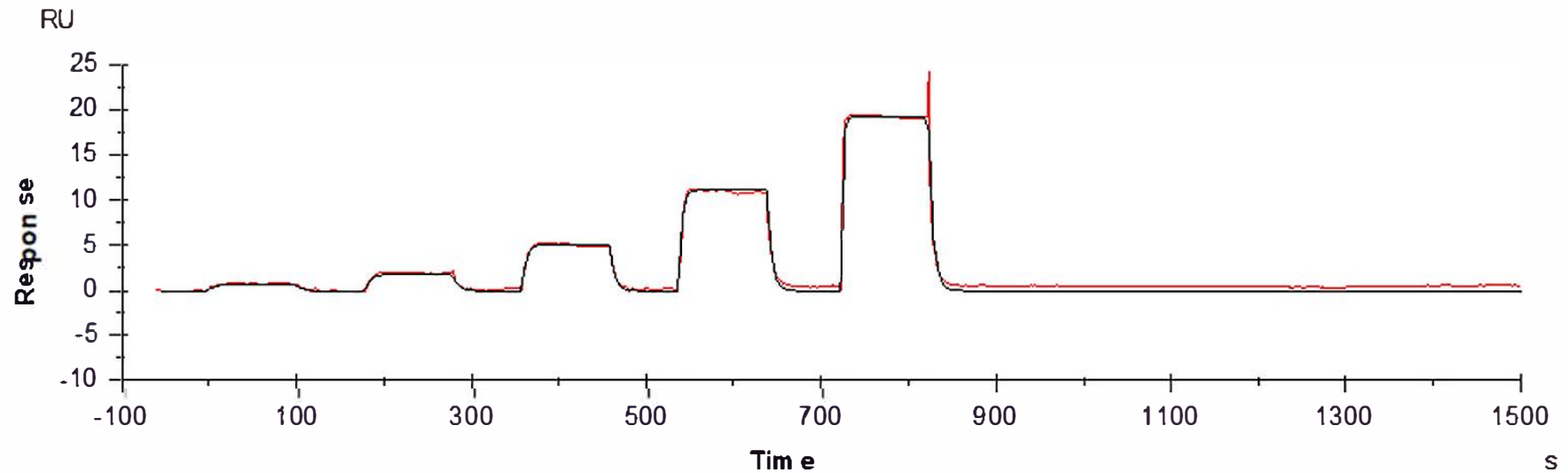


## Binding Kinetics of Axitinib on FGFR2[V564I](non-activated)



Data generated and provided by Carna Biosciences, Inc.