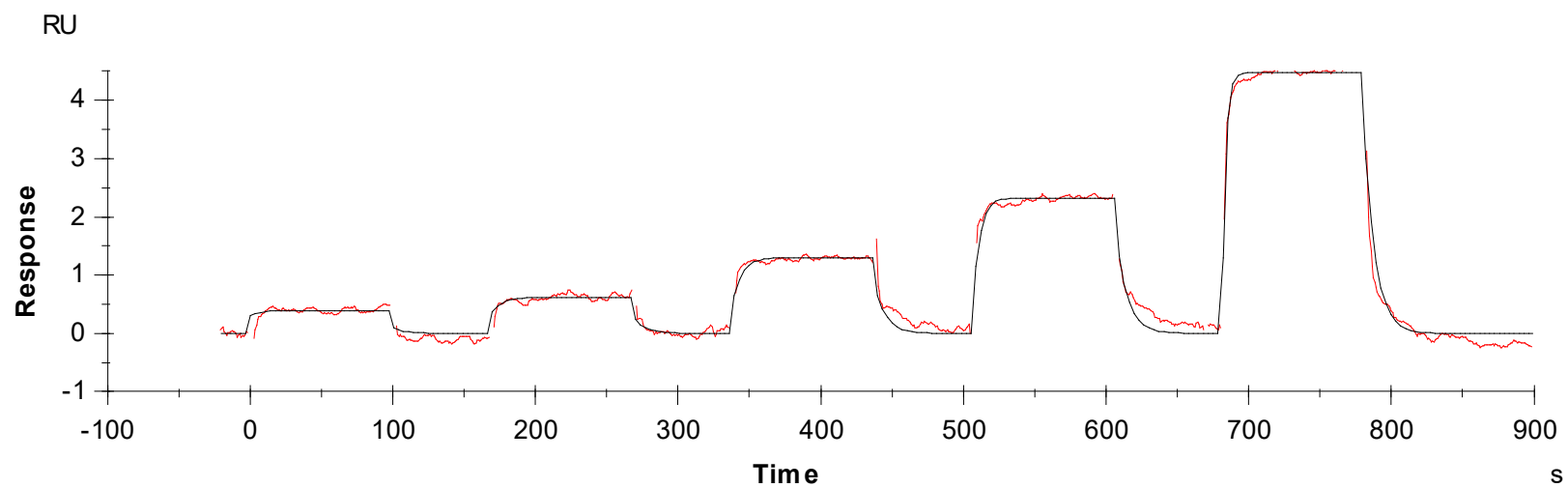


Binding Kinetics of dasatinib on STRADA



Data generated and provided by Carna Biosciences, Inc.