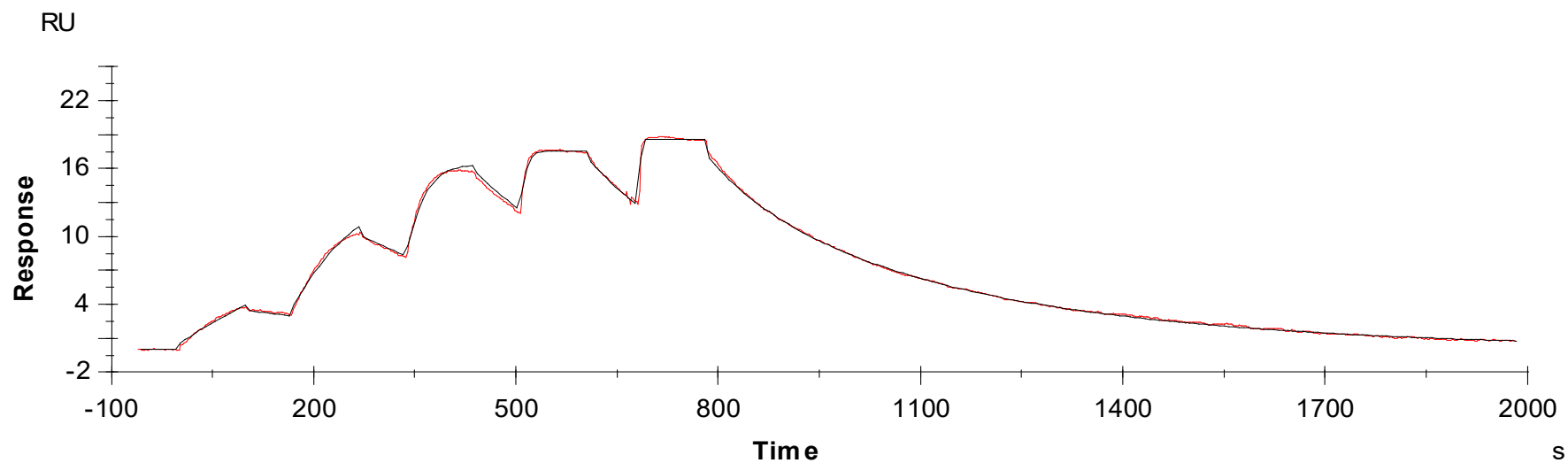




Binding Kinetics of Staurosporine on DYRK1A



Data generated and provided by Carna Biosciences, Inc.