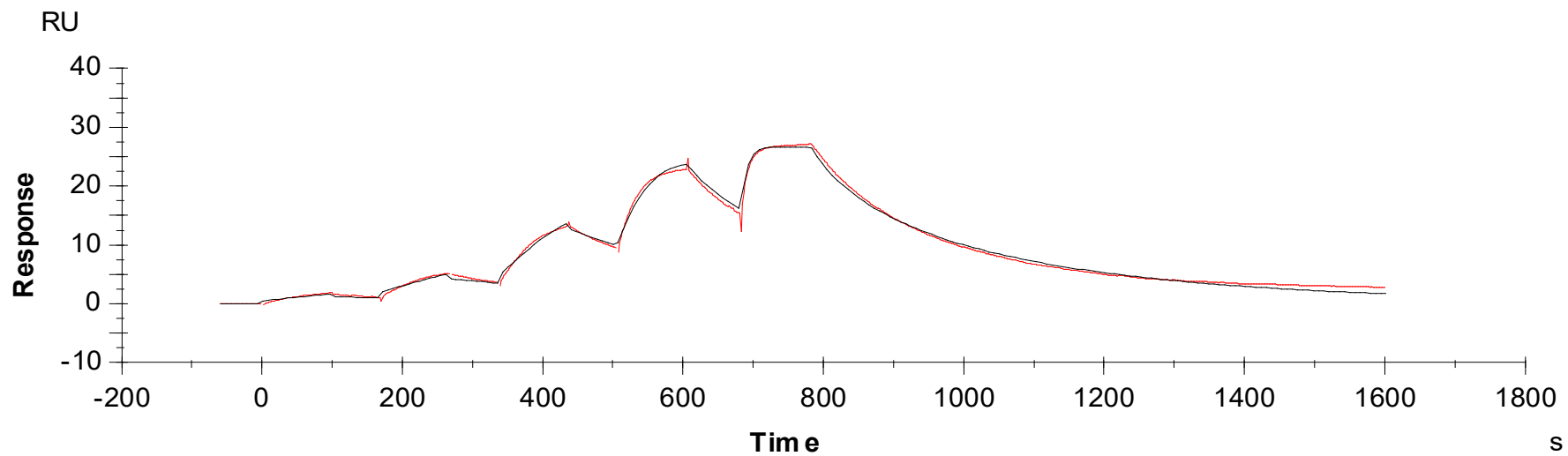


## Binding Kinetics of Cabozantinib on RET[Y791F]



Data generated and provided by Carna Biosciences, Inc.

# Binding kinetics of staurosporine on RET [Y791F]

