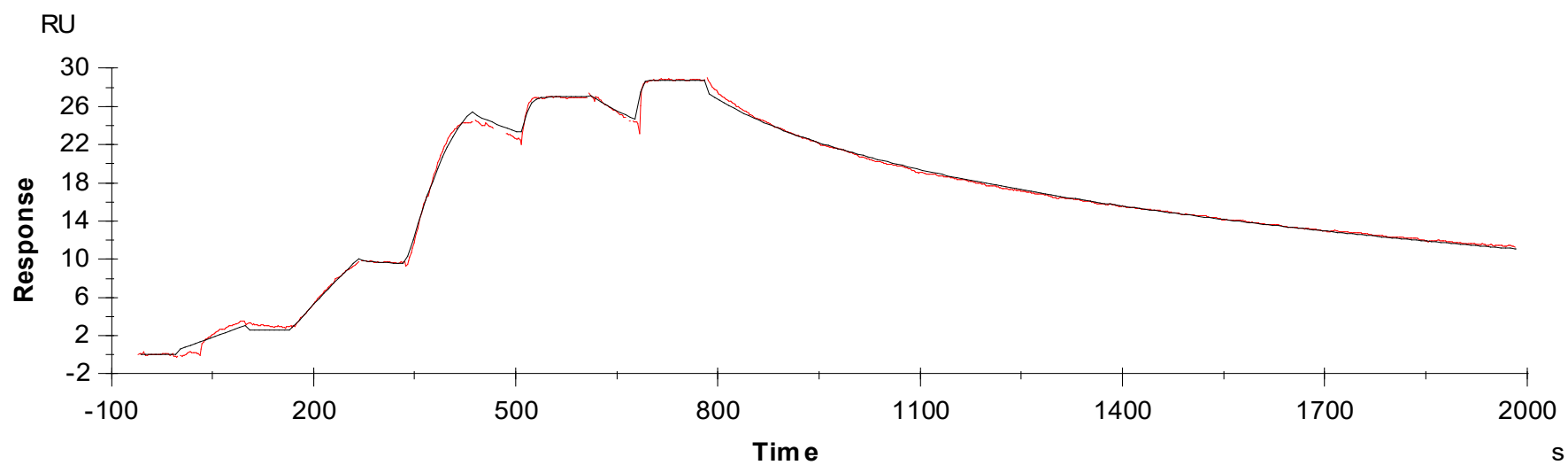


Binding Kinetics of Staurosporine on DAPK3



Data generated and provided by Carna Biosciences, Inc.