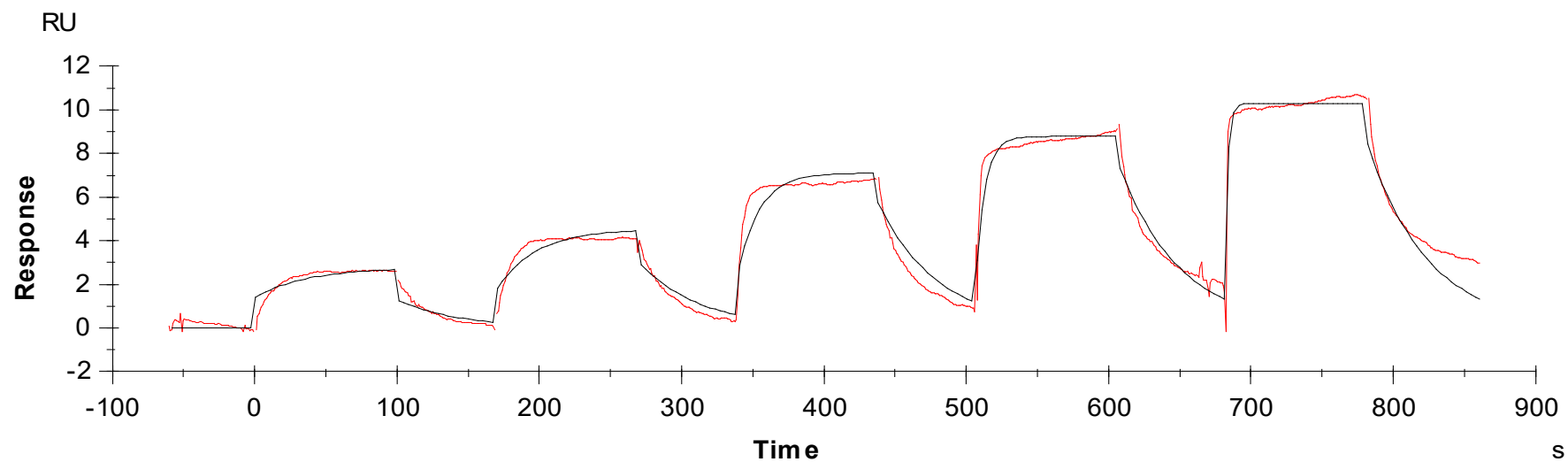


Binding Kinetics of K252A on MNK1(MKNK1)



Data generated and provided by Carna Biosciences, Inc.