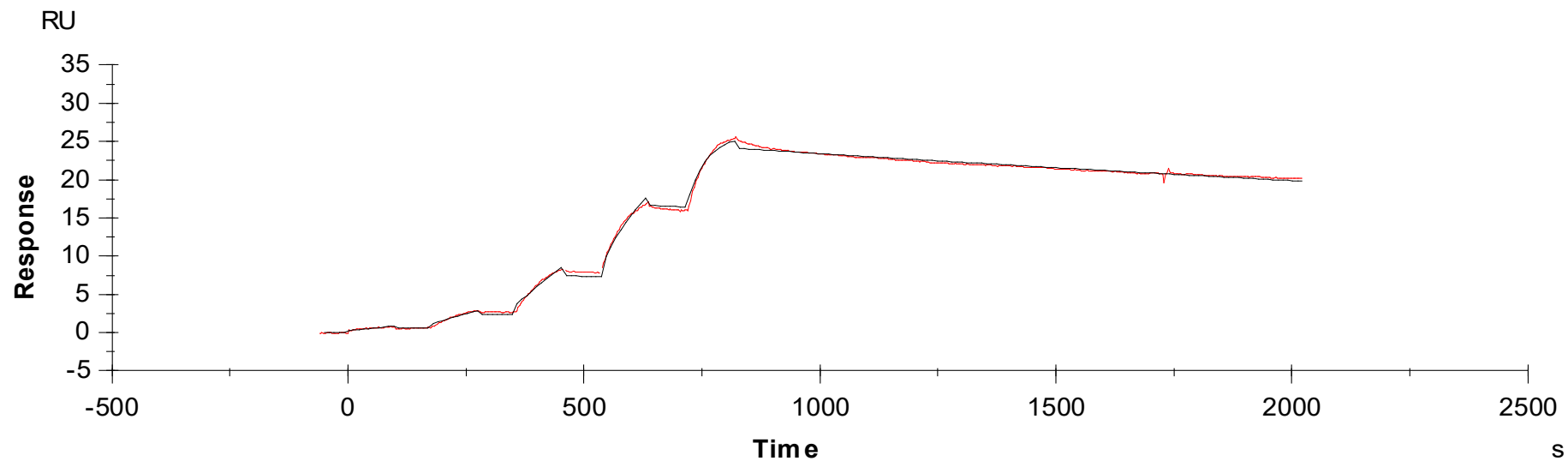


Binding Kinetics of ponatinib on FGFR2[C491A]



Data generated and provided by Carna Biosciences, Inc.