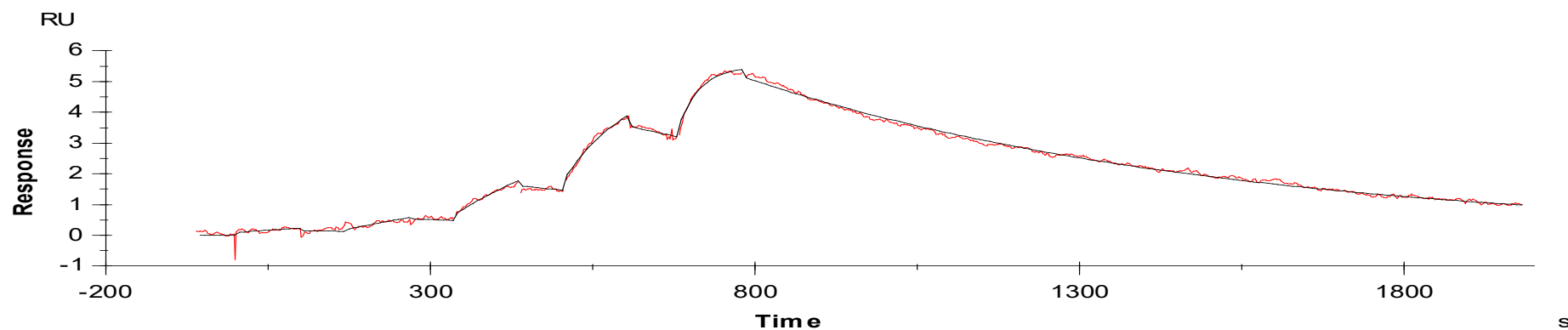


Binding Kinetics of Midostaurin on MST2



Data generated and provided by Carna Biosciences, Inc.