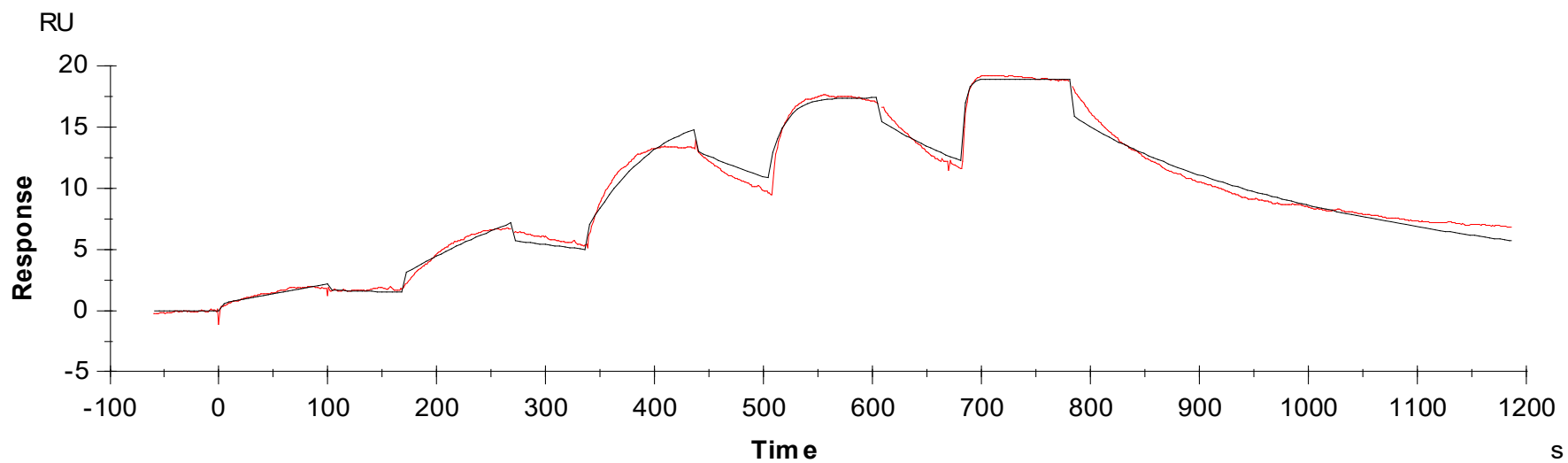


Binding Kinetics of Staurosporine on CLK2



Data generated and provided by Carna Biosciences, Inc.