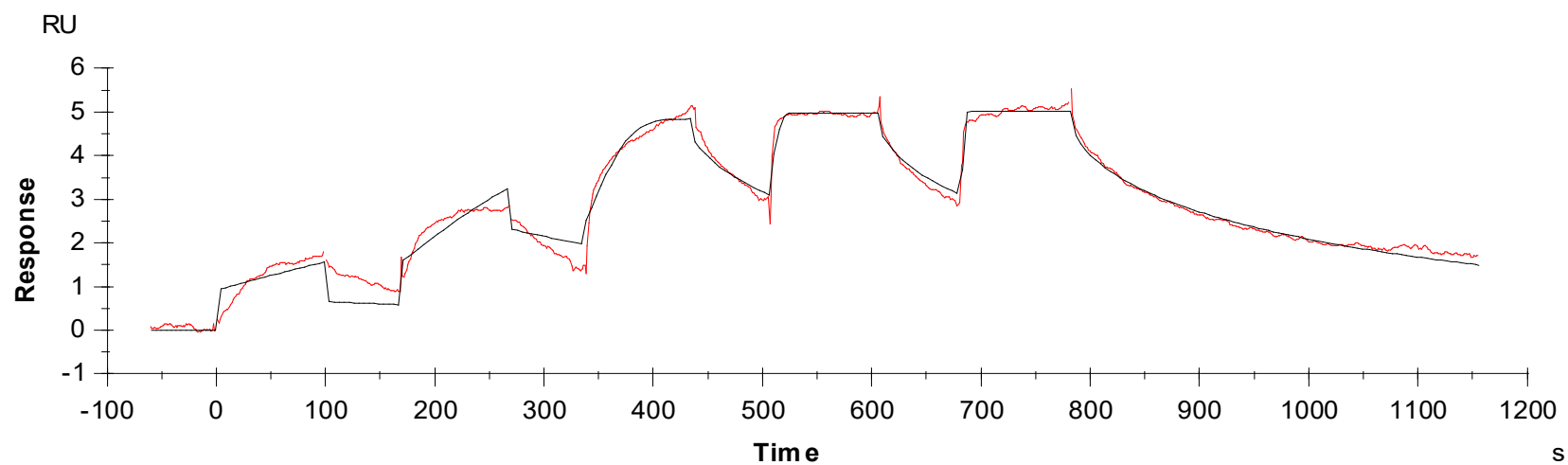




Binding Kinetics of Staurosporine on TSSK2



Data generated and provided by Carna Biosciences, Inc.