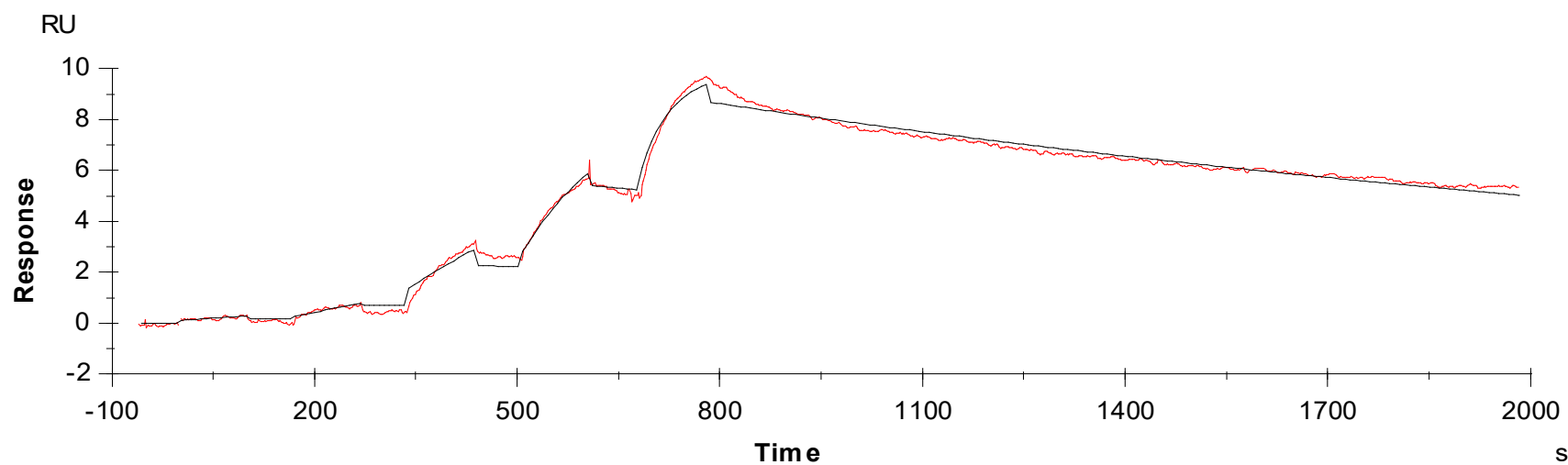




Binding Kinetics of Ponatinib on CDK19(CDC2L6)



Data generated and provided by Carna Biosciences, Inc.