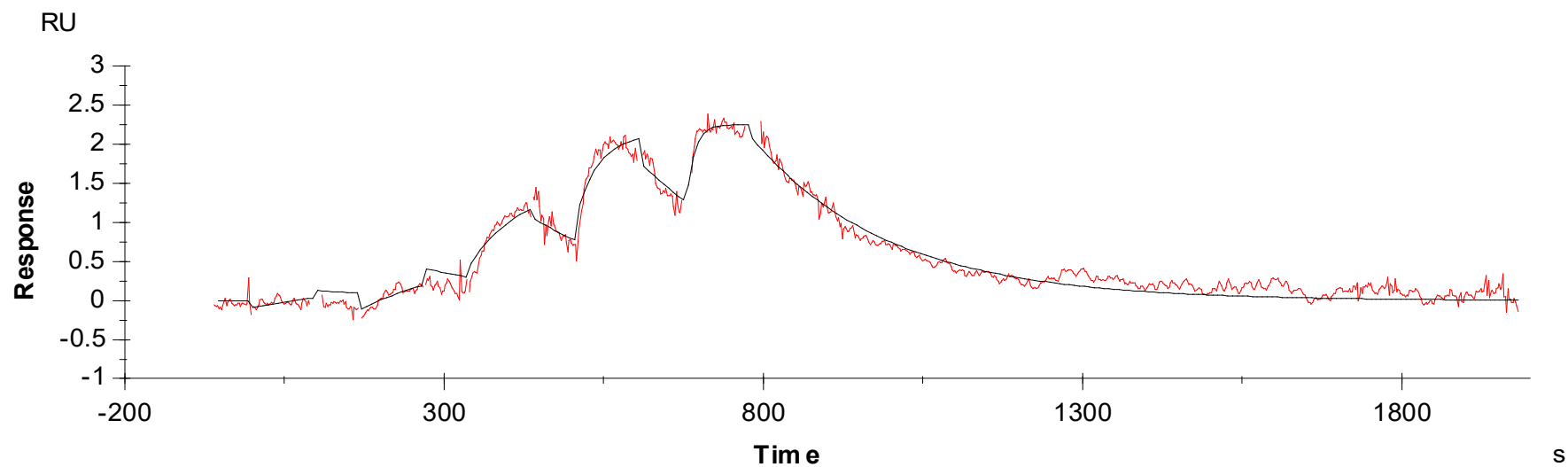


Binding Kinetics of TAE684 on ALK[L1196M/G1202R]



Data generated and provided by Carna Biosciences, Inc.