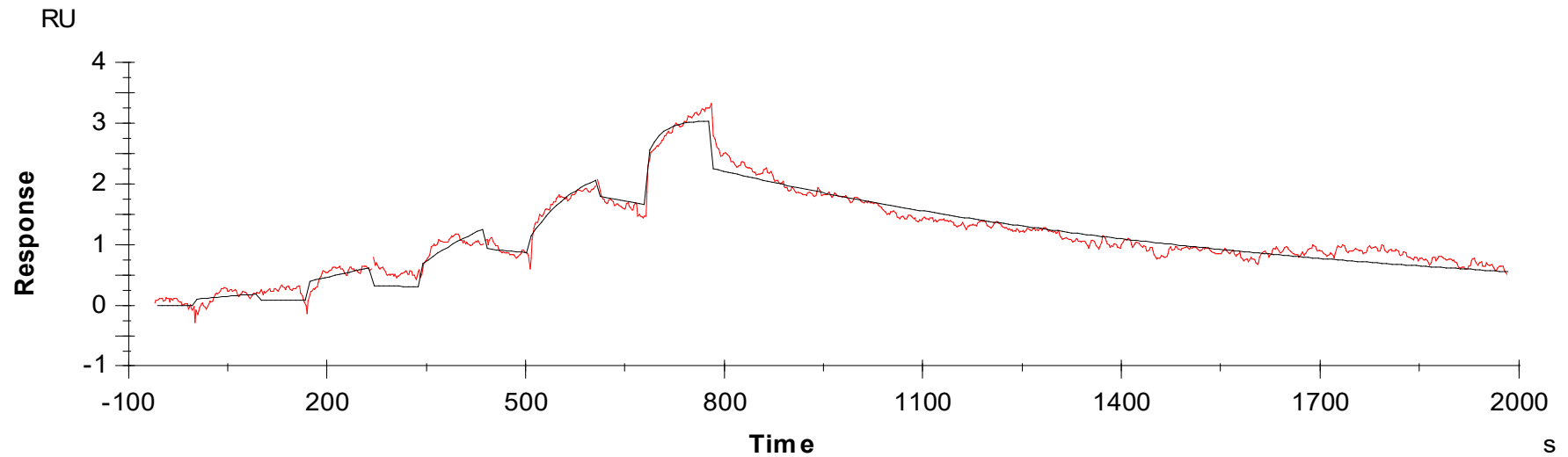


Binding Kinetics of TAE684 on ACVR1B(ALK4)



Data generated and provided by Carna Biosciences, Inc.