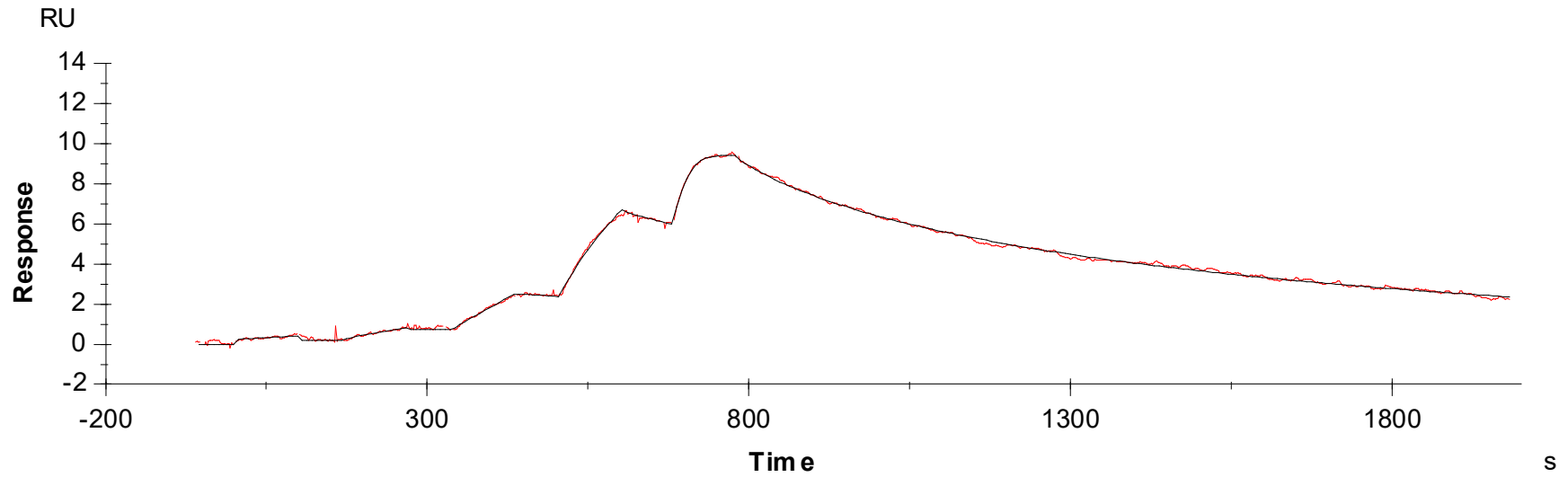


## Binding Kinetics of Staurosporine on RET[M918T](non-activated)



Data generated and provided by Carna Biosciences, Inc.