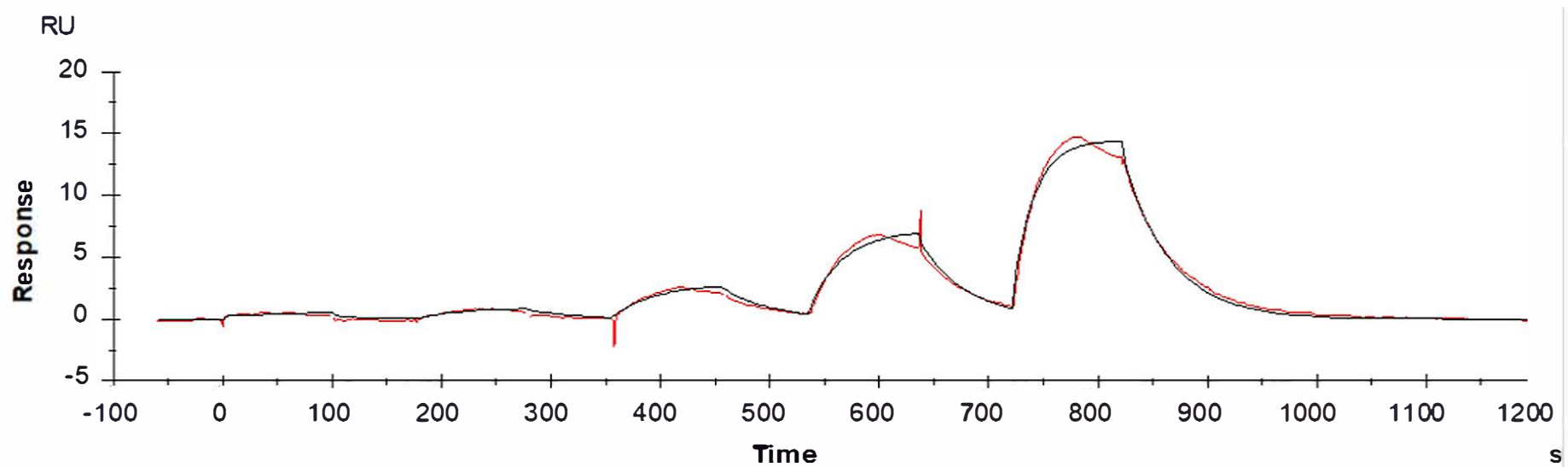


Binding Kinetics of ponatinib on FGFR2[C491A/V564I](non-activated)



Data generated and provided by Carna Biosciences, Inc.