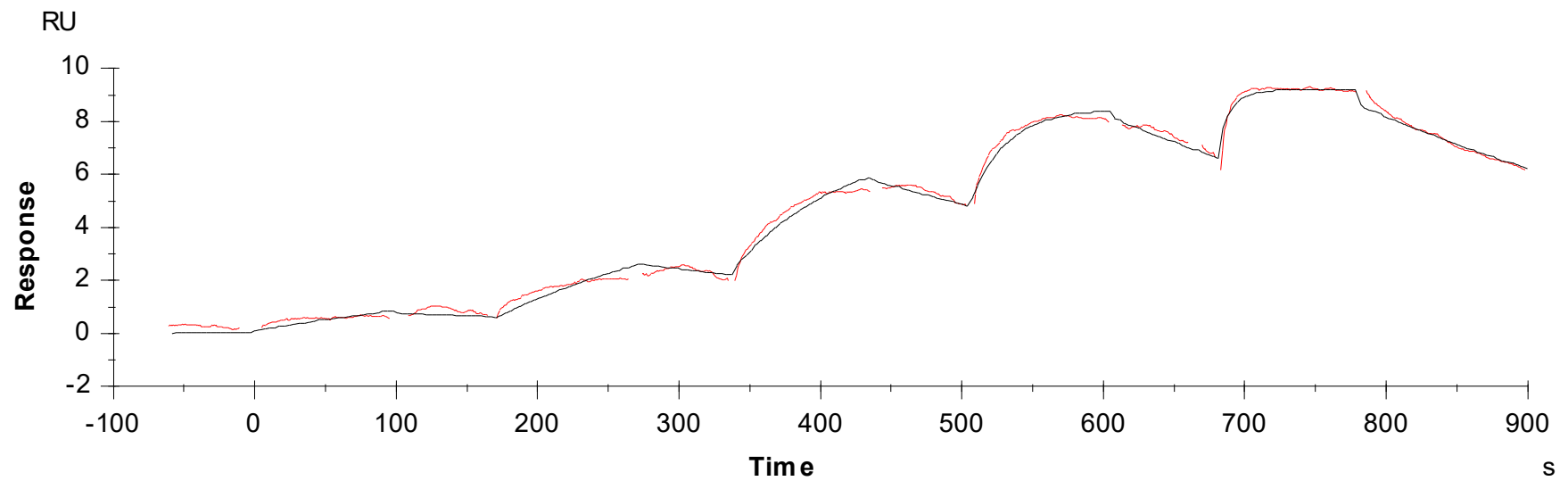


## Binding Kinetics of Staurosporine on PRKD3(PKD3)



Data generated and provided by Carna Biosciences, Inc.