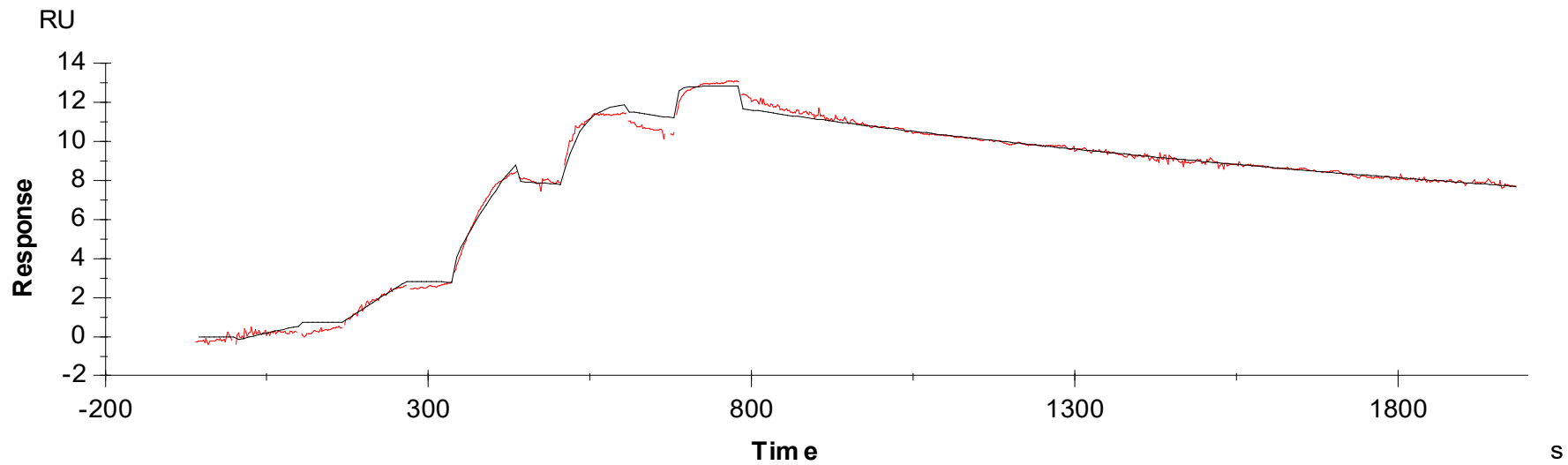


Binding Kinetics of TAE684 on ALK [V1180L]



Data generated and provided by Carna Biosciences, Inc.