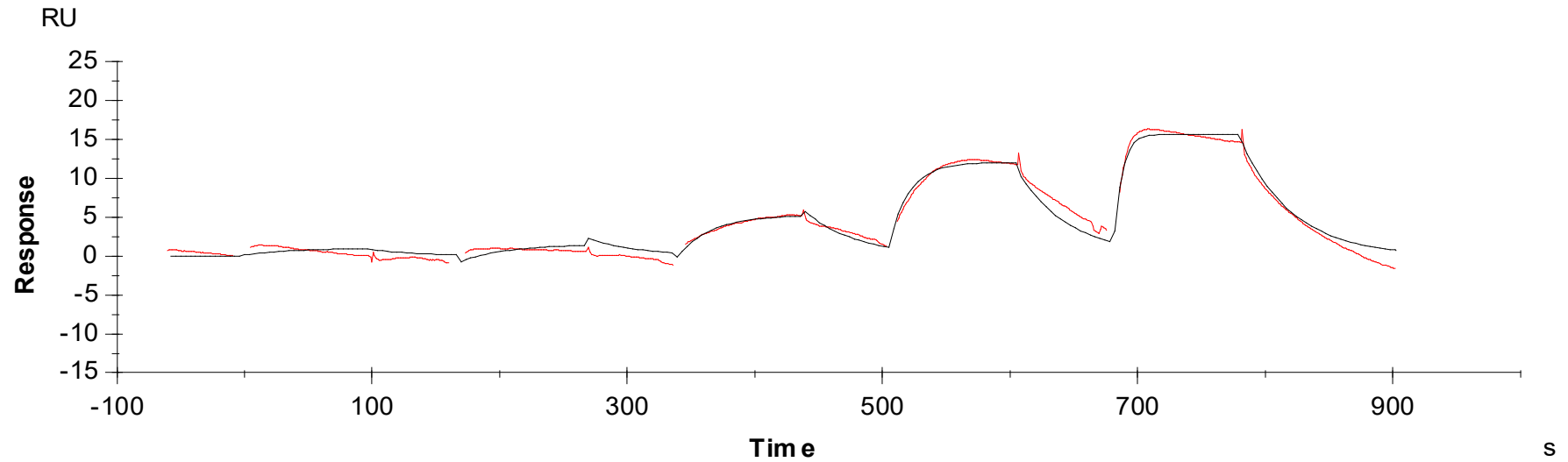


## Binding Kinetics of Crizotinib on MET[Y1230C]



Data generated and provided by Carna Biosciences, Inc.