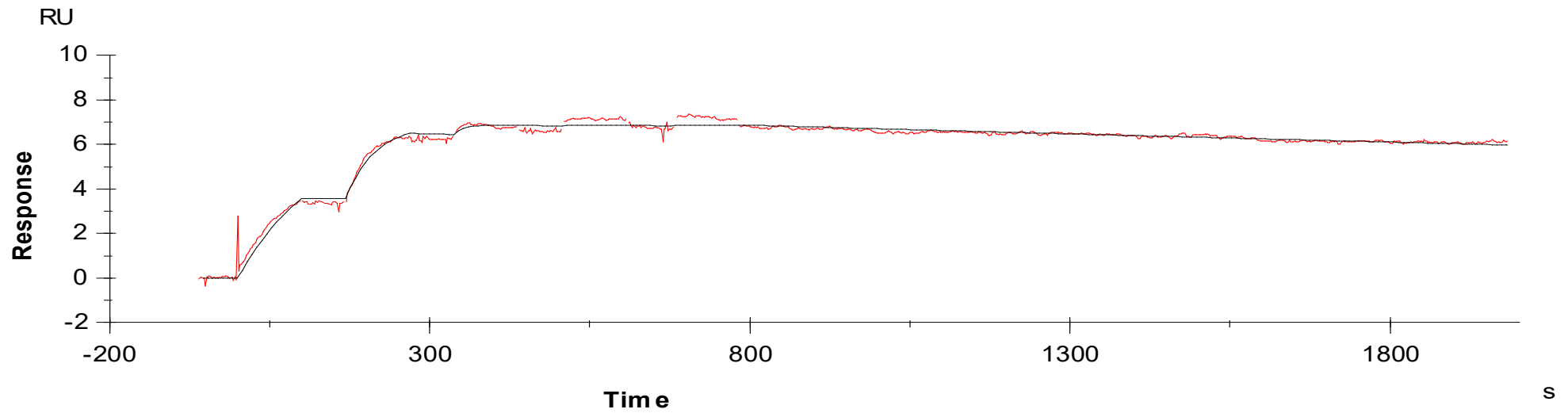


Binding Kinetics of Staurosporine on Nuak2



Data generated and provided by Carna Biosciences, Inc.