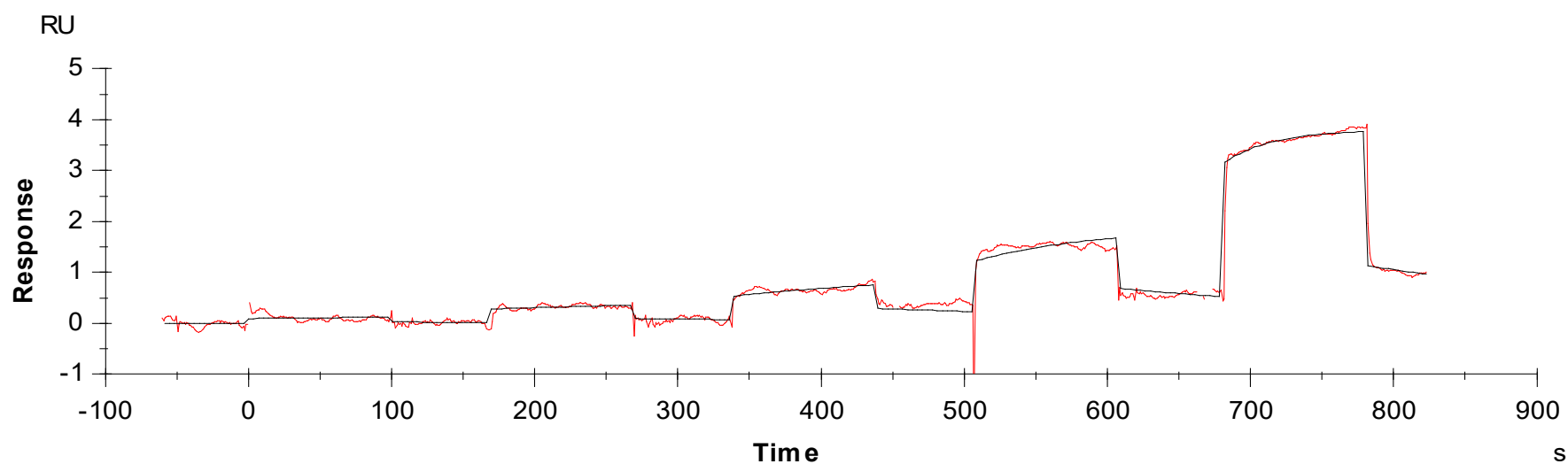




Binding Kinetics of fenebrutinib on BTK[A428D]



Data generated and provided by Carna Biosciences, Inc.