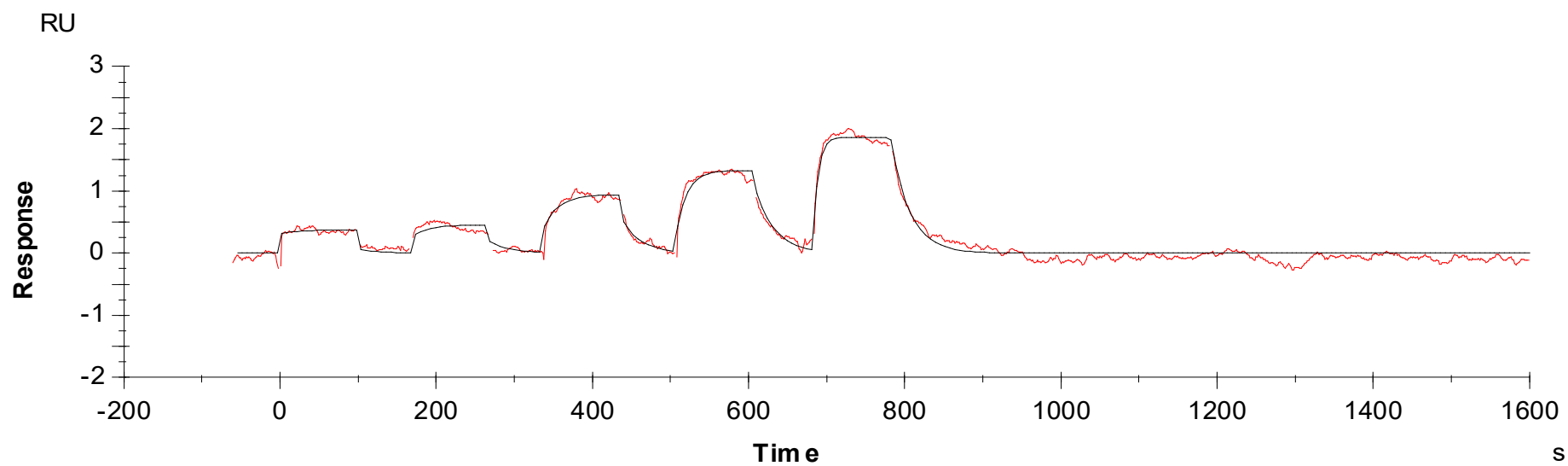


## Binding Kinetics of YM201636 on PIKFYVE



Data generated and provided by Carna Biosciences, Inc.