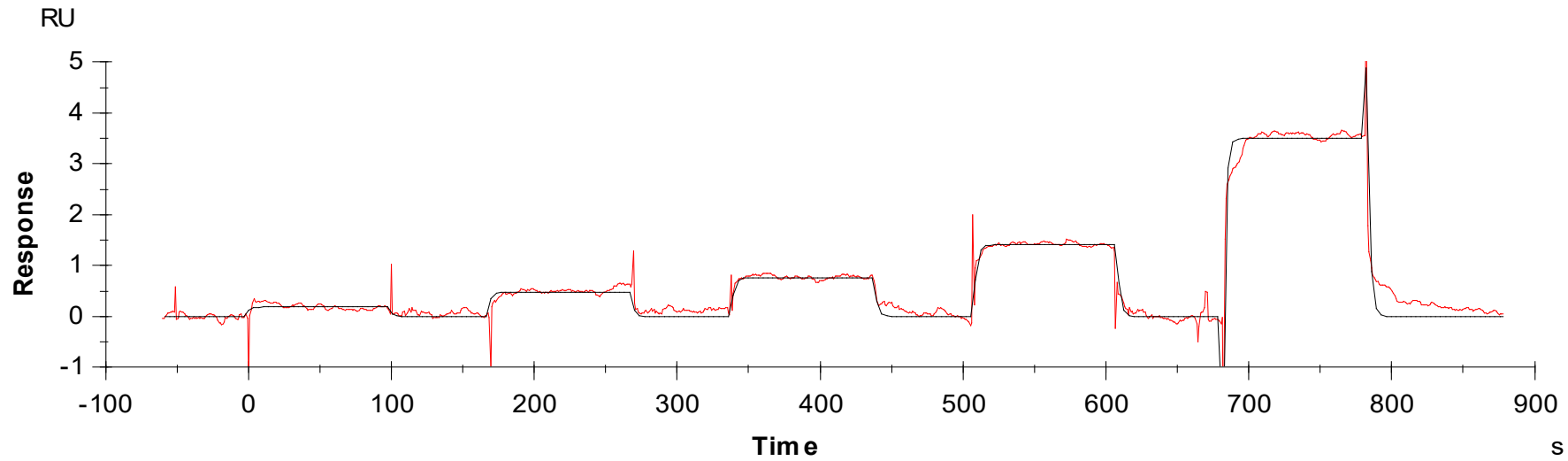




Binding Kinetics of Staurosporine on ERBB3(HER3)



Data generated and provided by Carna Biosciences, Inc.