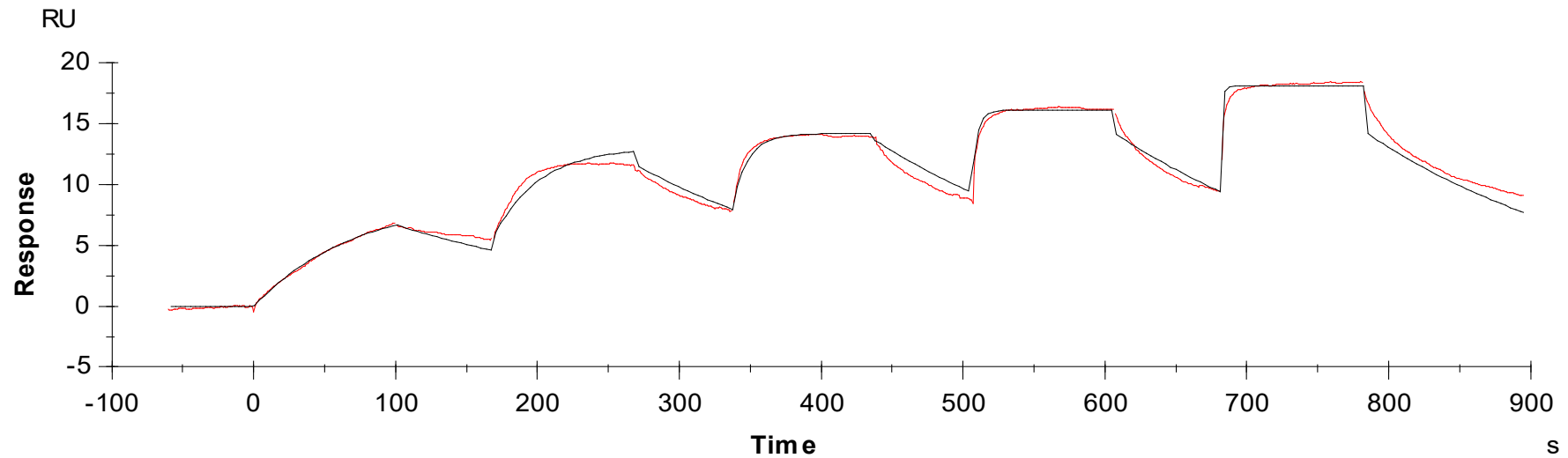


Binding Kinetics of crizotinib on ROS(ROS1)



Data generated and provided by Carna Biosciences, Inc.