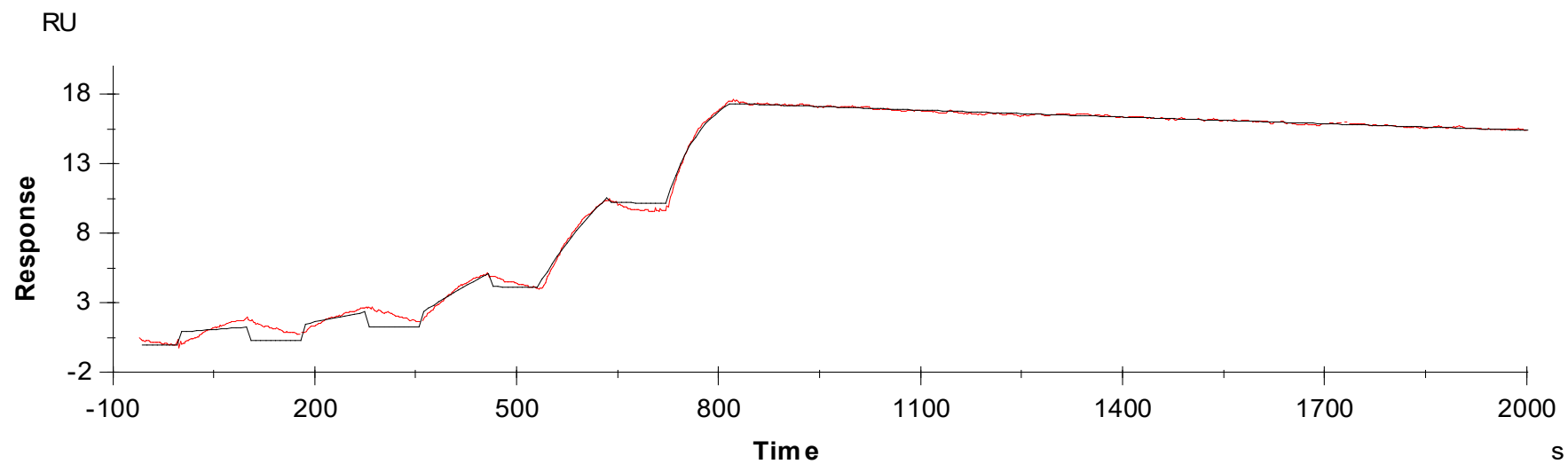


Binding Kinetics of Ponatinib on FGFR2[M537I]



Data generated and provided by Carna Biosciences, Inc.