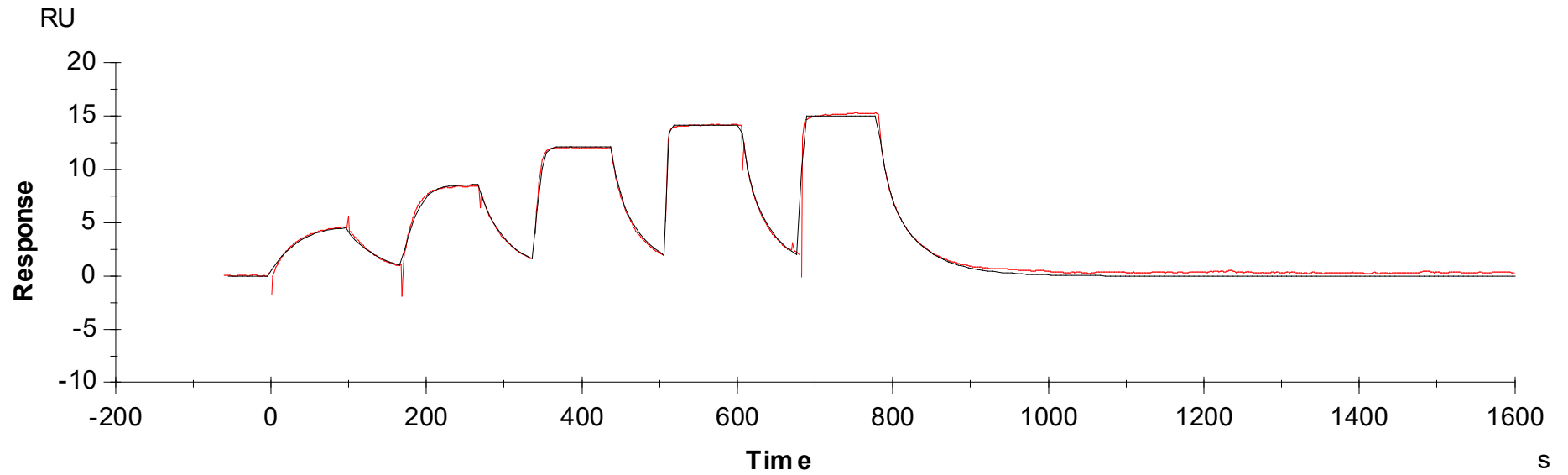




Binding Kinetics of AT7519 on CDK1(CDC2)/cyclinB1



Data generated and provided by Carna Biosciences, Inc.