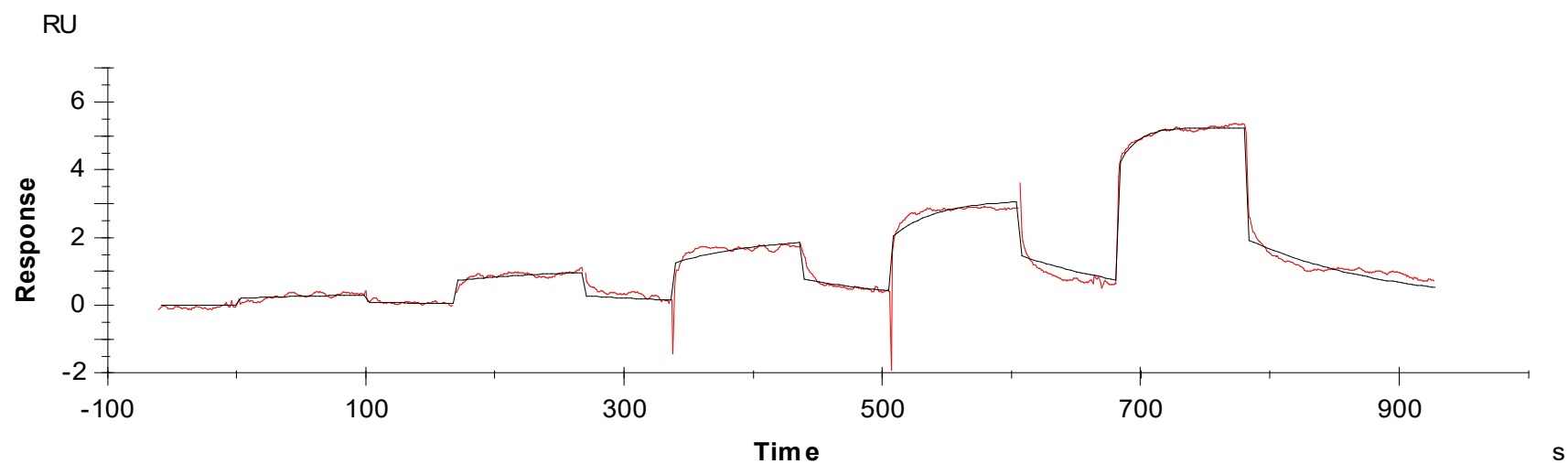




Binding Kinetics of Ibrutinib on BTK[V416L]



Data generated and provided by Carna Biosciences, Inc.