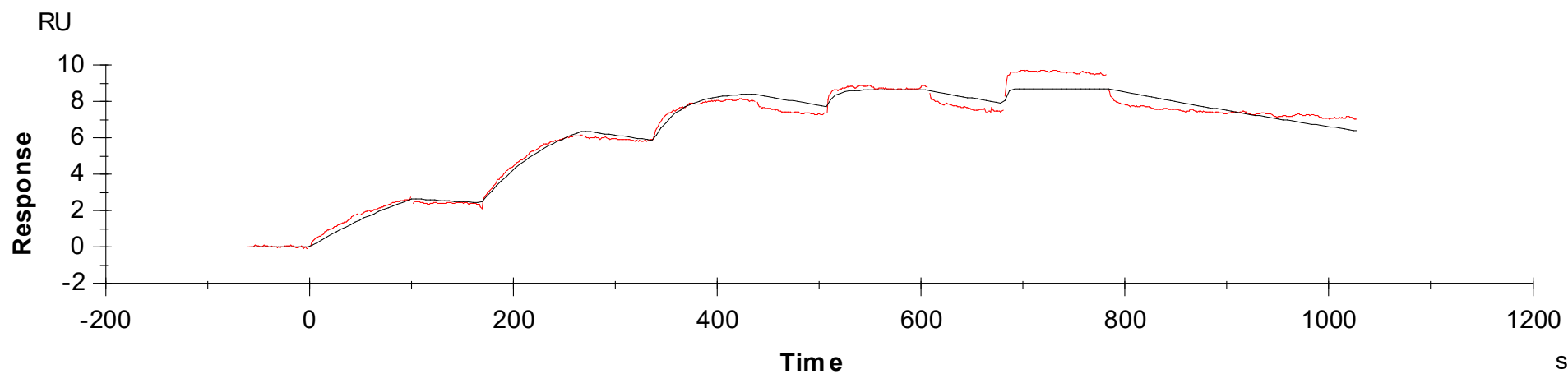


Binding Kinetics of Staurosporine on TBK1



Data generated and provided by Carna Biosciences, Inc.