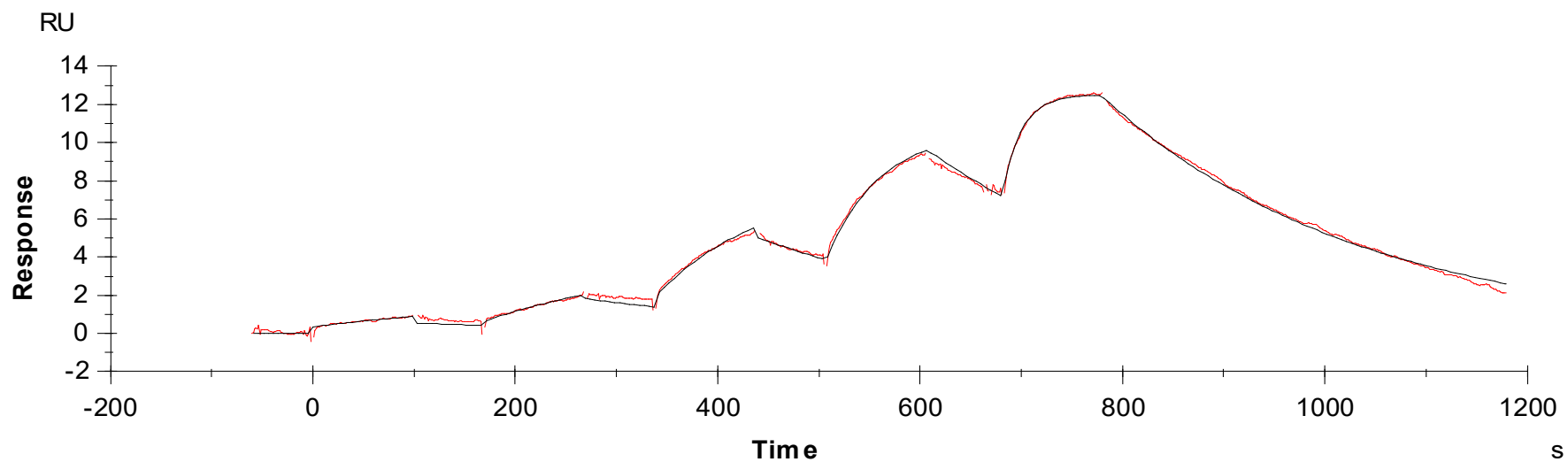


Binding Kinetics of Met/Ron inhibitor on MET[Y1230H]



Data generated and provided by Carna Biosciences, Inc.