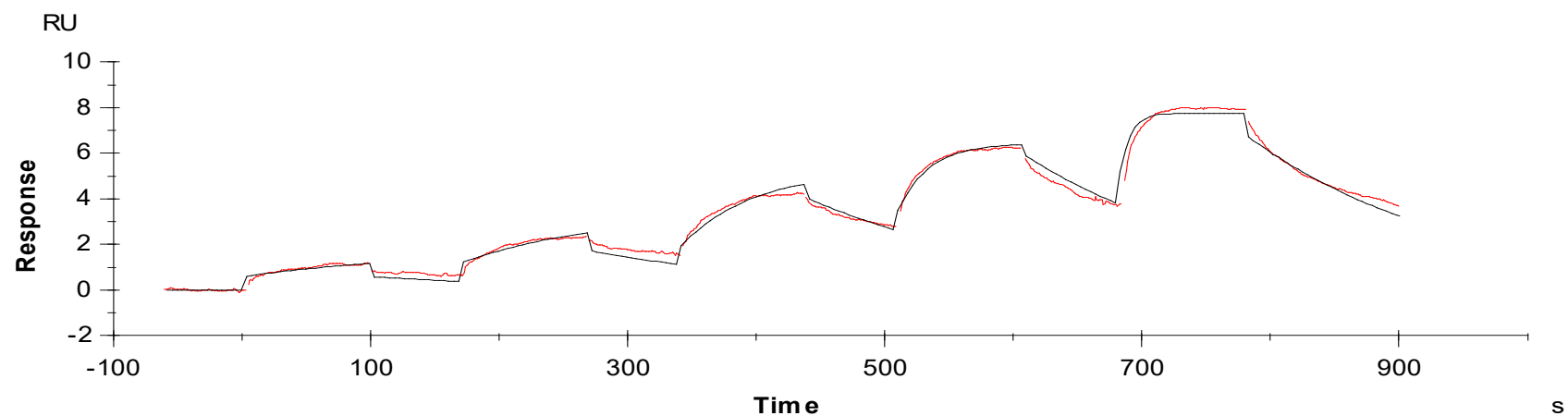


Binding Kinetics of Staurosporine on PKC η (PRKCH)



Data generated and provided by Carna Biosciences, Inc.