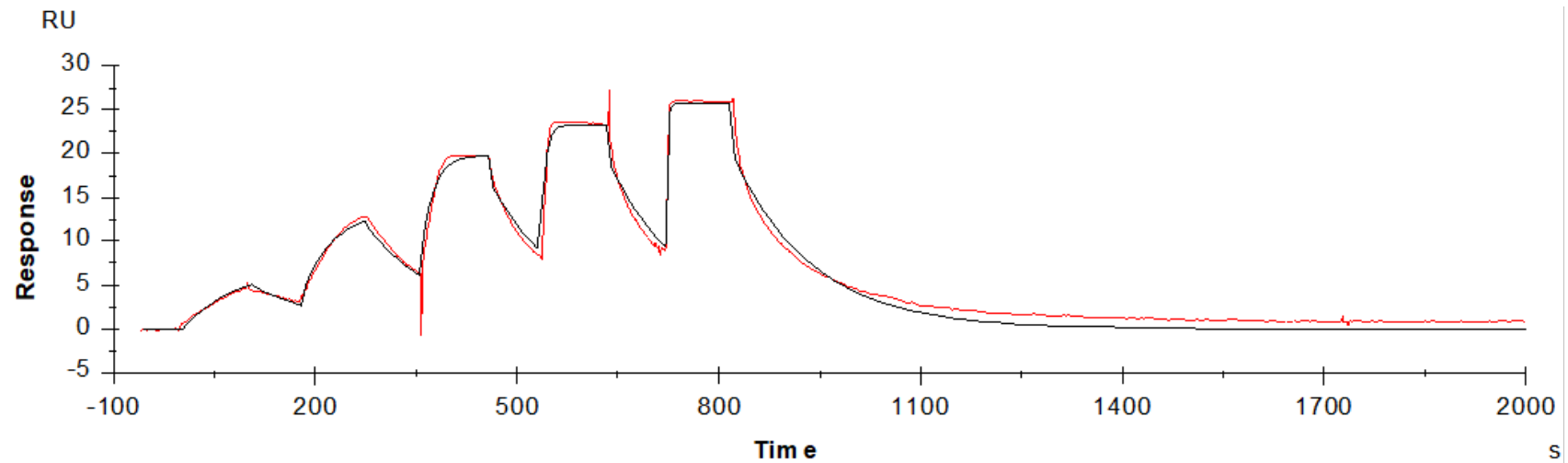


Binding Kinetics of axitinib on FGFR2[N549H](non-activated)



Data generated and provided by Carna Biosciences, Inc.