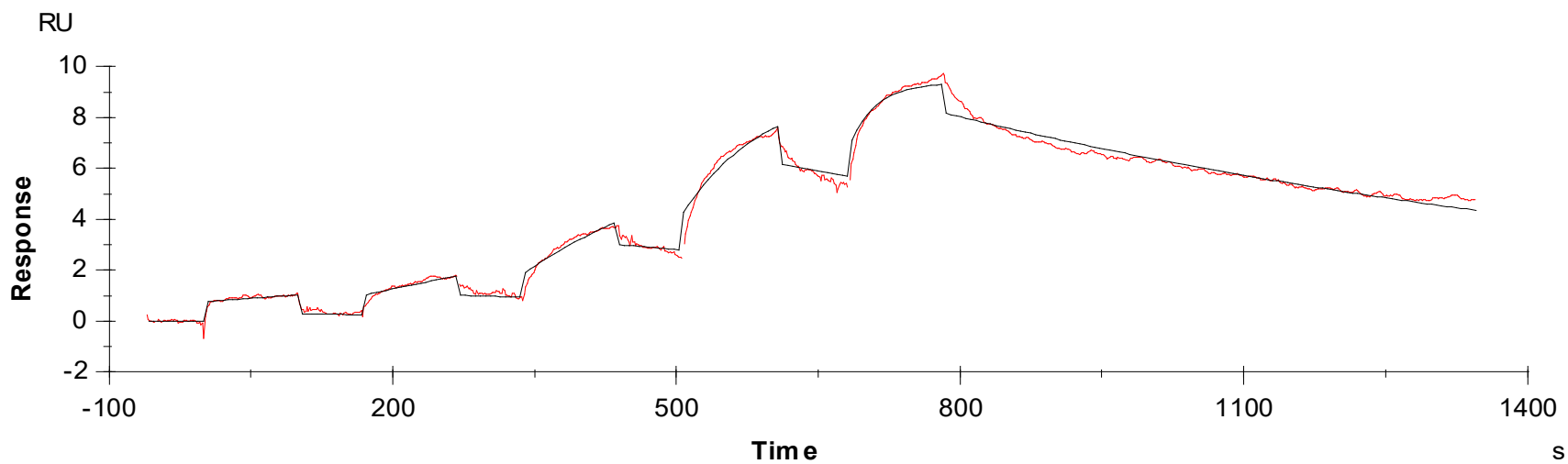




Binding Kinetics of Ponatinib on CDK8



Data generated and provided by Carna Biosciences, Inc.