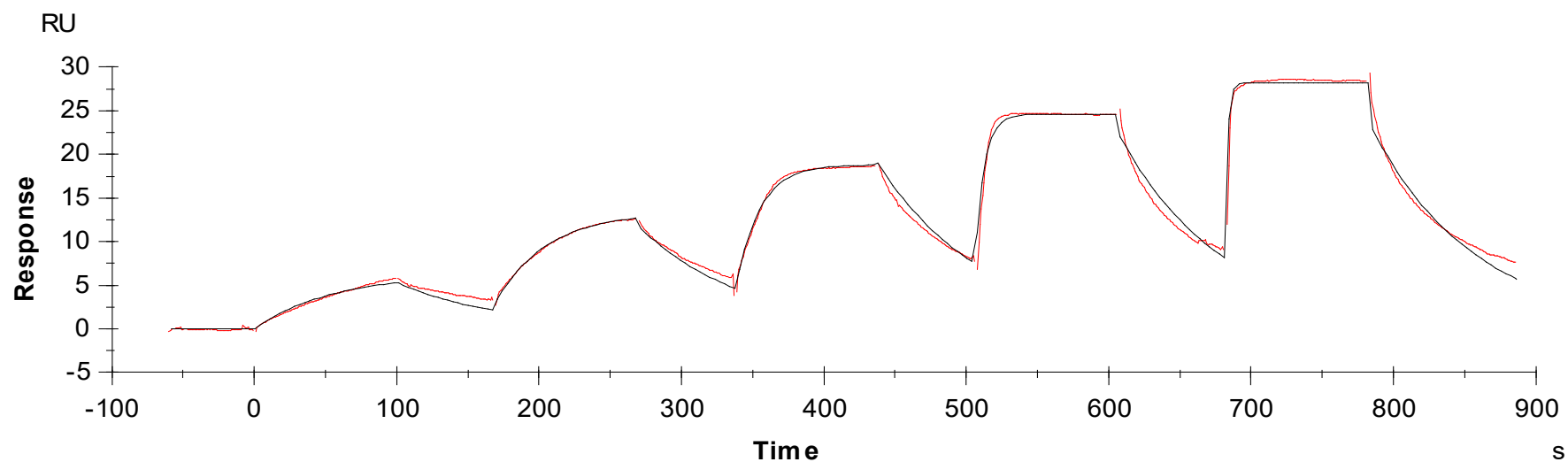


## Binding Kinetics of sunitinib on KIT[non-activated]



Data generated and provided by Carna Biosciences, Inc.