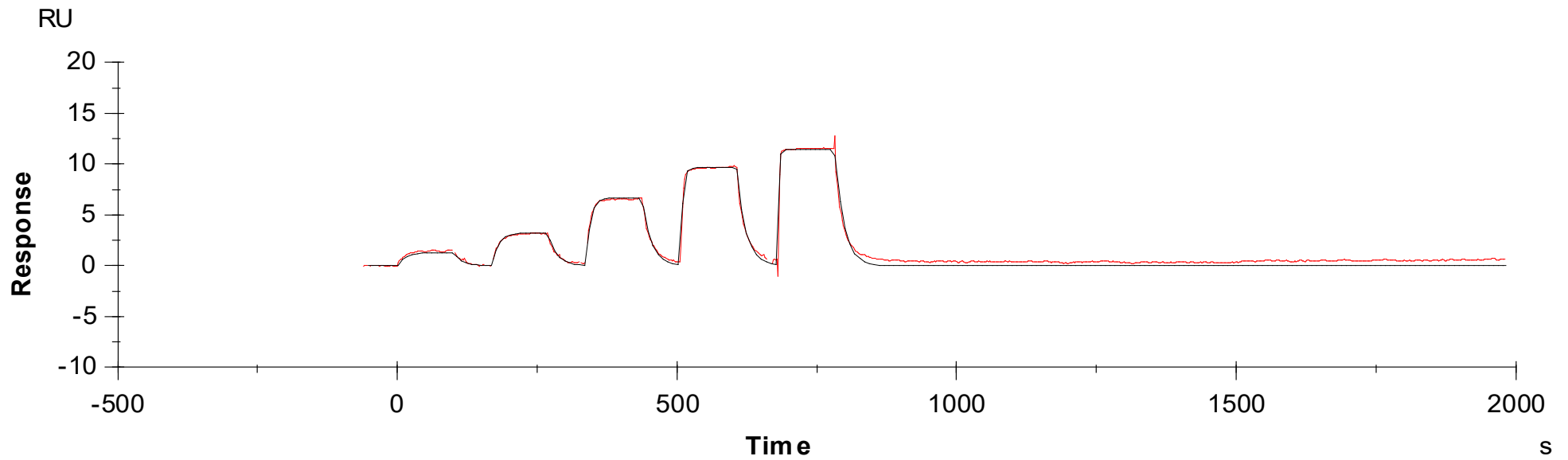


Binding Kinetics of AT7519 on CDK3/cyclinE1



Data generated and provided by Carna Biosciences, Inc.