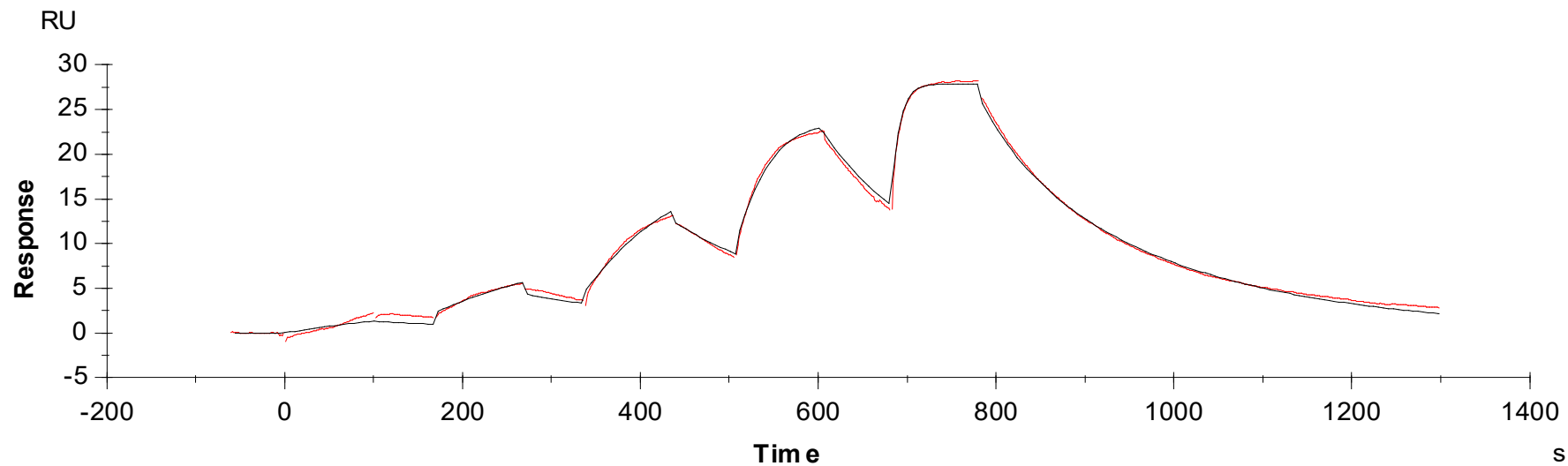


## Binding Kinetics of Cabozantinib on RET[G691S]



Data generated and provided by Carna Biosciences, Inc.

# Binding kinetics of staurosporine on RET [G691S]

