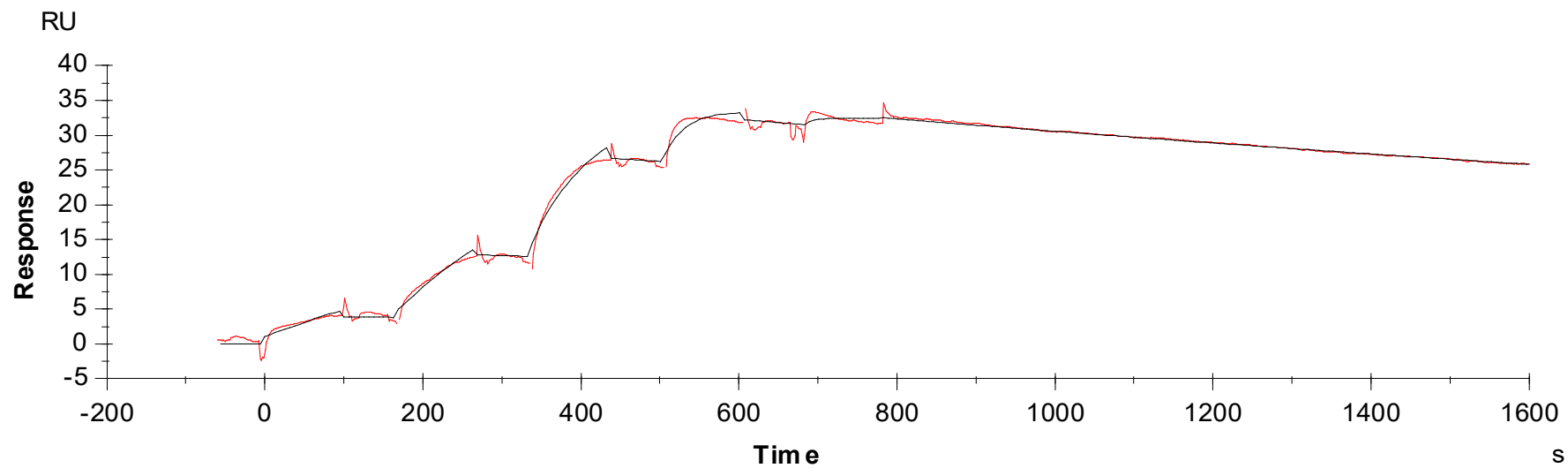


Binding Kinetics of Foretinib on MET[D1228H]



Data generated and provided by Carna Biosciences, Inc.