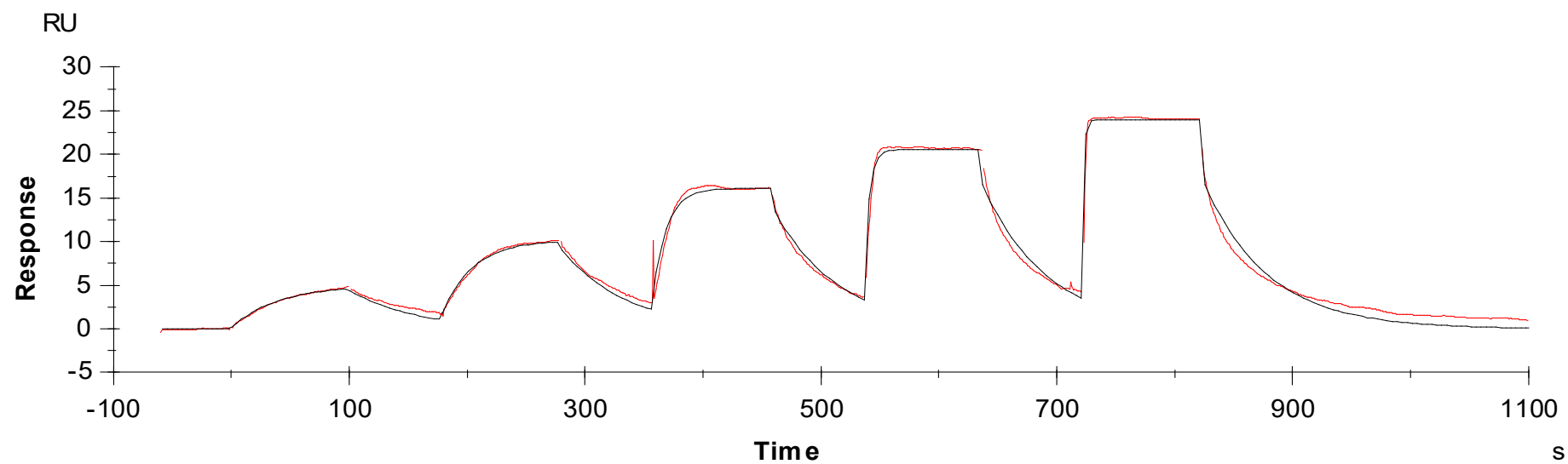


Binding Kinetics of Axitinib on FGFR2[N549H]



Data generated and provided by Carna Biosciences, Inc.