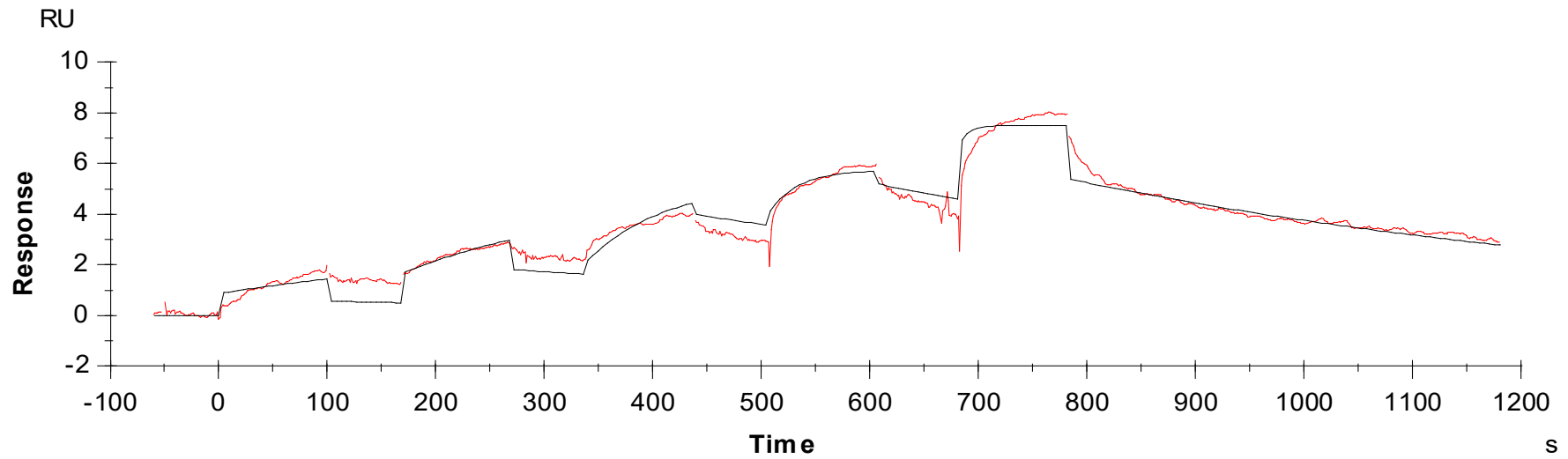




Binding Kinetics of Dasatinib on TEC(non-activated)



Data generated and provided by Carna Biosciences, Inc.