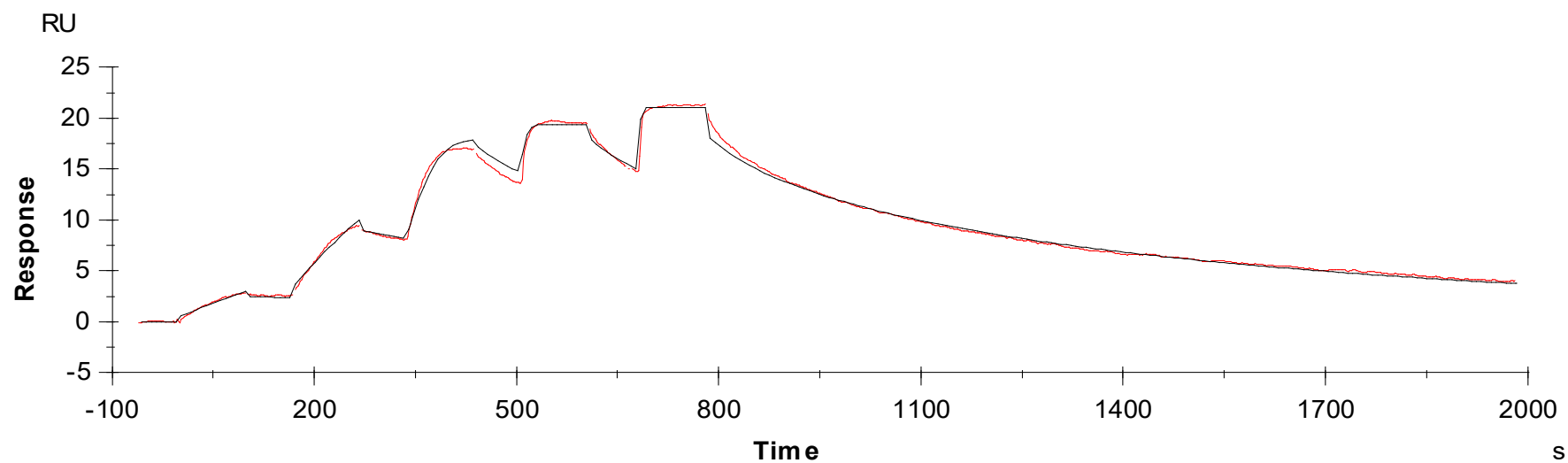




Binding Kinetics of Staurosporine on PRKD2(PKD2)



Data generated and provided by Carna Biosciences, Inc.