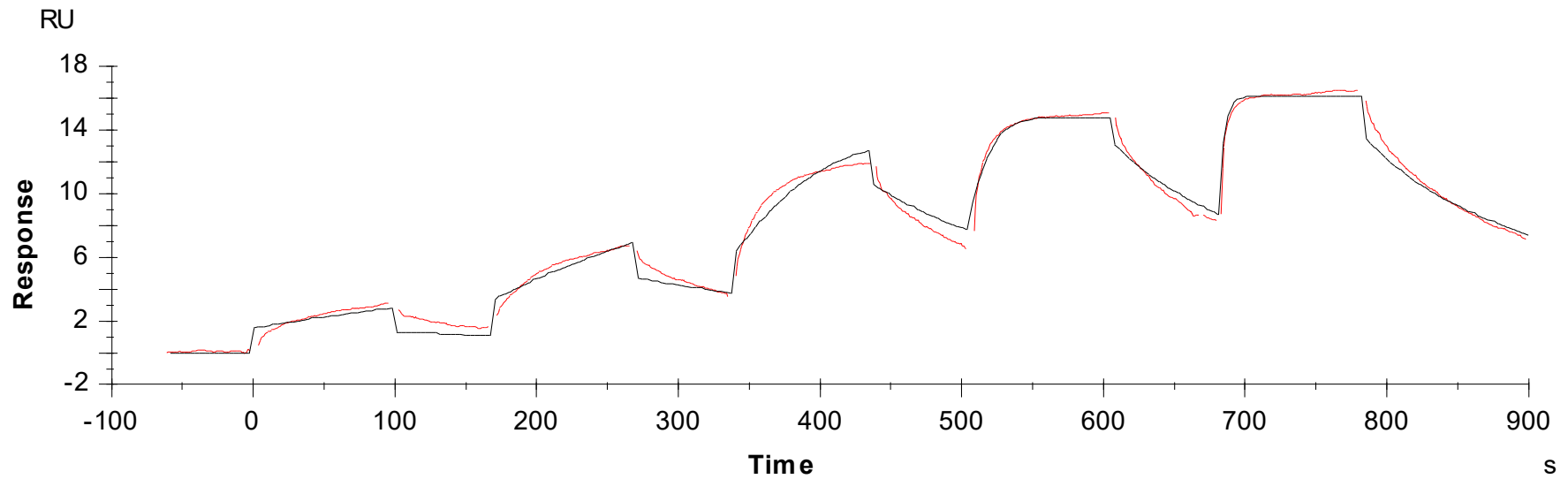


## Binding Kinetics of WZ4003 on NuaK1



Data generated and provided by Carna Biosciences, Inc.