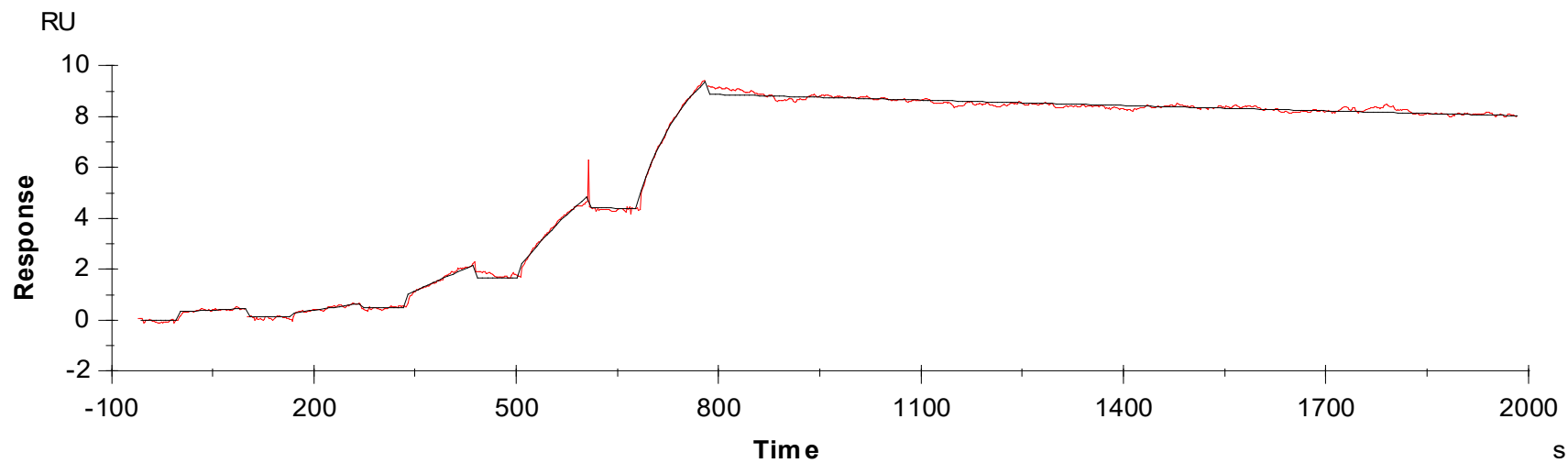




## Binding Kinetics of Ponatinib on CDK19(CDC2L6) / CycC



Data generated and provided by Carna Biosciences, Inc.