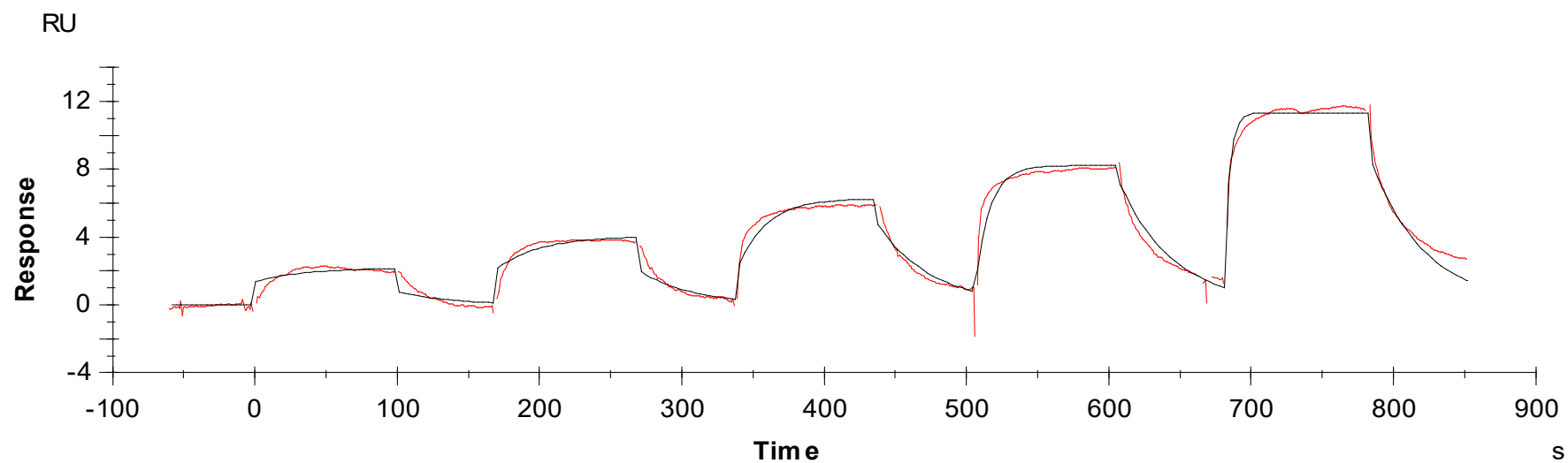


Binding Kinetics of K252A on PKC γ (PRKCG)



Data generated and provided by Carna Biosciences, Inc.