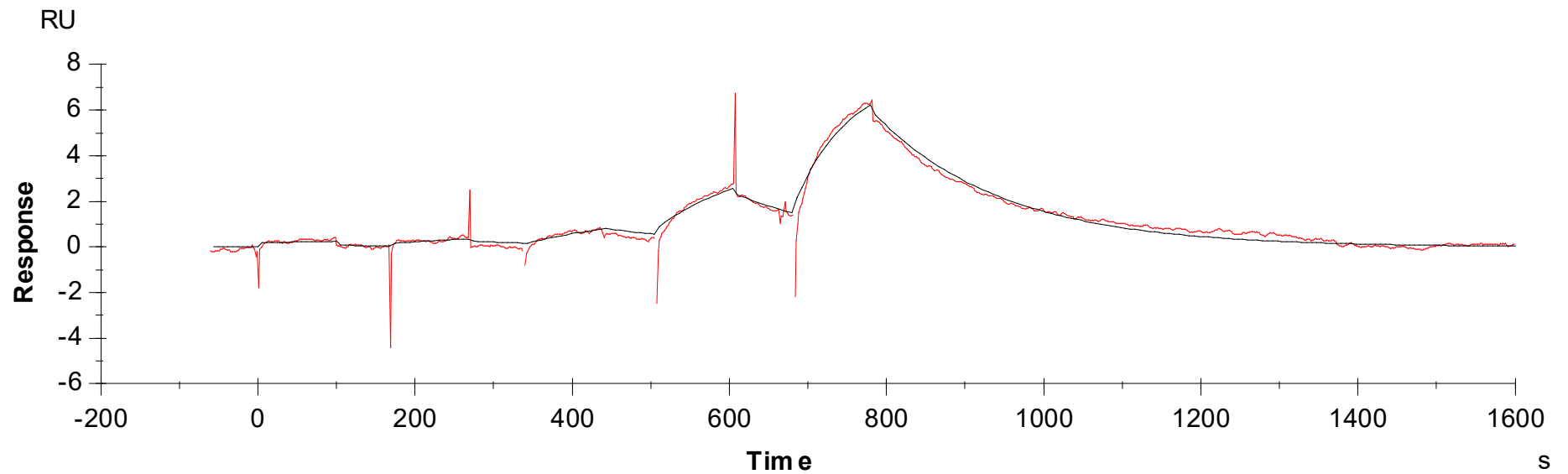


Binding Kinetics of Met/Ron inhibitor on MET[M1250T]



Data generated and provided by Carna Biosciences, Inc.