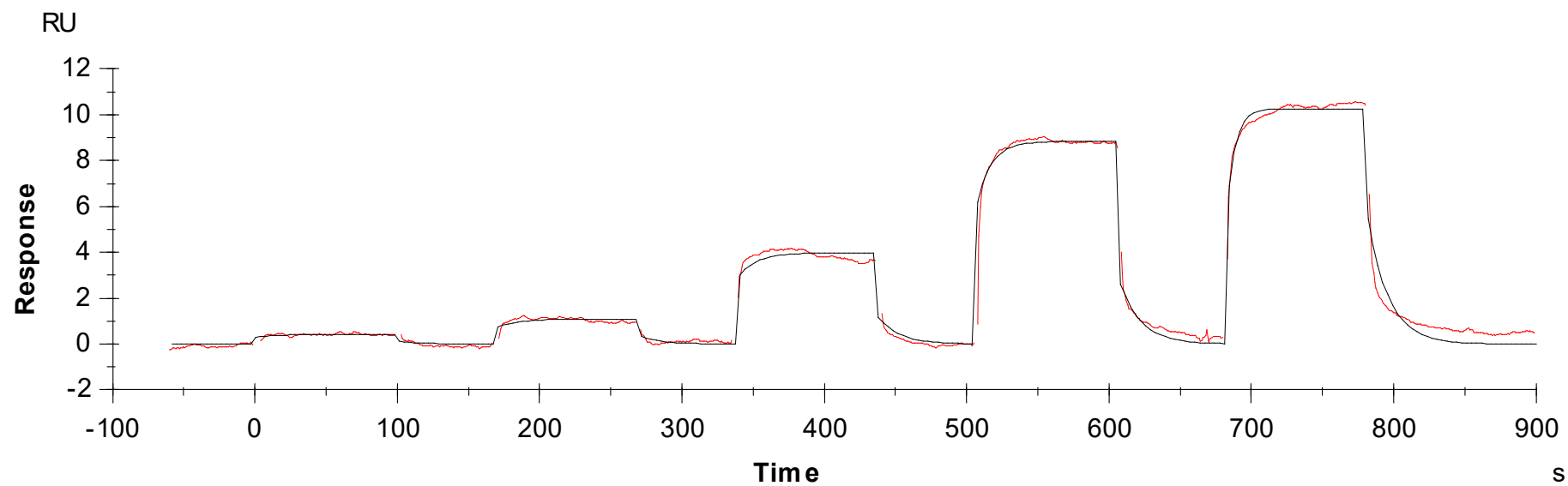


Binding Kinetics of midostaurine on PAN3



Data generated and provided by Carna Biosciences, Inc.