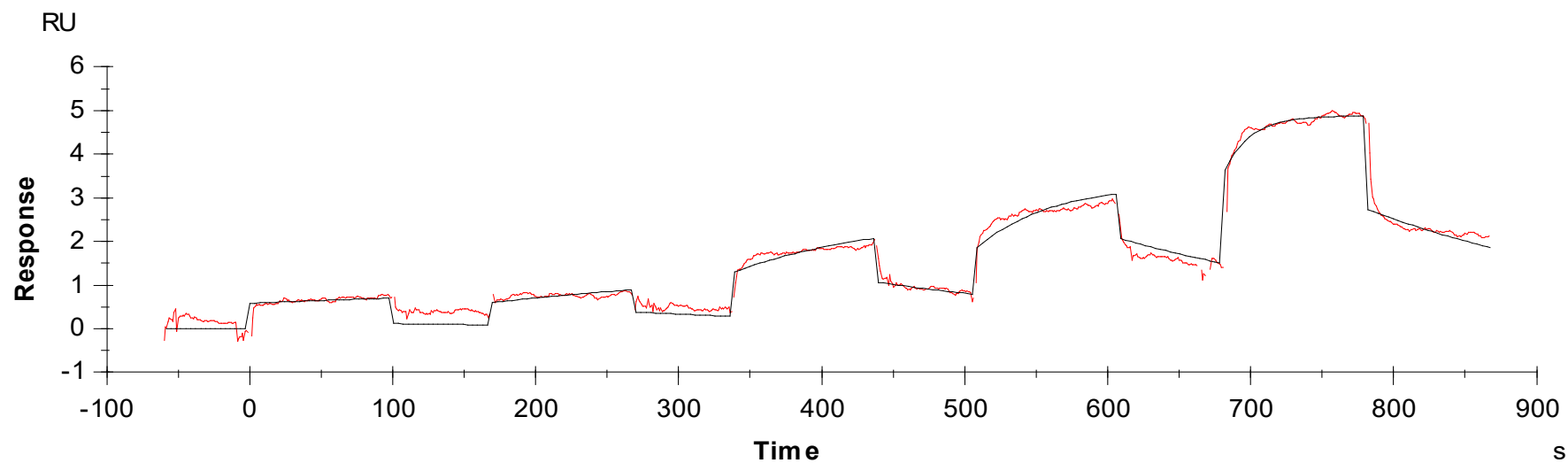




Binding Kinetics of K252A on DLK(MAP3K12)



Data generated and provided by Carna Biosciences, Inc.