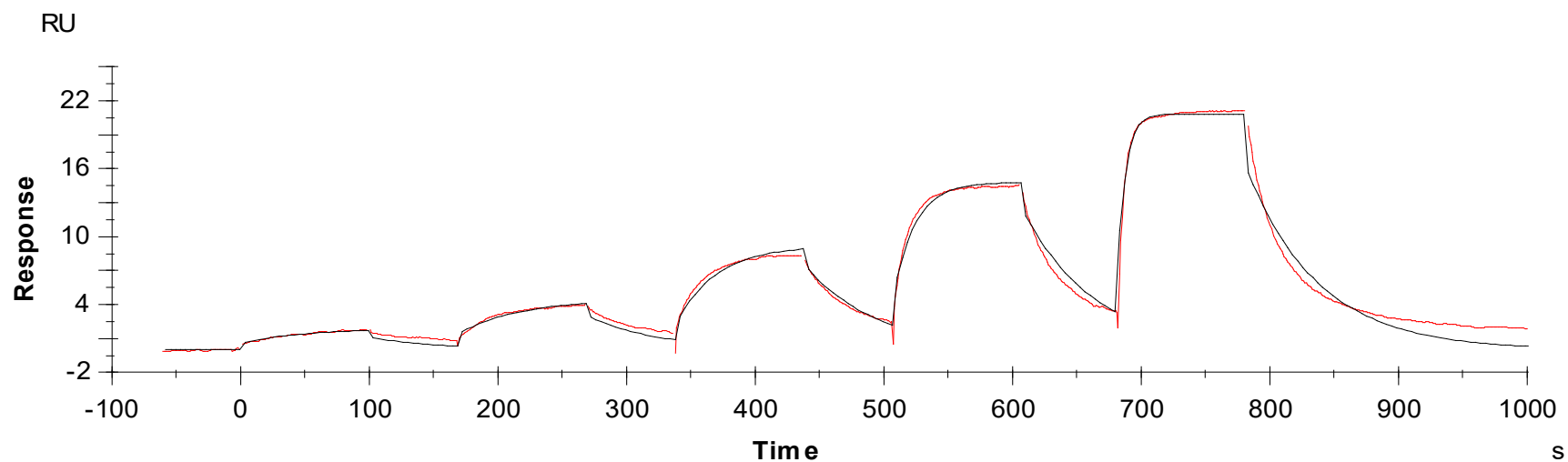




Binding Kinetics of Ribociclib on CDK6/CycD3



Data generated and provided by Carna Biosciences, Inc.