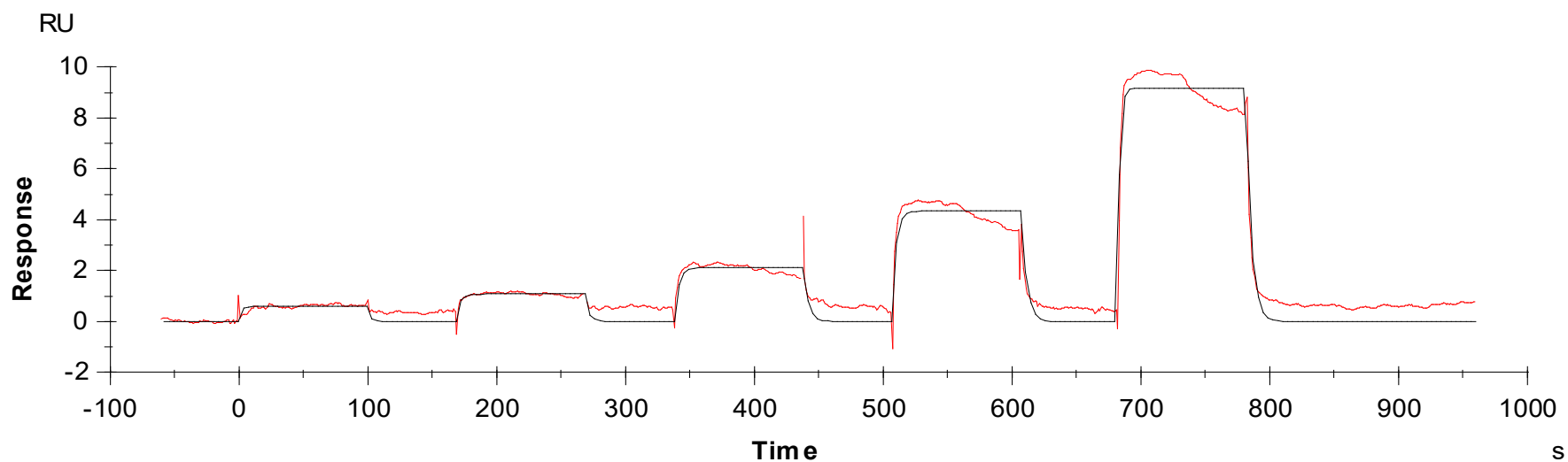




Binding Kinetics of Staurosporine on BARK1(ADRBK1)



Data generated and provided by Carna Biosciences, Inc.