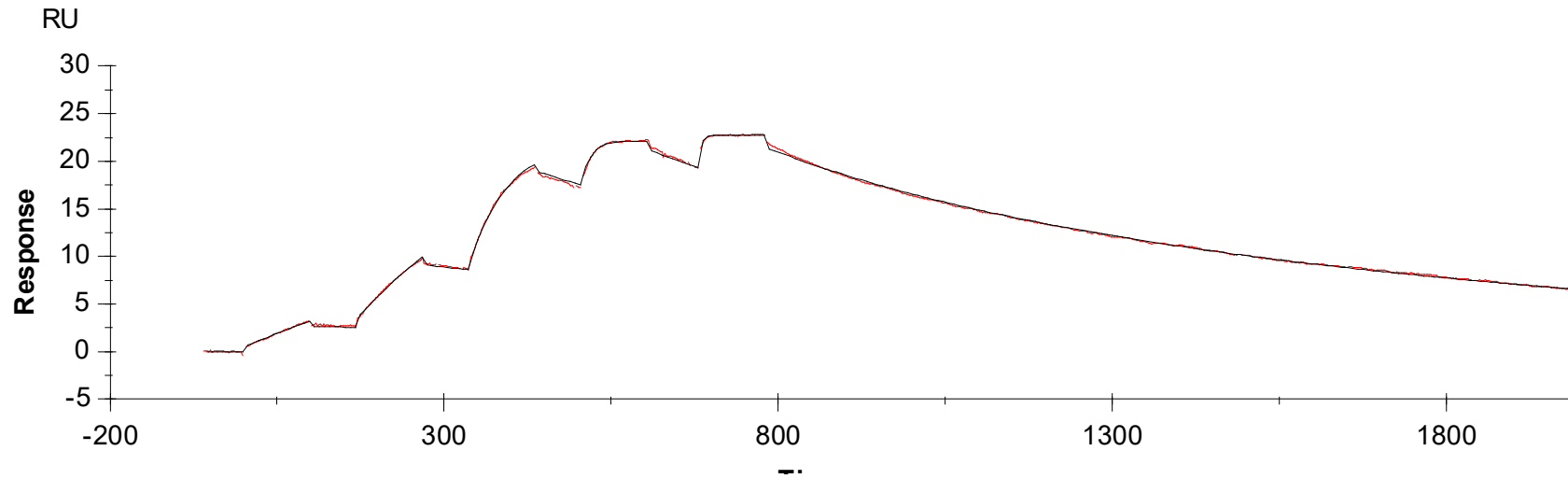


## Binding Kinetics of Cabozantinib on RET[G691S](non-activated)



Data generated and provided by Carna Biosciences, Inc.