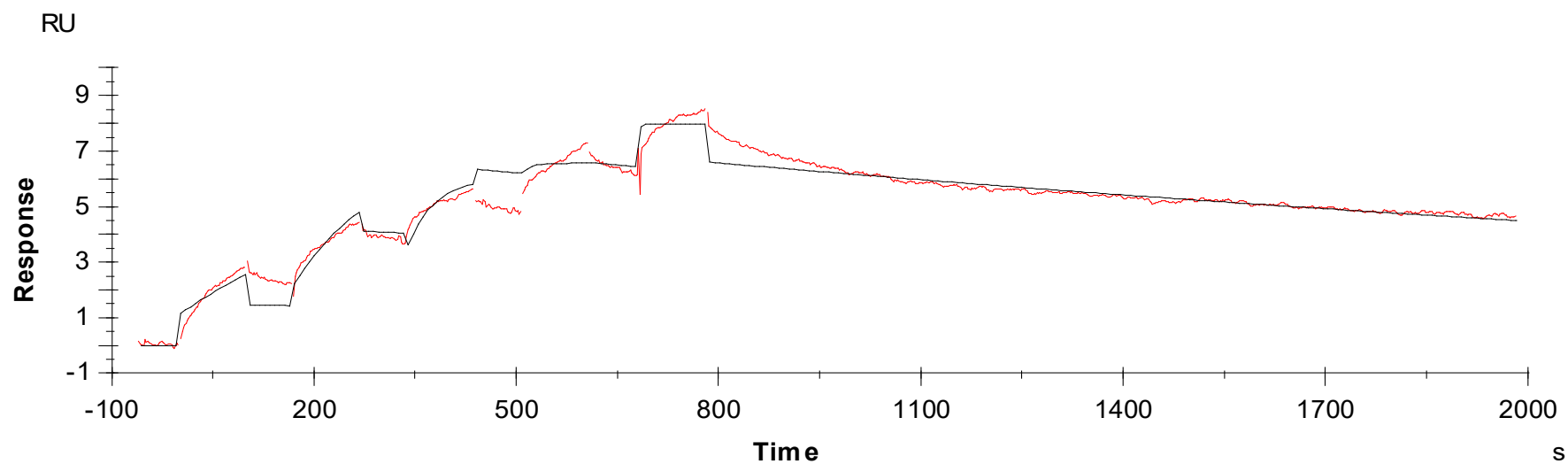




Binding Kinetics of Ibrutinib on BTK[M437R]



Data generated and provided by Carna Biosciences, Inc.