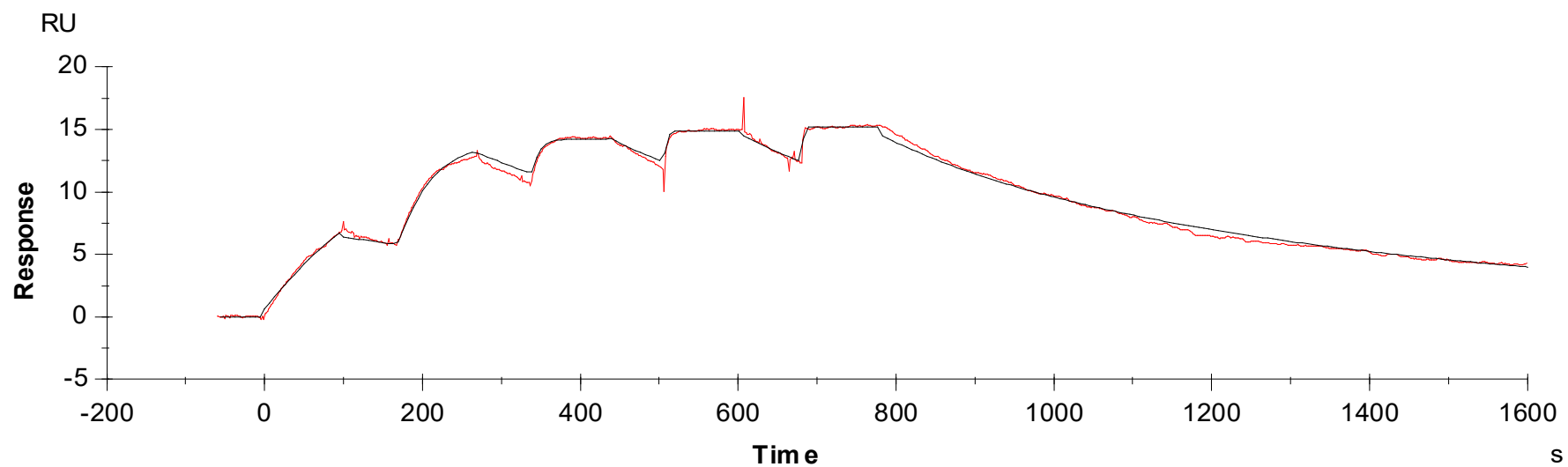




## Binding Kinetics of Dasatinib on BRK



Data generated and provided by Carna Biosciences, Inc.