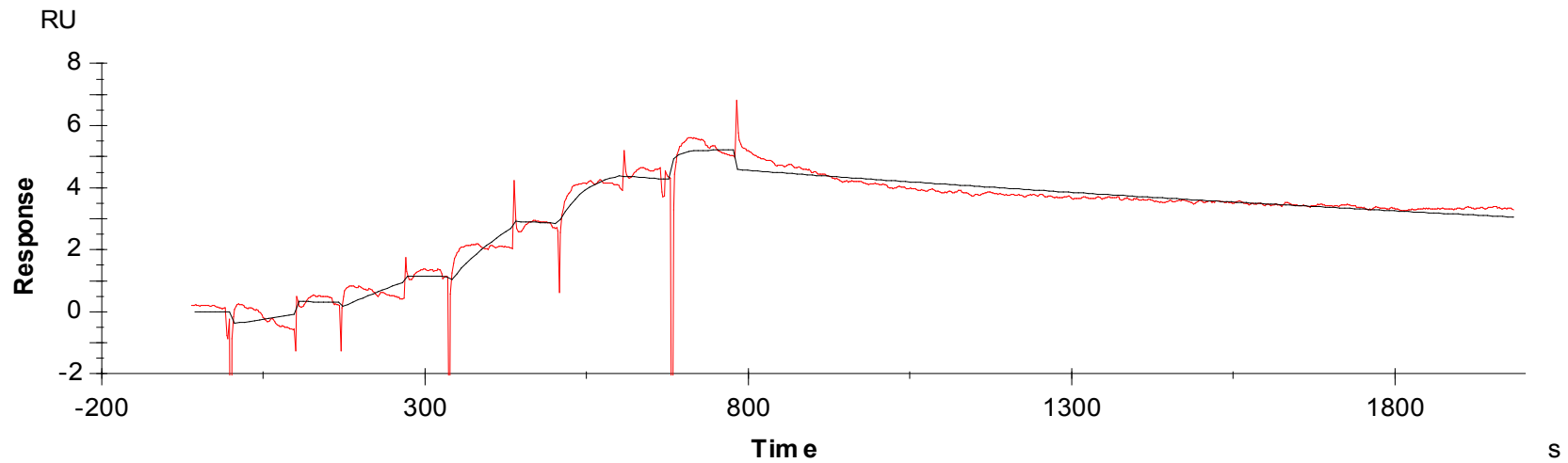


Binding Kinetics of staurosporine on FLT3



Data generated and provided by Carna Biosciences, Inc.