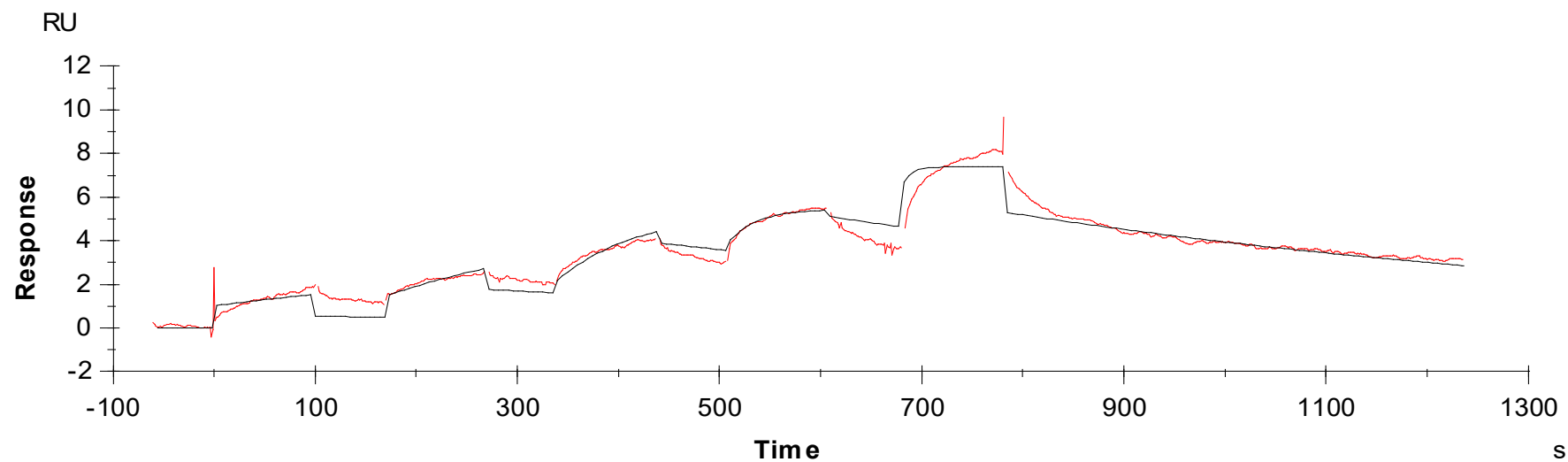




Binding Kinetics of Dasatinib on TEC



Data generated and provided by Carna Biosciences, Inc.