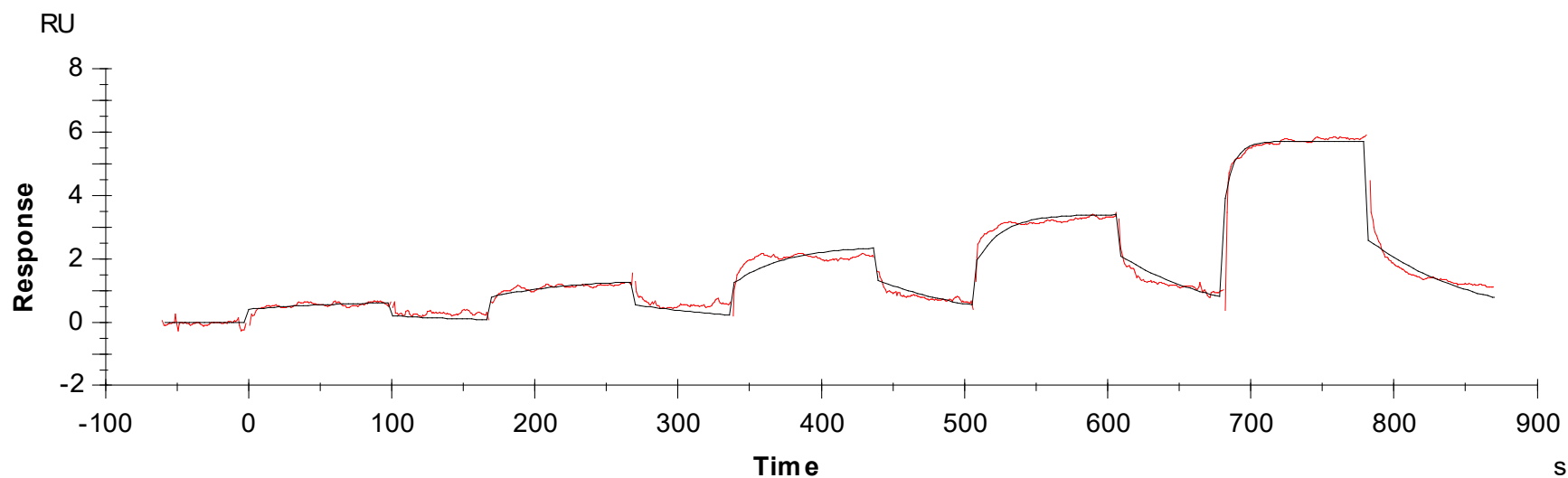




Binding Kinetics of Ibrutinib on BTK[V416L][non-activated]]



Data generated and provided by Carna Biosciences, Inc.