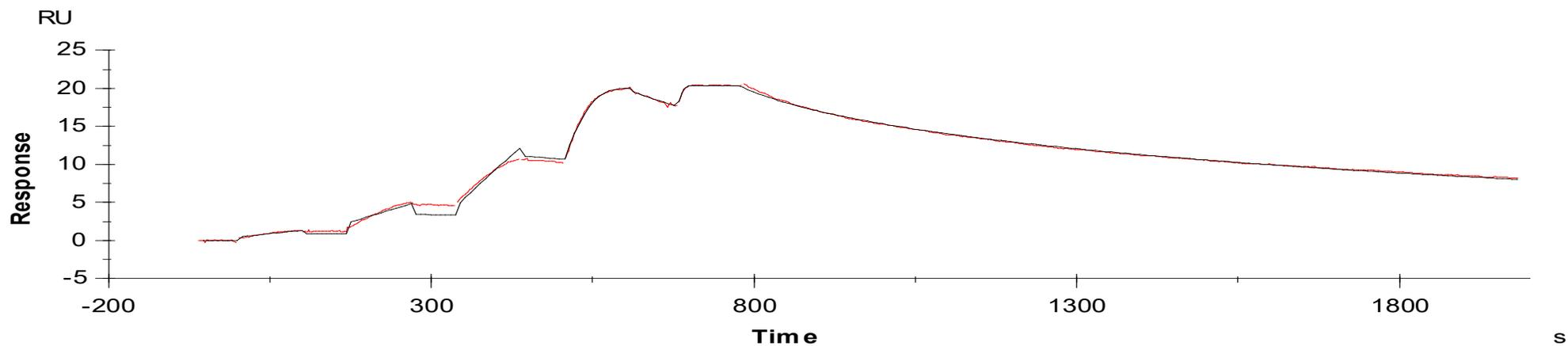


Binding Kinetics of Staurosporine on PAK6



Data generated and provided by Carna Biosciences, Inc.