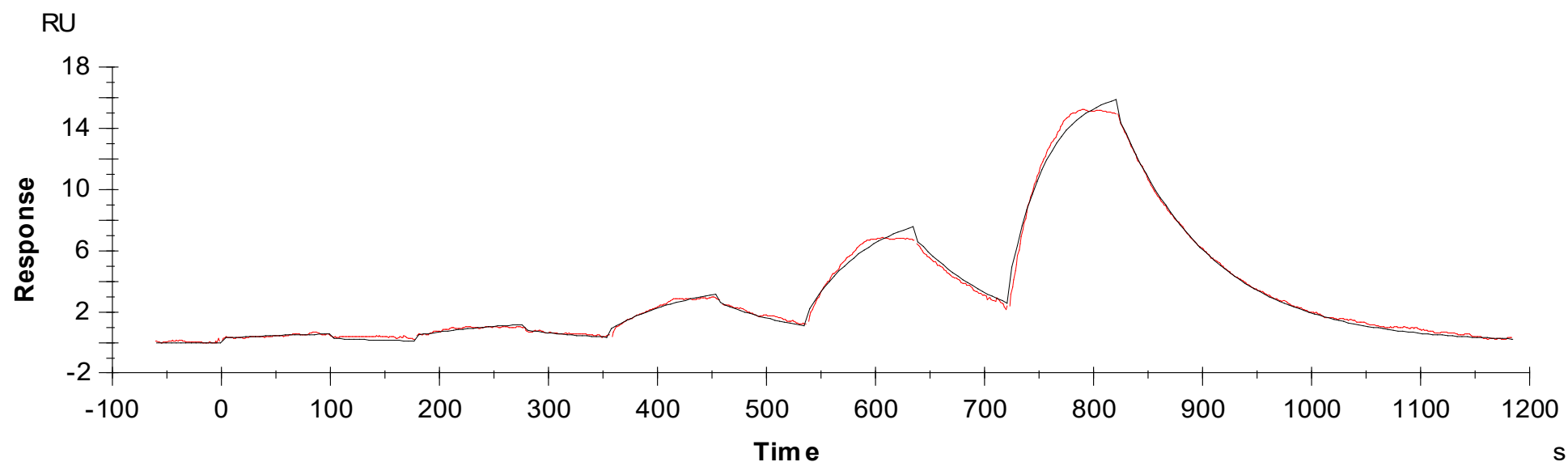


Binding Kinetics of Ponatinib on FGFR2[C491A/V564L][non-activated]



Data generated and provided by Carna Biosciences, Inc.