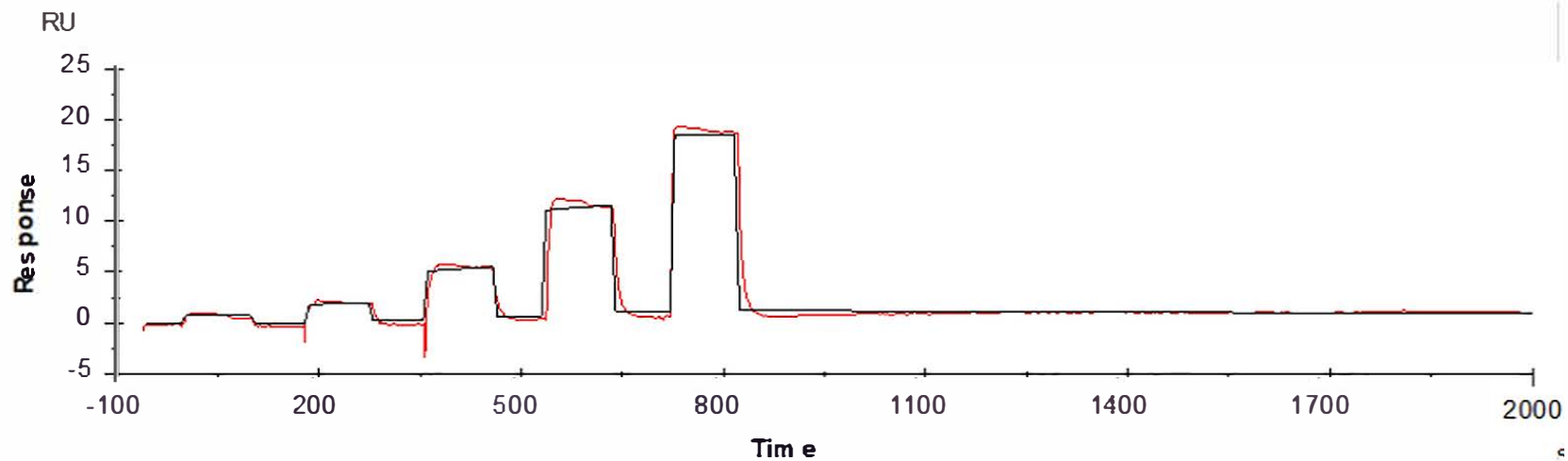


## Binding Kinetics of axitinib on FGFR2[V564L](non-activated)



Data generated and provided by Carna Biosciences, Inc.