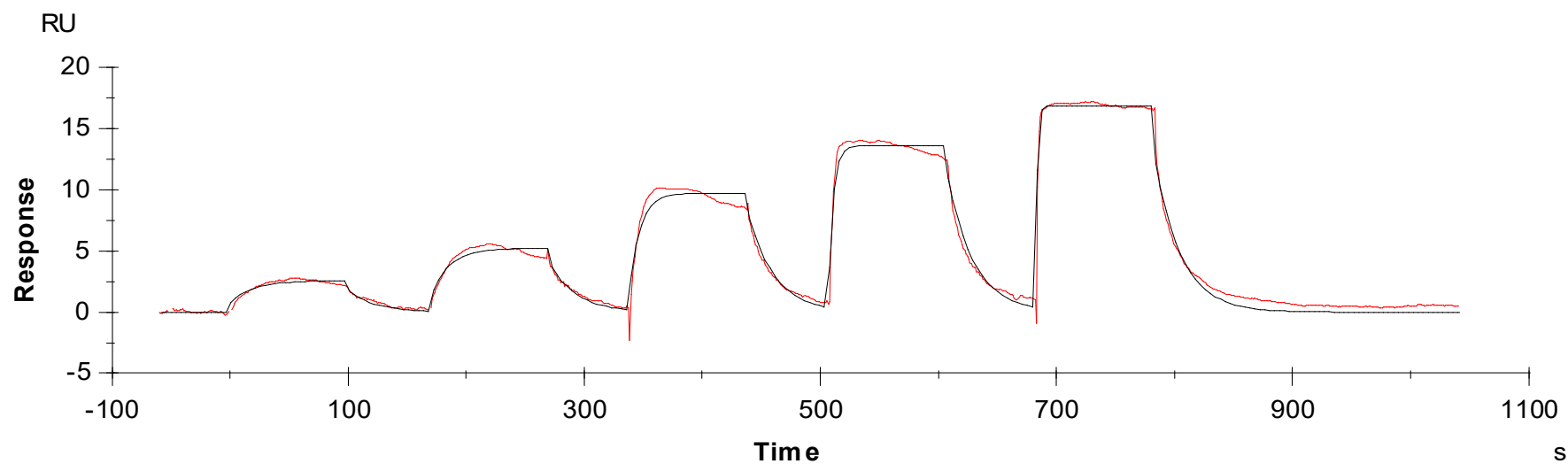




Binding Kinetics of Staurosporine on CDK7



Data generated and provided by Carna Biosciences, Inc.