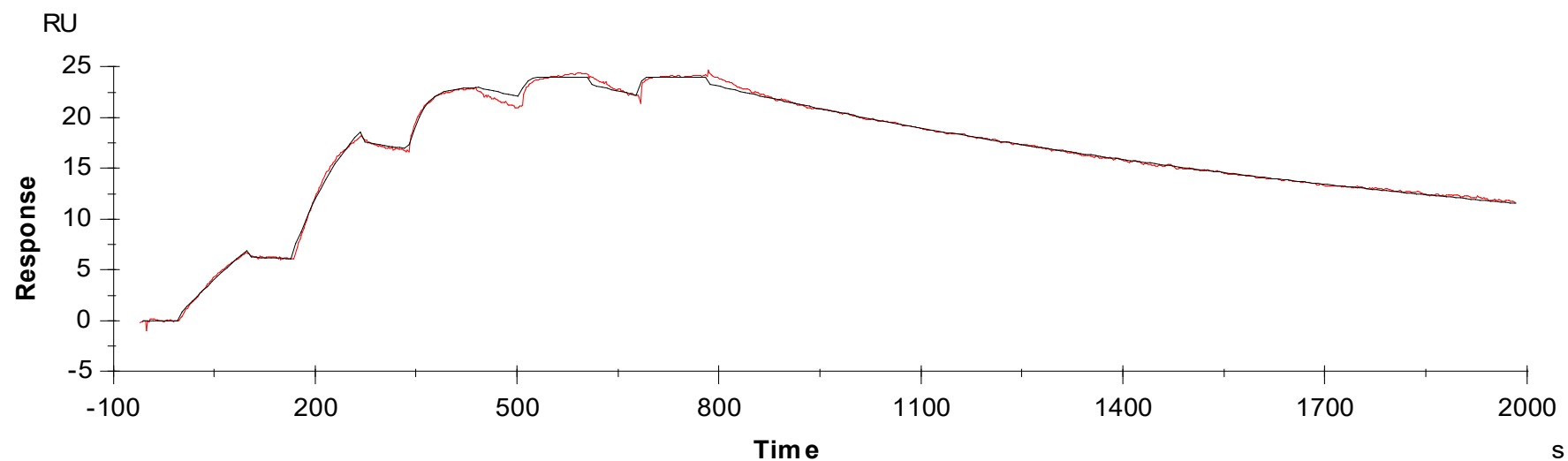


Binding Kinetics of Staurosporine on MUSK



Data generated and provided by Carna Biosciences, Inc.