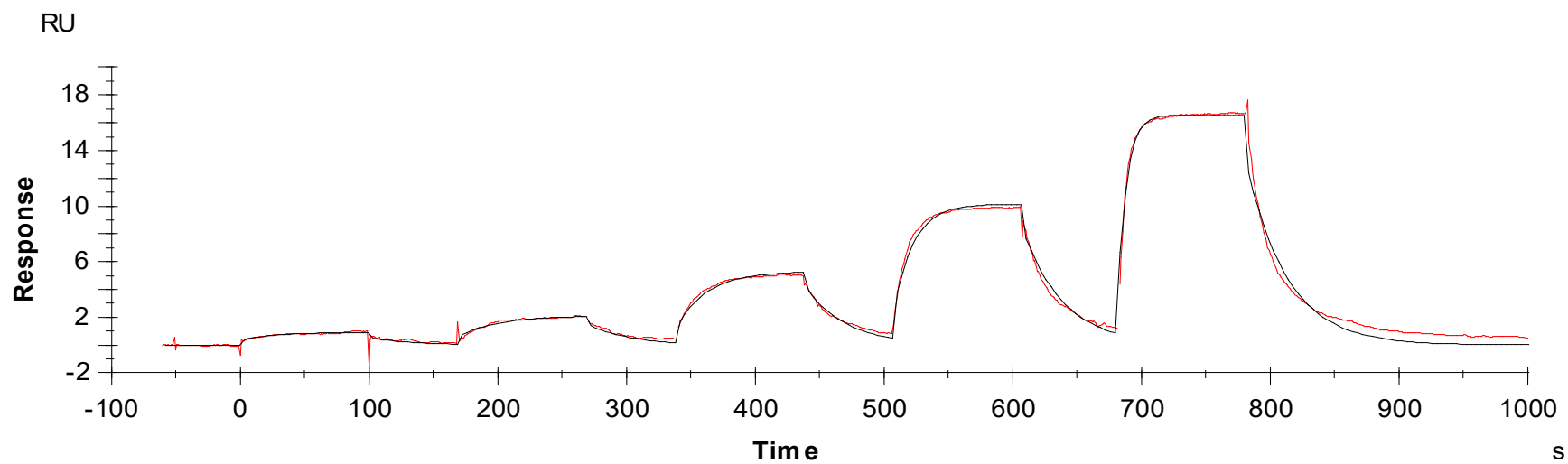




Binding Kinetics of Ribociclib on CDK6/CycD3[non-activated]



Data generated and provided by Carna Biosciences, Inc.