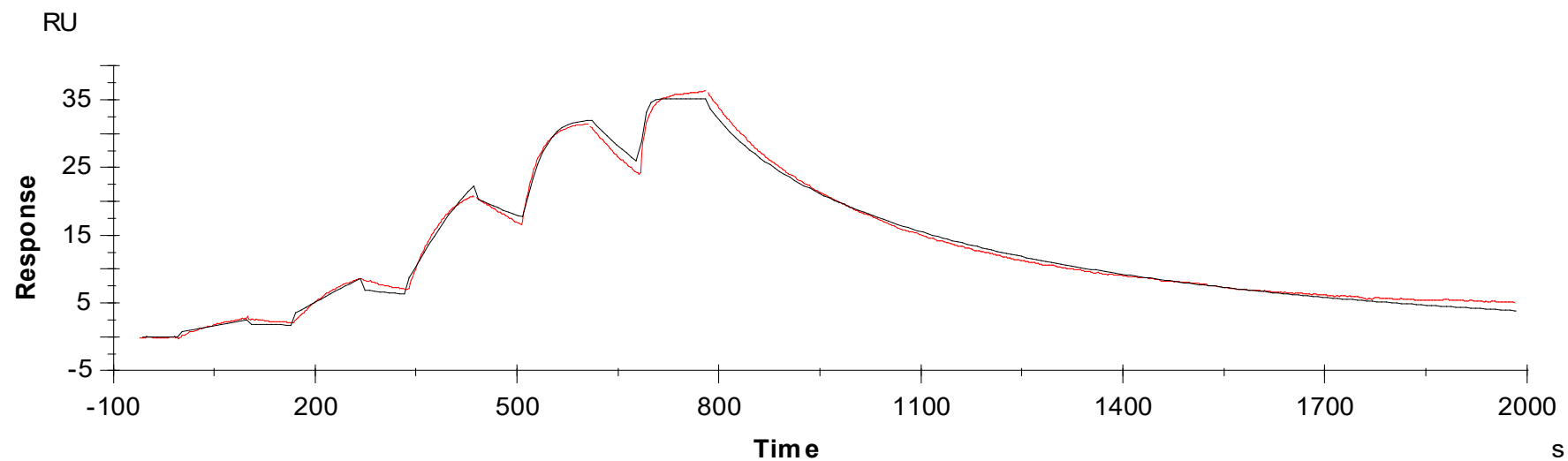


Binding Kinetics of Staurosporine on MUSK(on-activated)



Data generated and provided by Carna Biosciences, Inc.