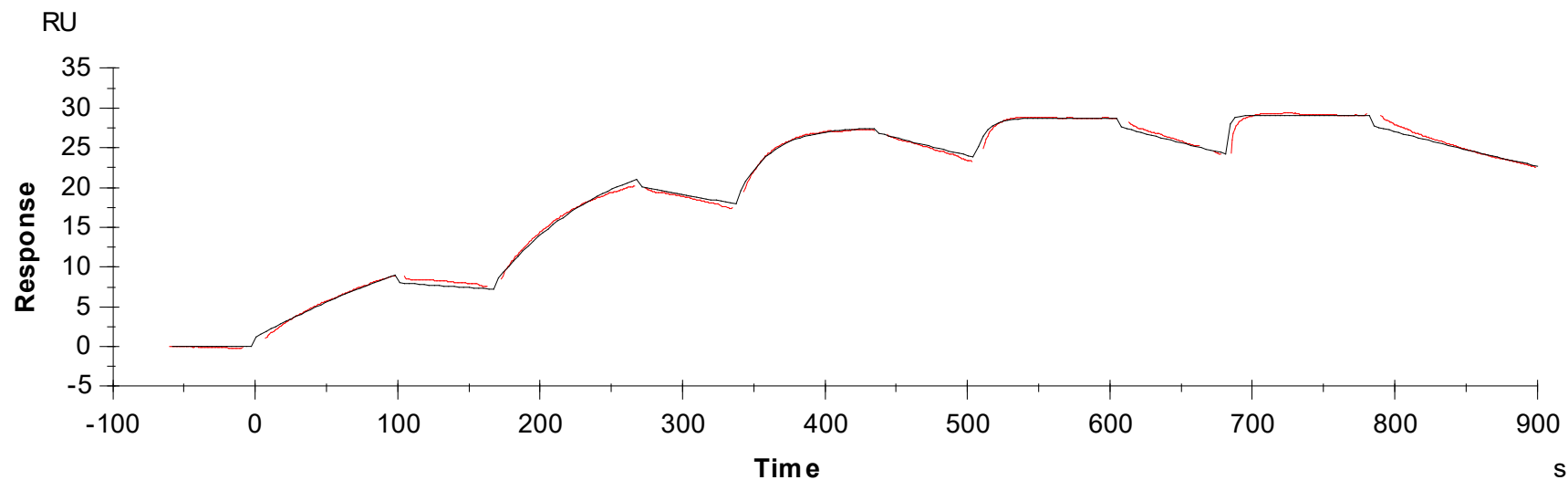


Binding Kinetics of Midostaurin on CAMK2 α (CAMK2A)



Data generated and provided by Carna Biosciences, Inc.