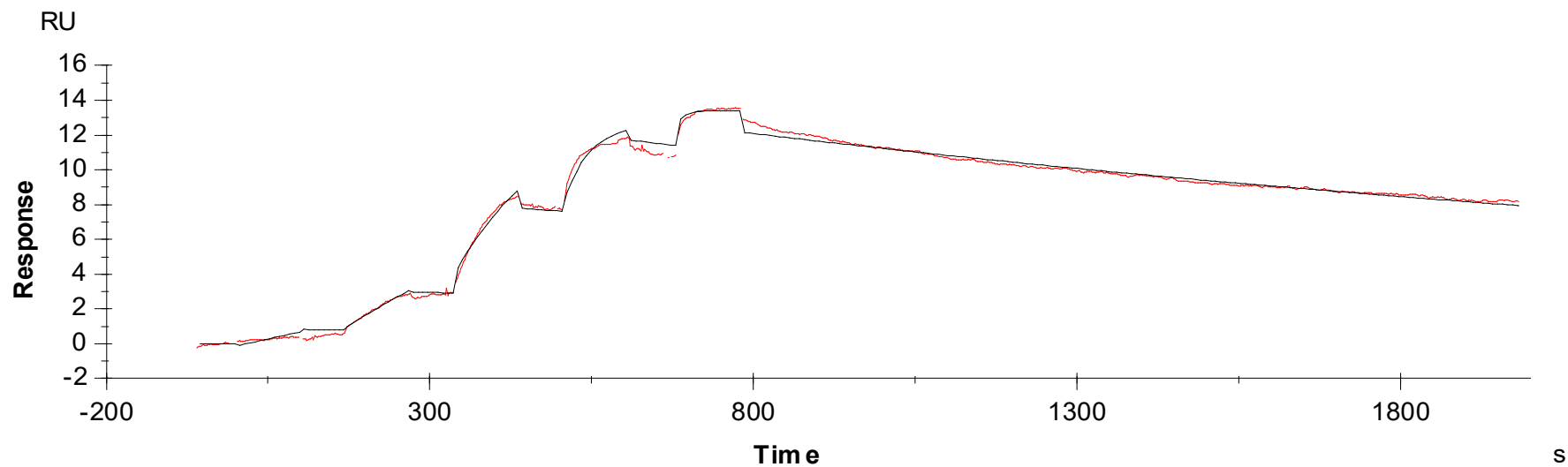


Binding Kinetics of TAE684 on ALK[G1269A]



Data generated and provided by Carna Biosciences, Inc.