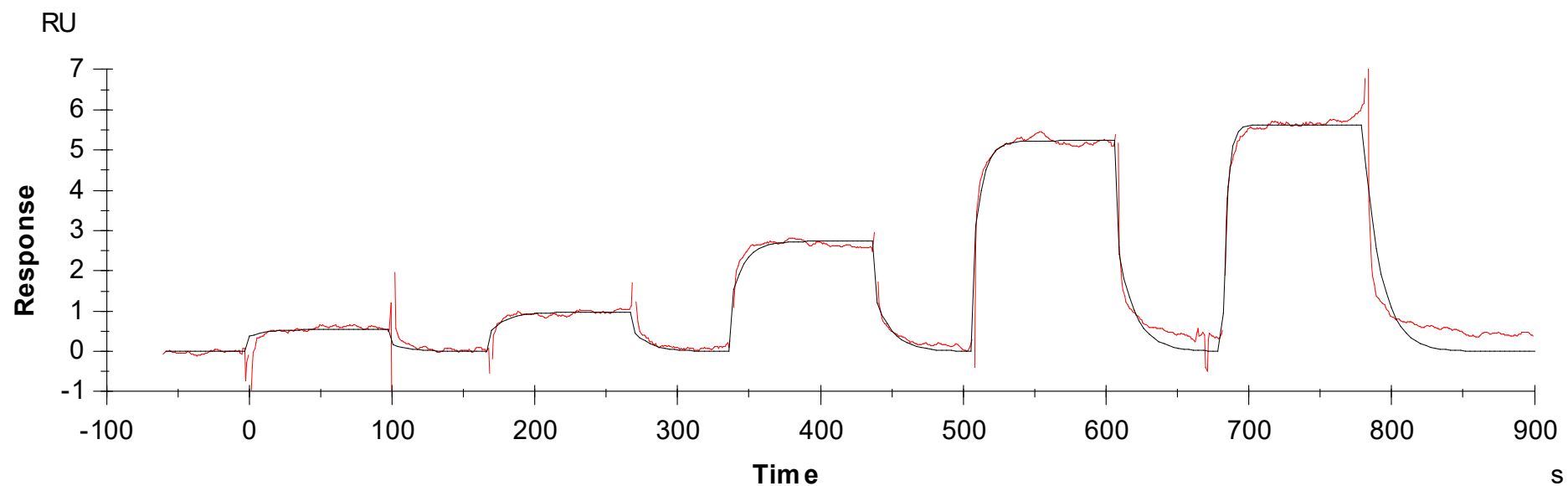


Binding Kinetics of midostaurine on JAK1(JH2)



Data generated and provided by Carna Biosciences, Inc.