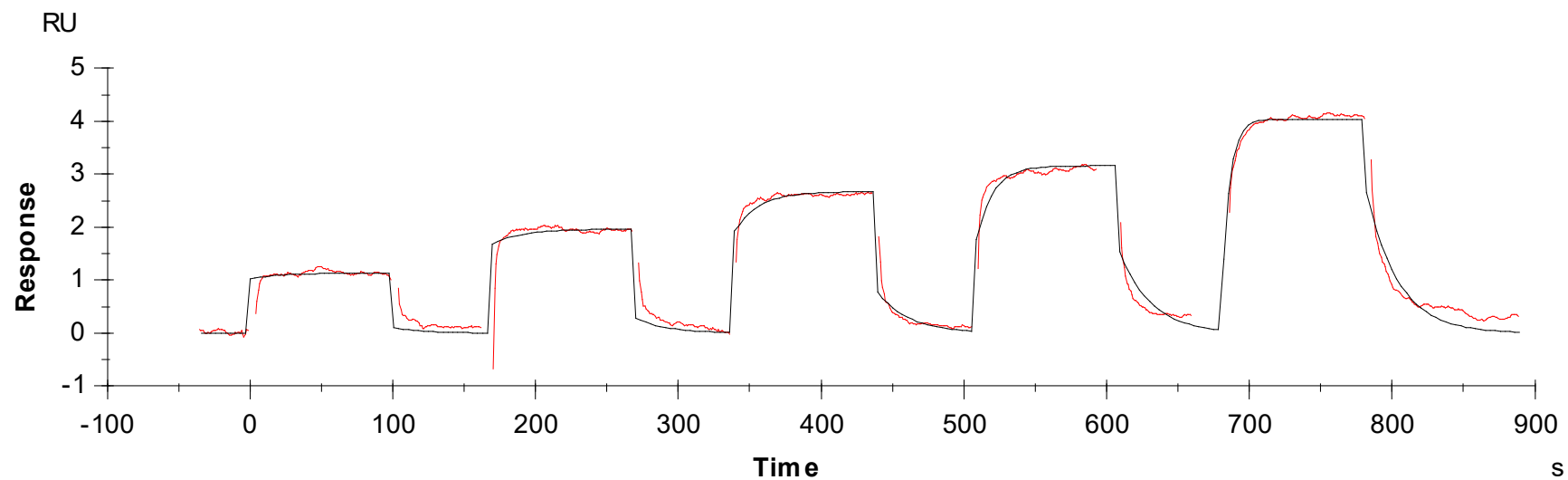


Binding Kinetics of CTx-0294885 on NEK1



Data generated and provided by Carna Biosciences, Inc.