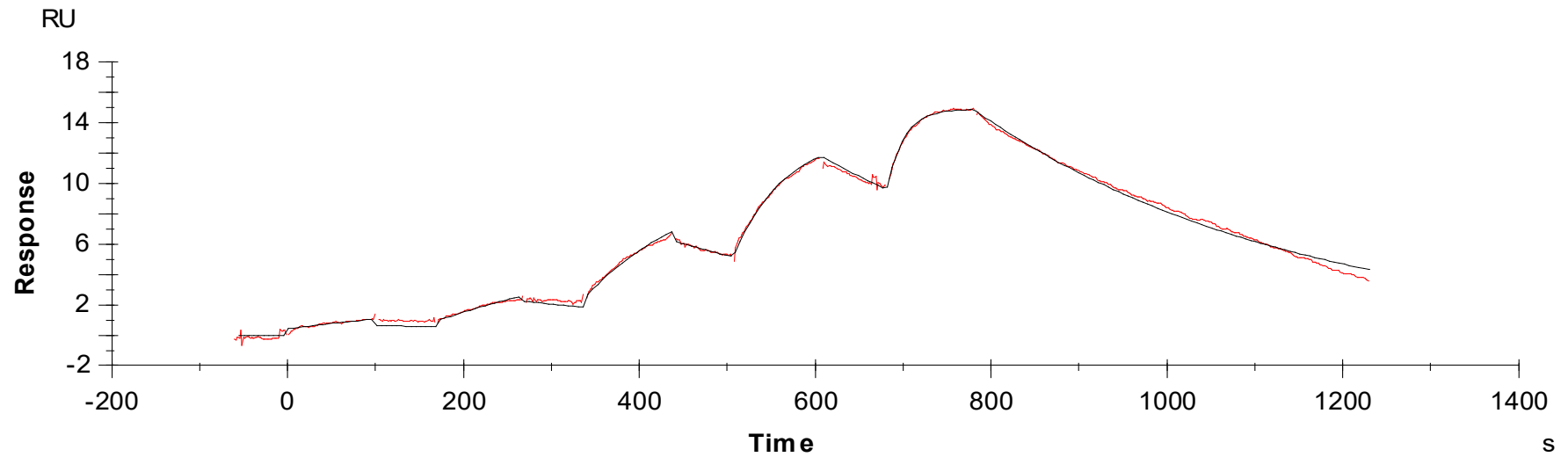


## Binding Kinetics of Met/Ron inhibitor on MET



Data generated and provided by Carna Biosciences, Inc.

# Binding kinetics of crizotinib on MET

