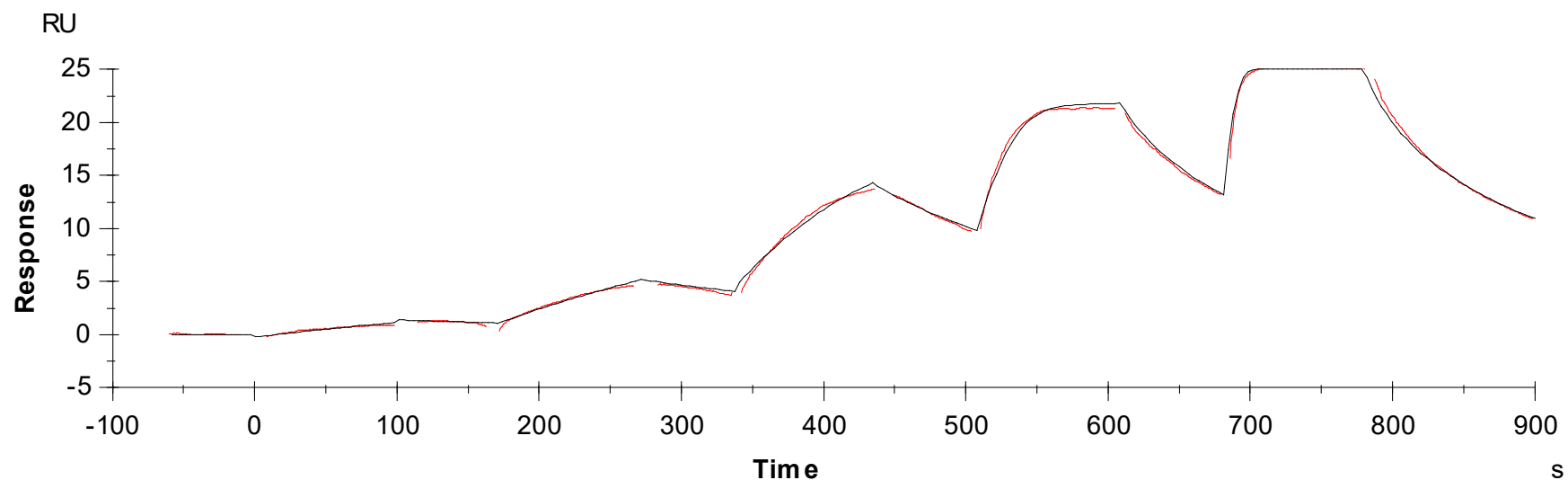


Binding Kinetics of MK-2 inhibitor III on MAPKAPK2



Data generated and provided by Carna Biosciences, Inc.