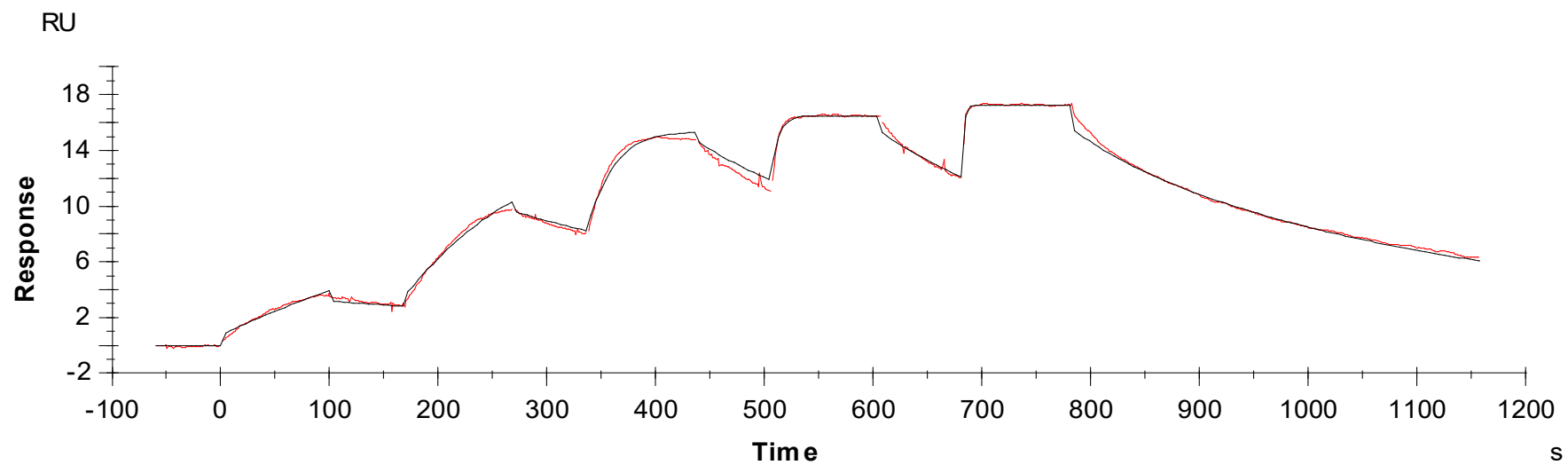




Binding Kinetics of Staurosporine on CDK10(PISSLRE)/CycQ



Data generated and provided by Carna Biosciences, Inc.