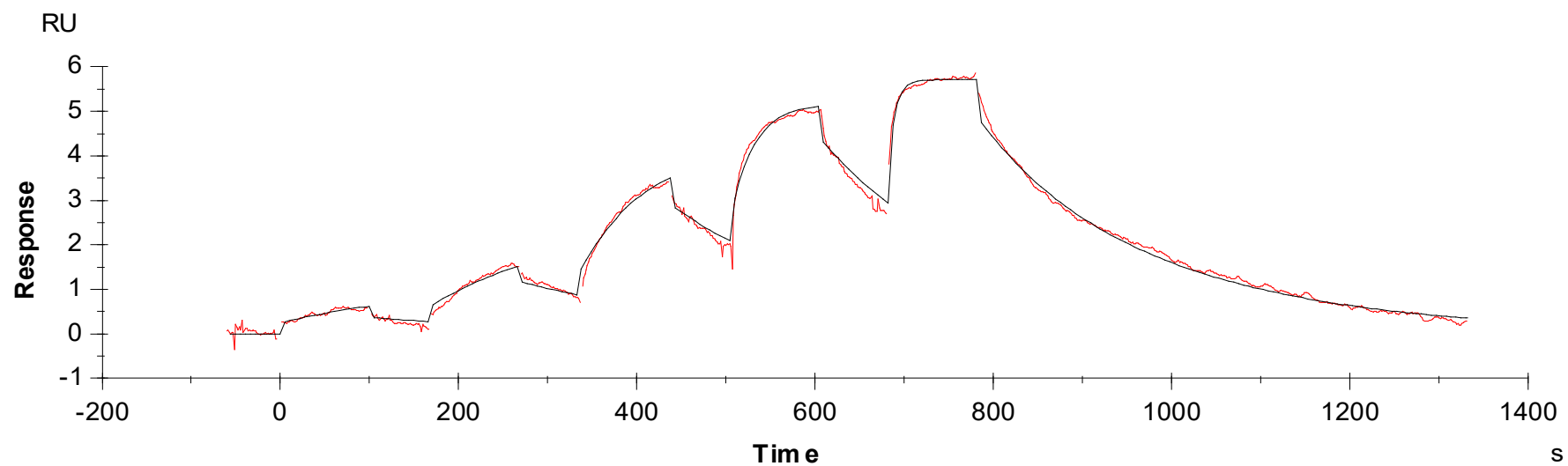


## Binding Kinetics of Brigatinib on ALK[F1174L/G1202R]



Data generated and provided by Carna Biosciences, Inc.