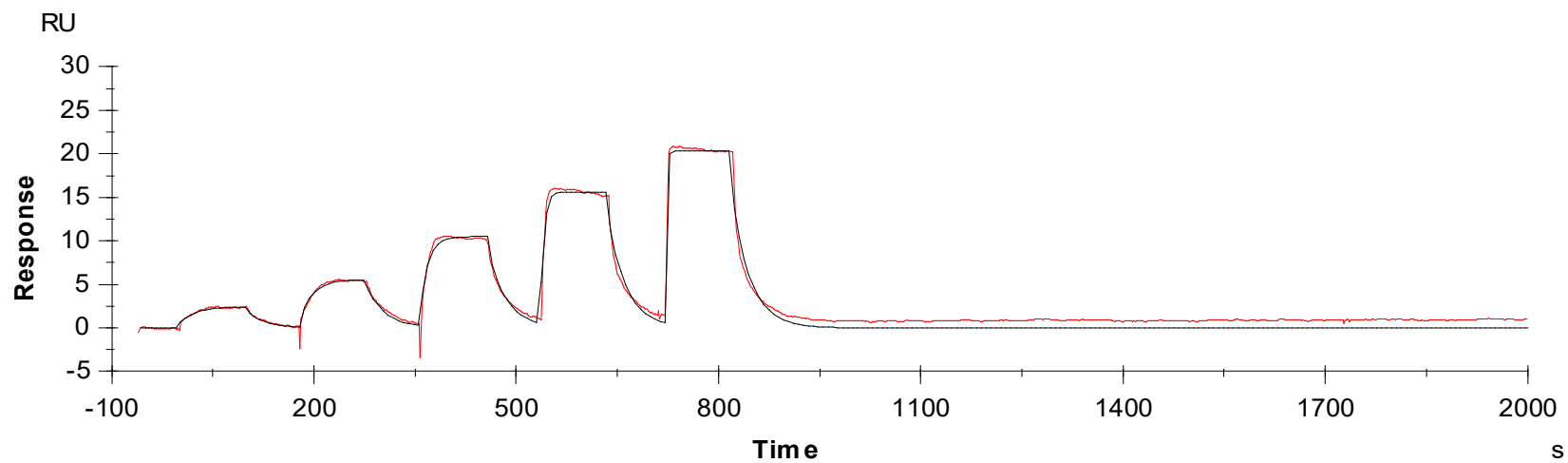


Binding Kinetics of axitinib on FGFR2[V564L]



Data generated and provided by Carna Biosciences, Inc.