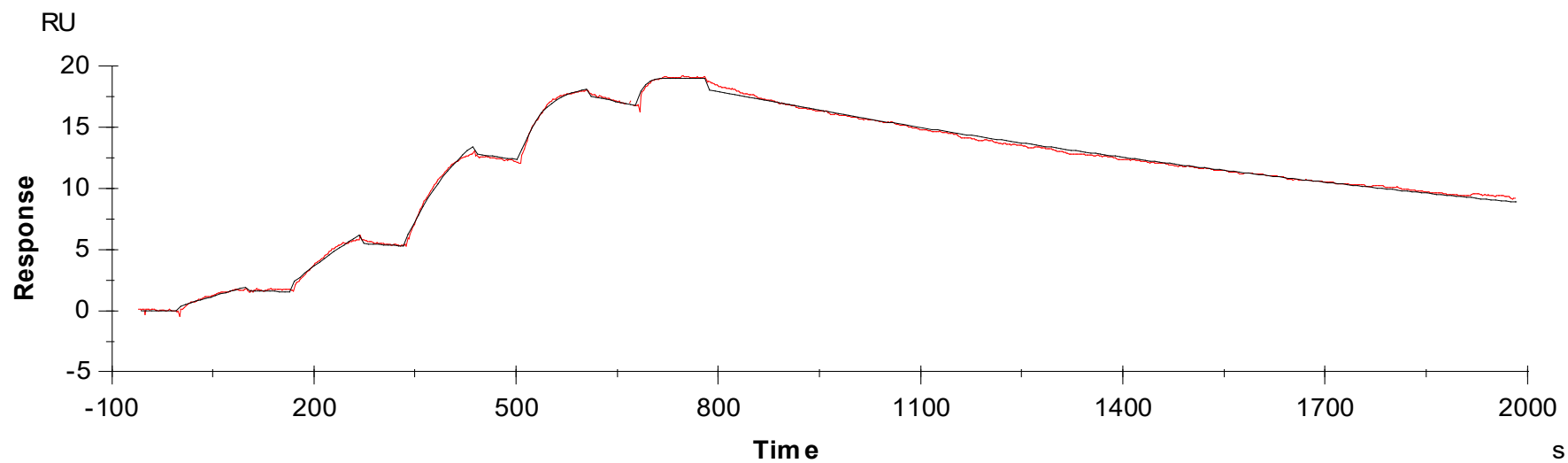




Binding Kinetics of Staurosporine on ROCK2



Data generated and provided by Carna Biosciences, Inc.