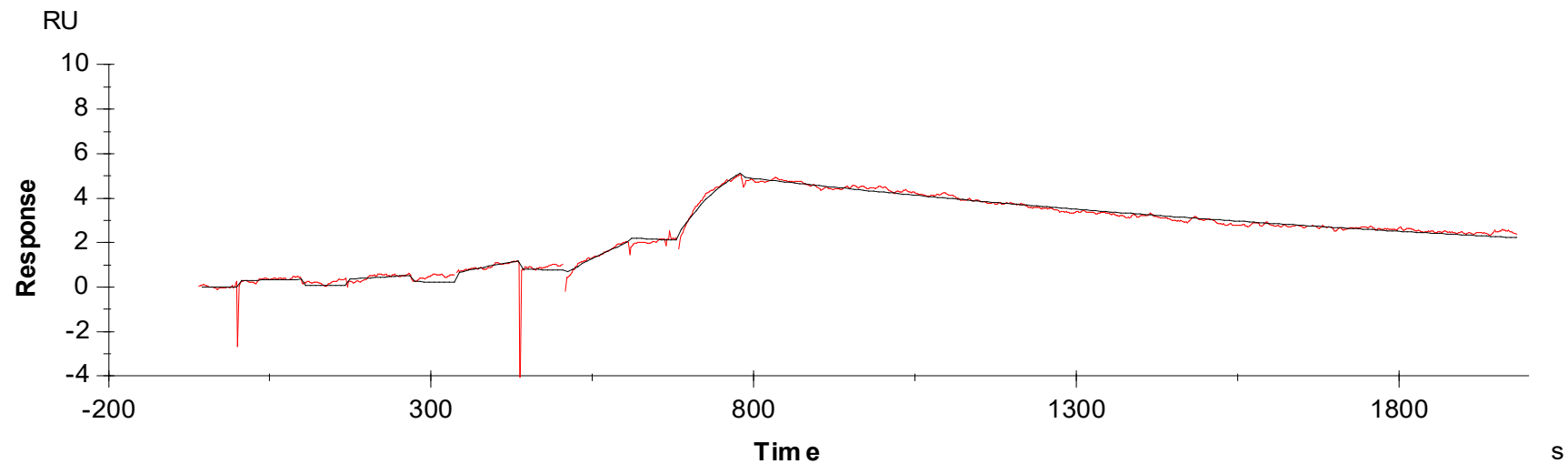




Binding Kinetics of Foretinib on MET[D1228N]



Data generated and provided by Carna Biosciences, Inc.