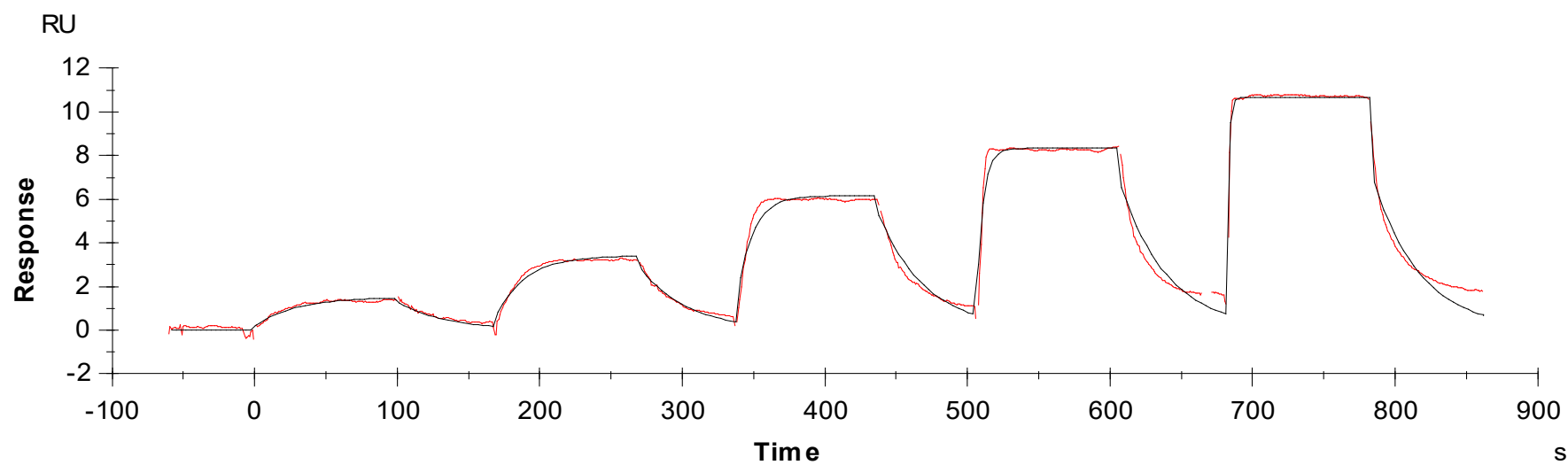


Binding Kinetics of sunitinib on DDR1(SRC treated)



Data generated and provided by Carna Biosciences, Inc.