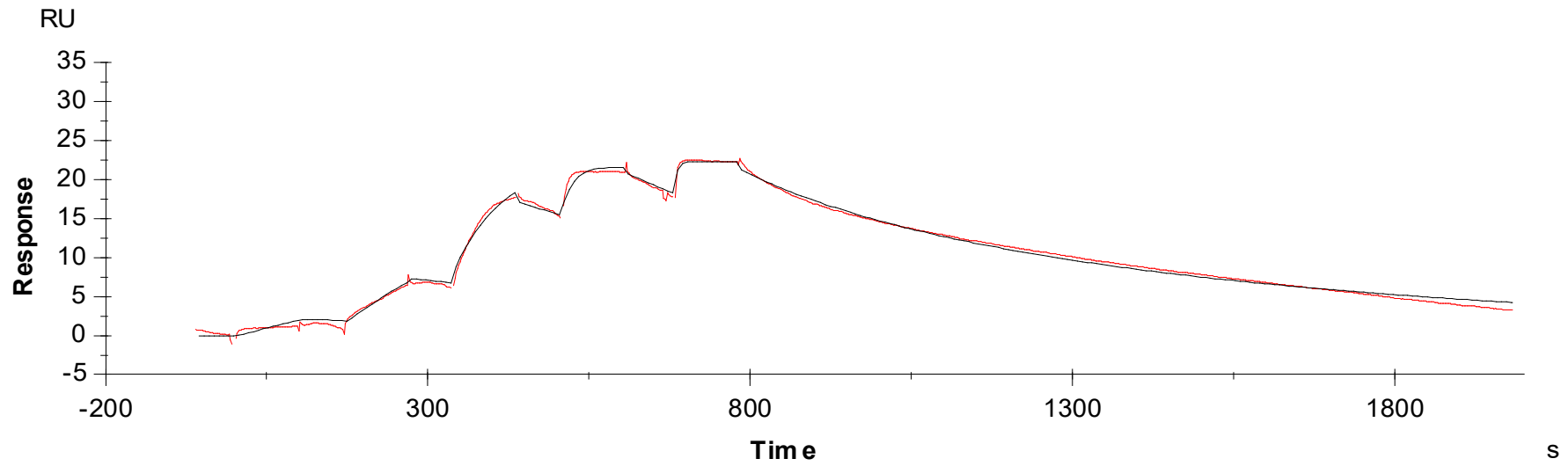


Binding Kinetics of Afatinib on EGFR[C797S]



Data generated and provided by Carna Biosciences, Inc.