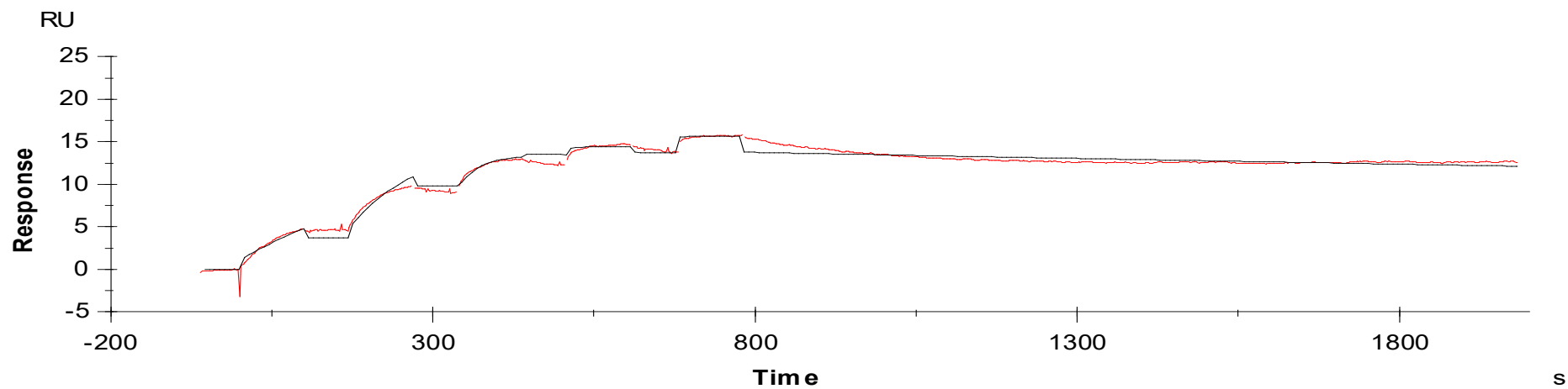


Binding Kinetics of Staurosporine on LATS2/MOBKL1A



Data generated and provided by Carna Biosciences, Inc.