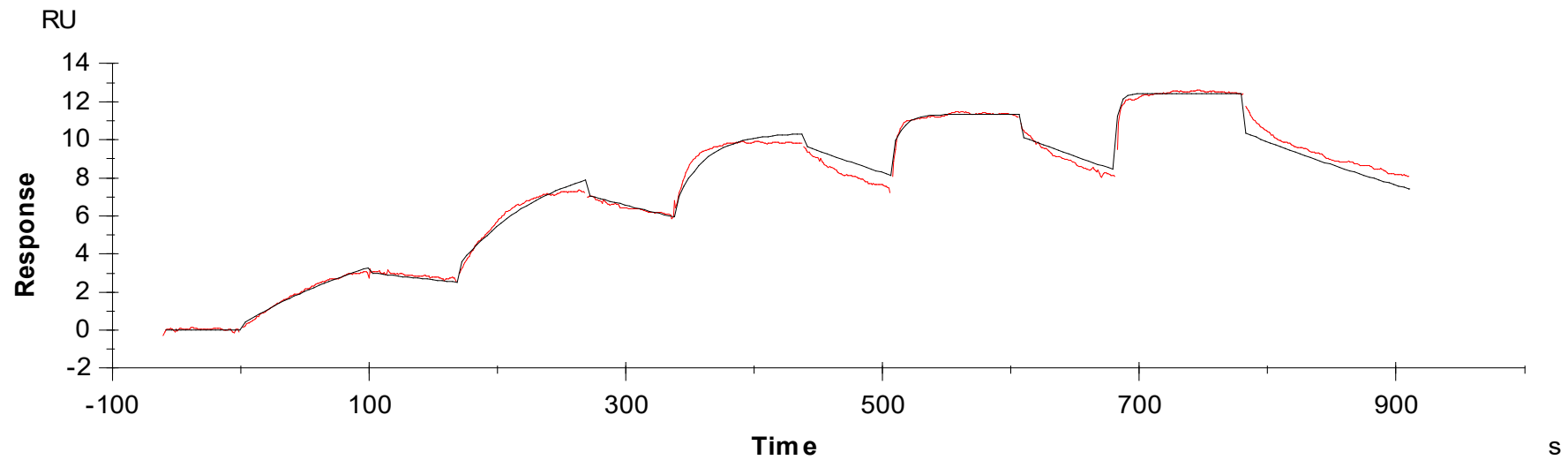




Binding Kinetics of Staurosporine on IRAK3



Data generated and provided by Carna Biosciences, Inc.