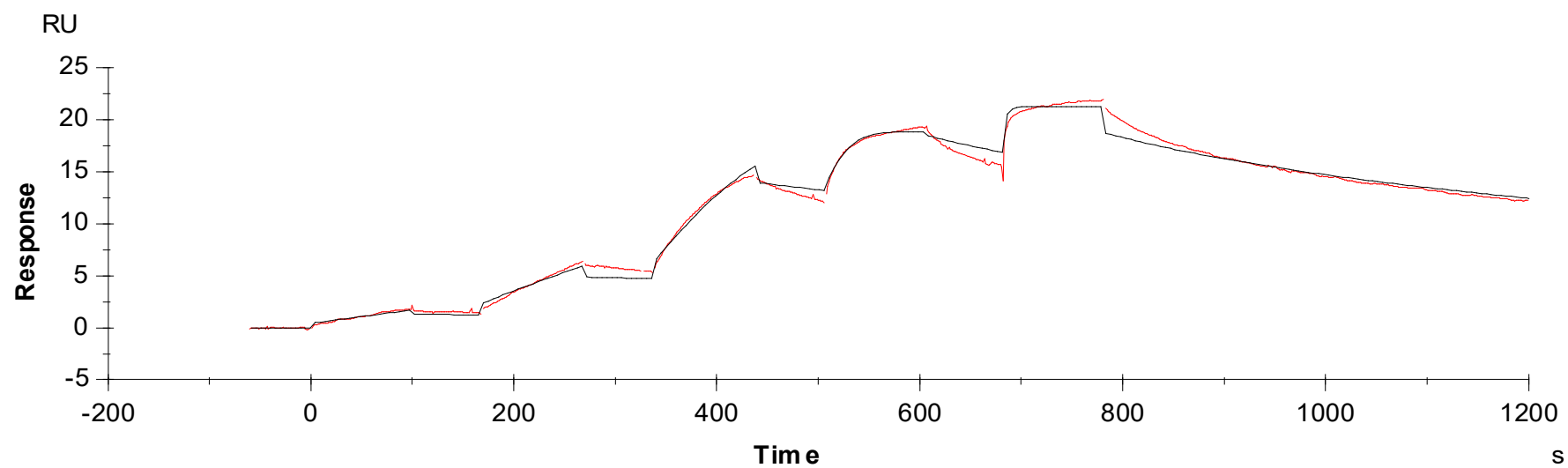


## Binding Kinetics of Brigatinib on ALK[R1275Q]



Data generated and provided by Carna Biosciences, Inc.