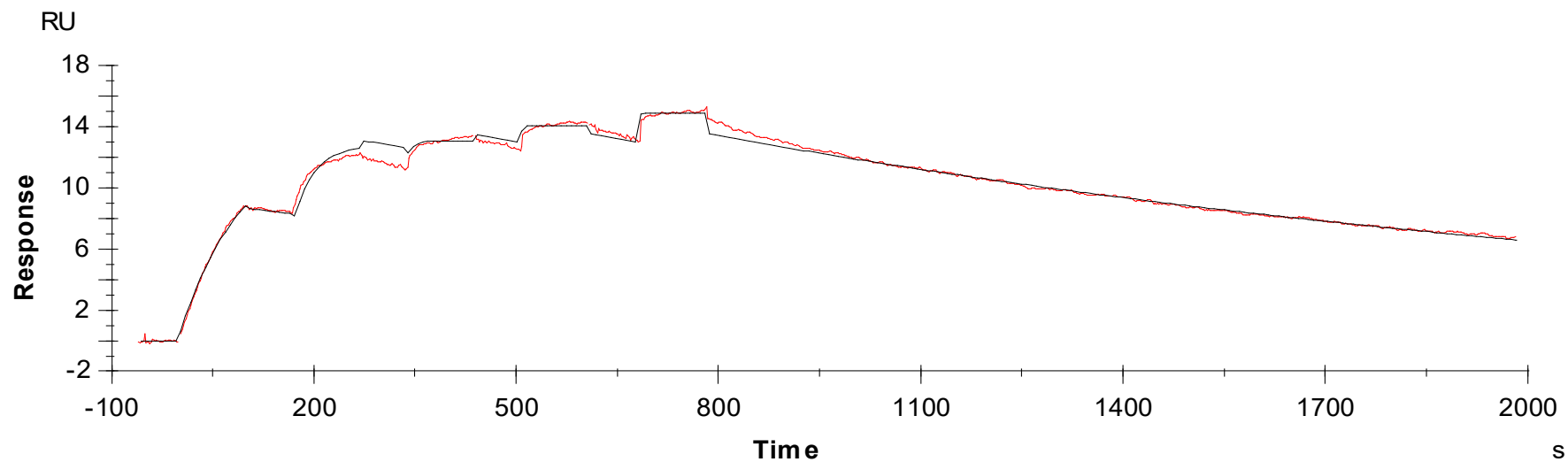




Binding Kinetics of Dasatinib on TXK



Data generated and provided by Carna Biosciences, Inc.