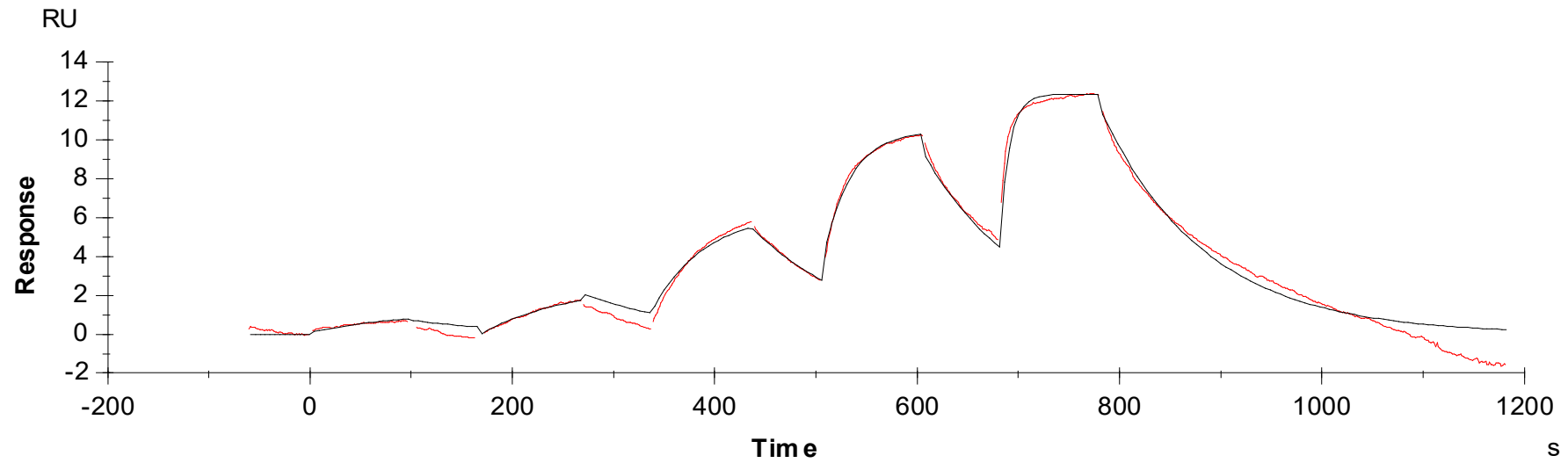


Binding Kinetics of Brigatinib on ALK[I1171N/L1198F]



Data generated and provided by Carna Biosciences, Inc.