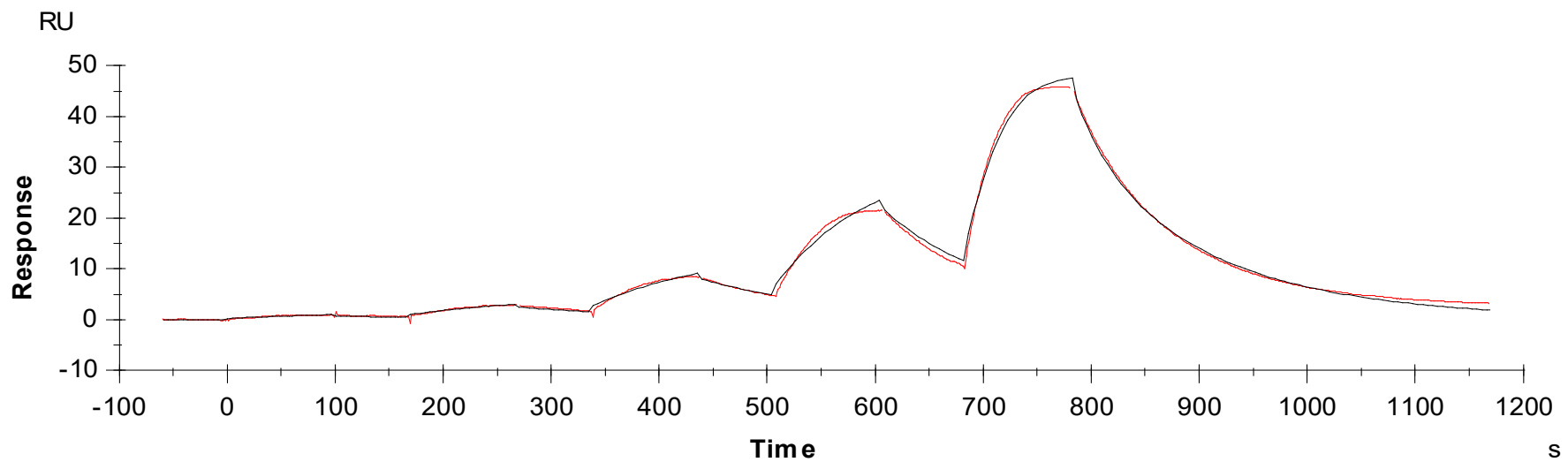




Binding Kinetics of Staurosporine on MLKL



Data generated and provided by Carna Biosciences, Inc.