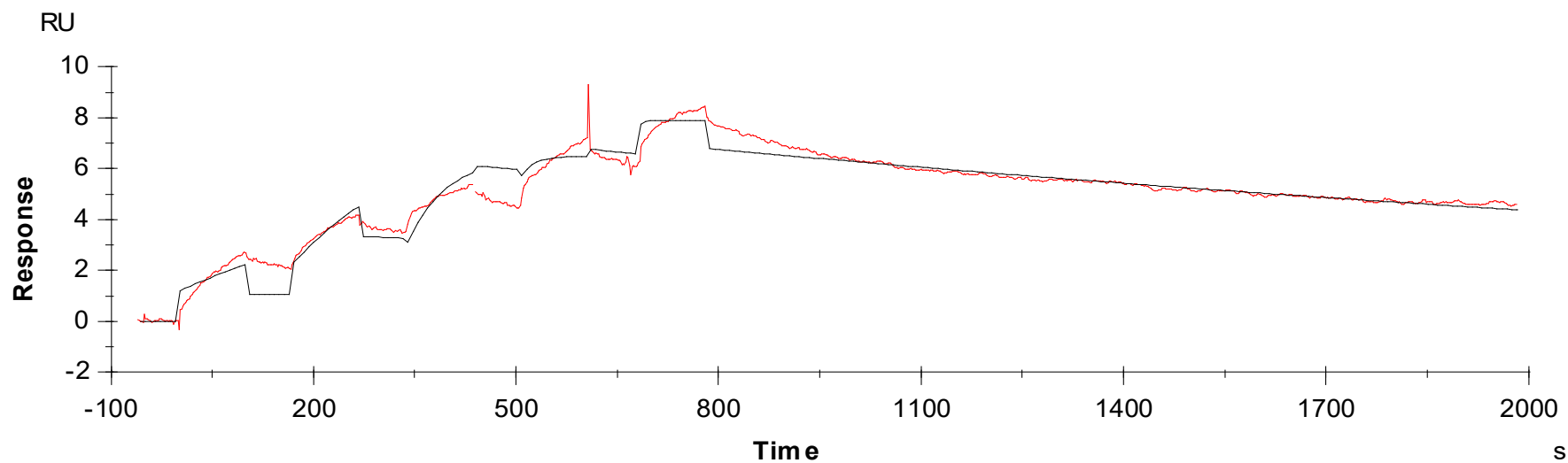




Binding Kinetics of Ibrutinib on BTK[M437R][non-activated]



Data generated and provided by Carna Biosciences, Inc.