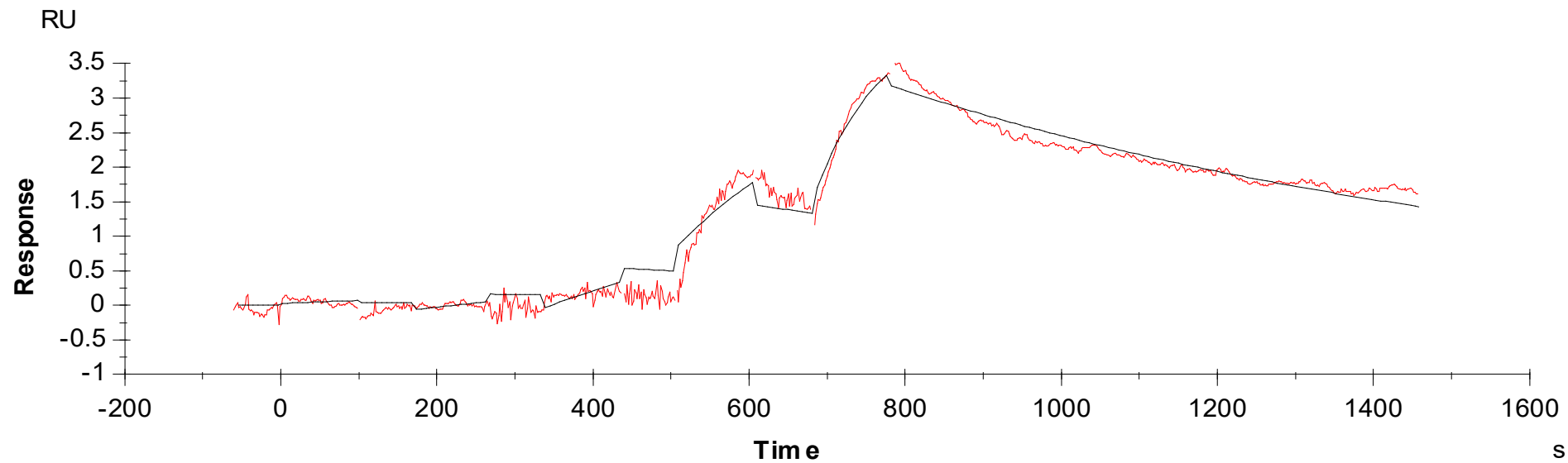


## Binding Kinetics of Compound X (Carna's proprietary) on DGK $\alpha$ (196-735aa)



Data generated and provided by Carna Biosciences, Inc.