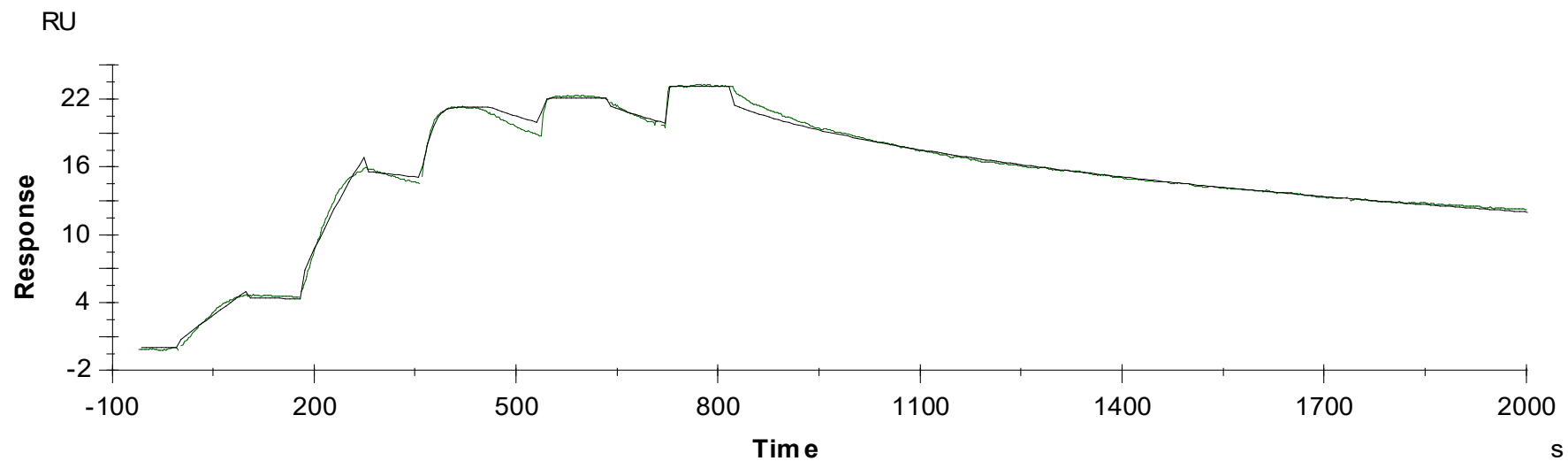


Binding Kinetics of Staurosporine on FGFR2[N549K/V564F]



Data generated and provided by Carna Biosciences, Inc.