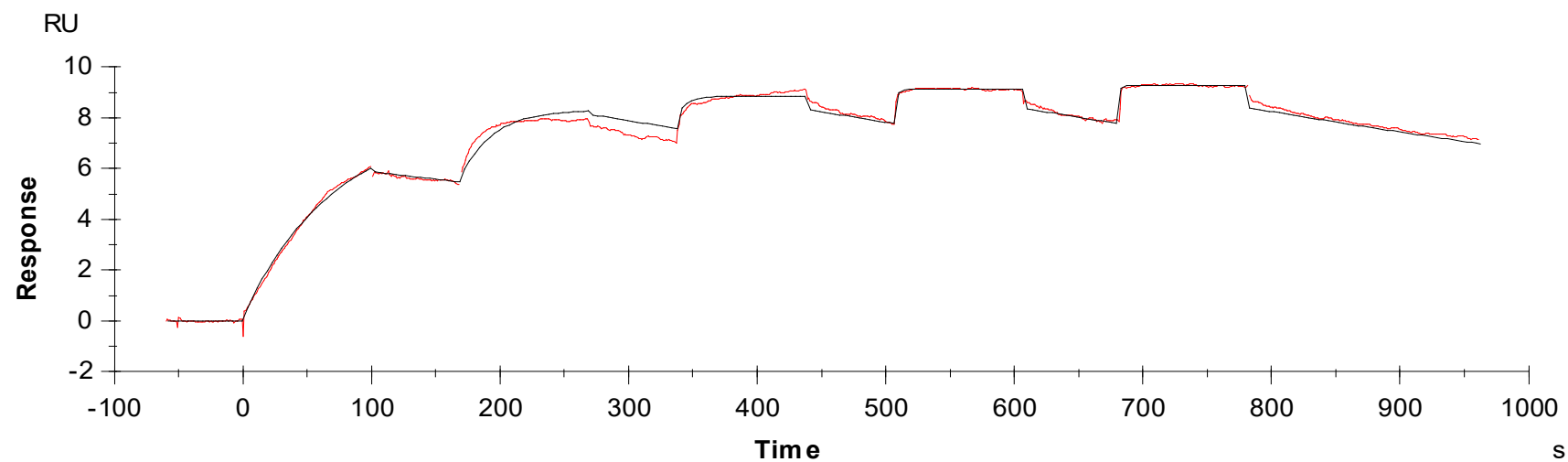




Binding Kinetics of Staurosporine on TSSK1B(TSSK1)



Data generated and provided by Carna Biosciences, Inc.