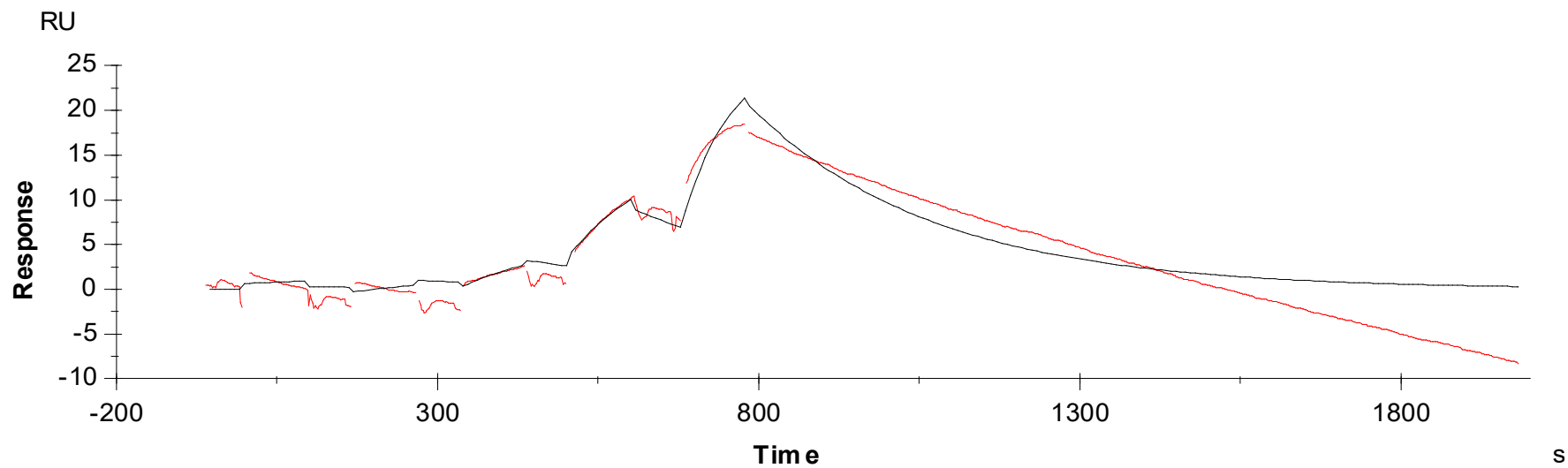


## Binding Kinetics of imatinib on DDR1



Data generated and provided by Carna Biosciences, Inc.