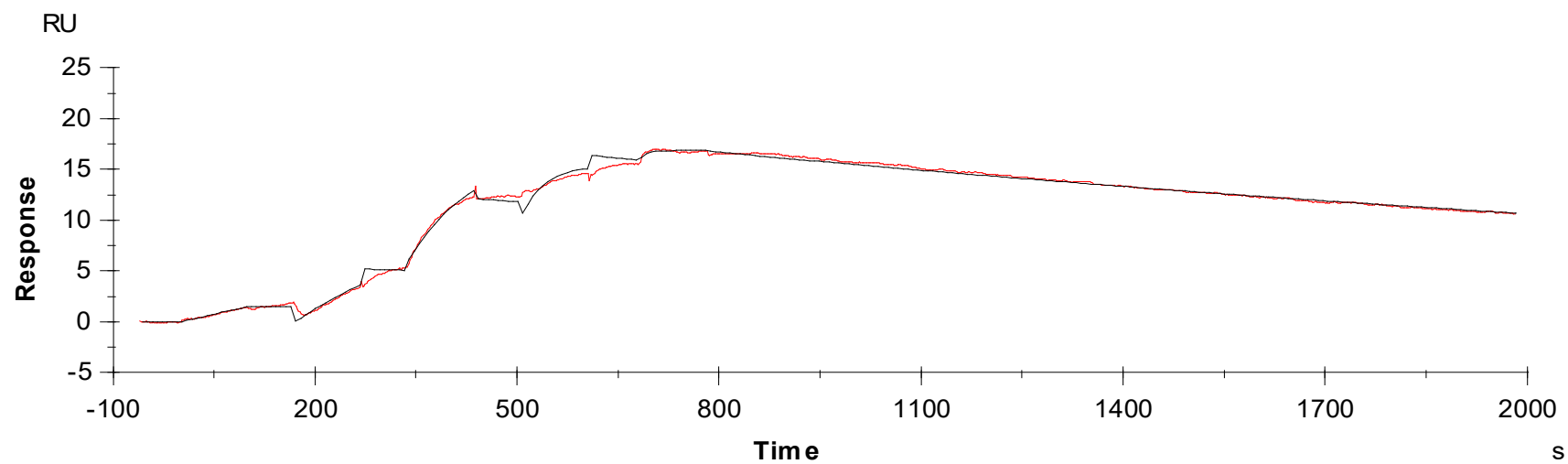




## Binding Kinetics of Staurosporine on CLK4



**Data generated and provided by Carna Biosciences, Inc.**