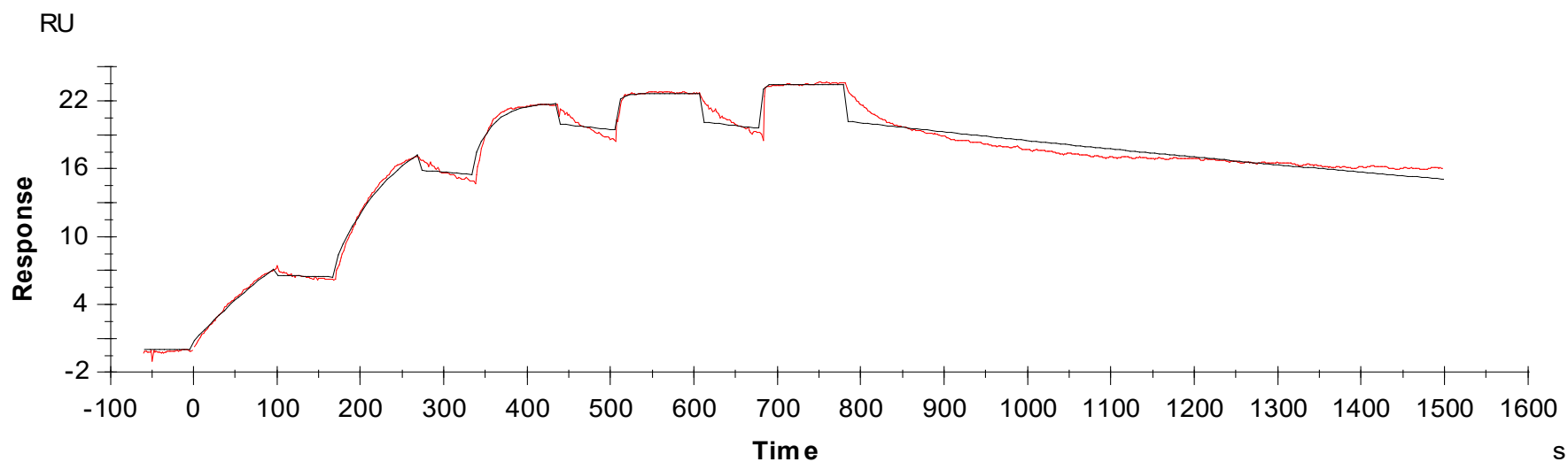




Binding Kinetics of Dinaciclib on CDK2/CycE1



Data generated and provided by Carna Biosciences, Inc.

Binding kinetics of SNS-032 on CDK2/CycE1

