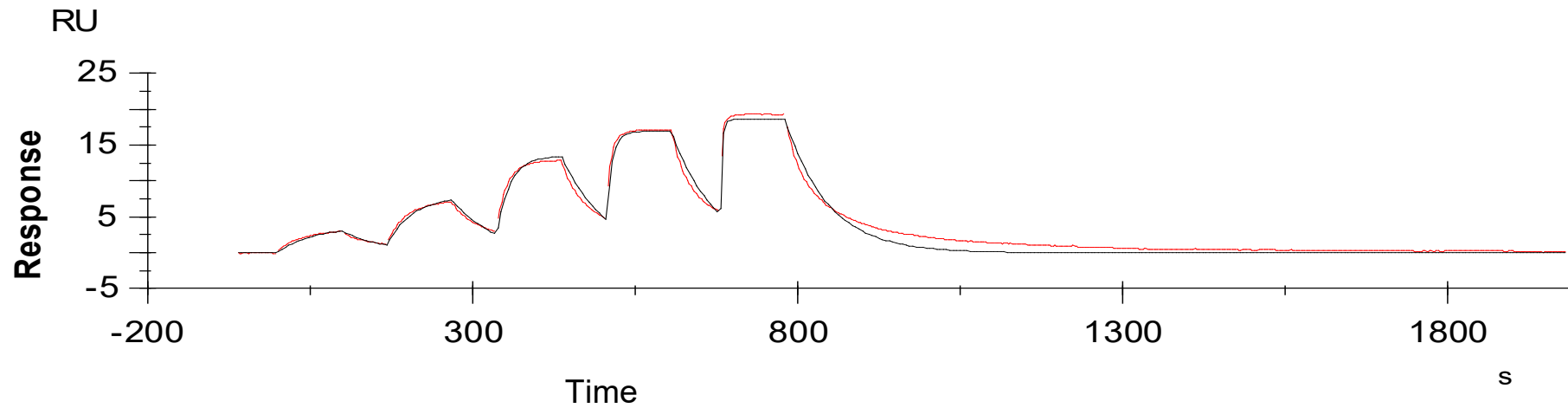


Binding Kinetics of Dasatinib on ACVR2A



Data generated and provided by Carna Biosciences, Inc.