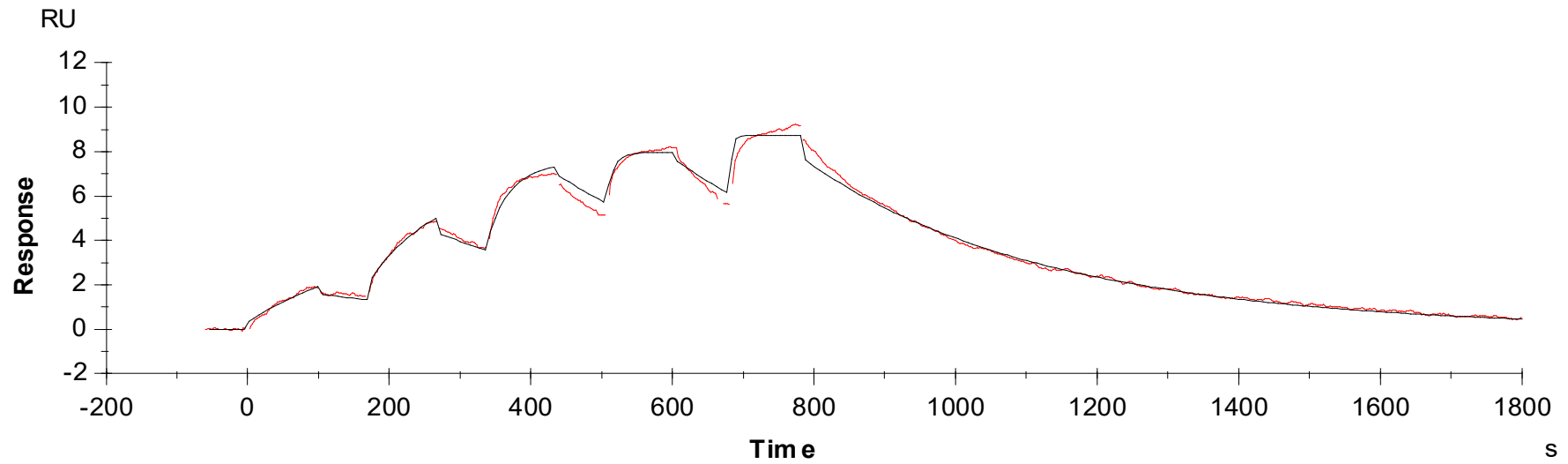
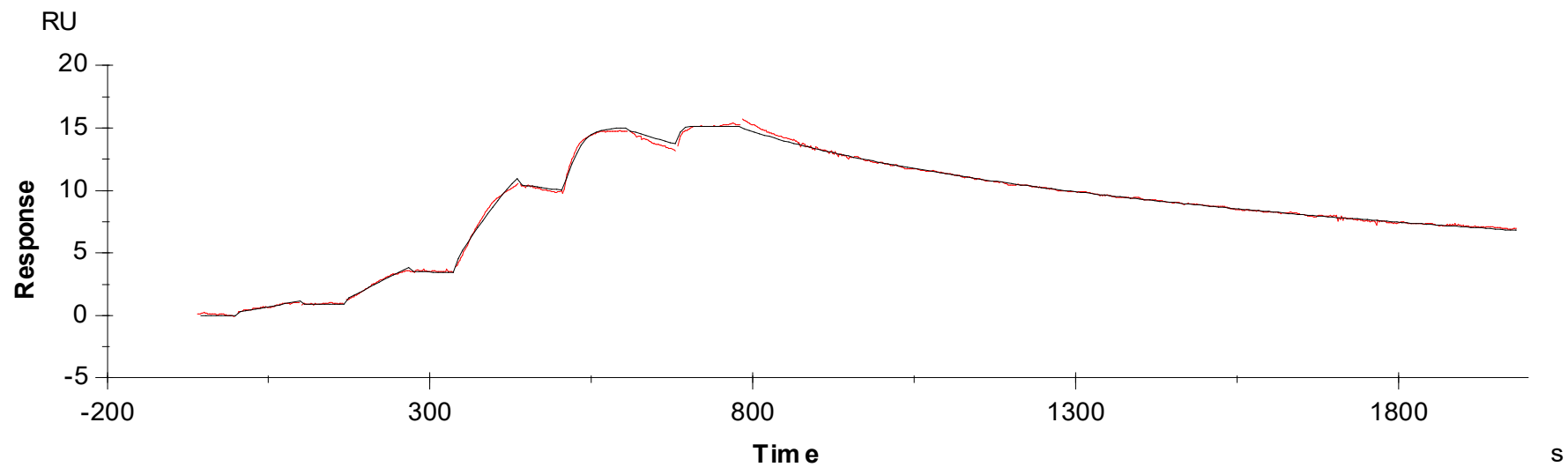


Binding Kinetics of GSK1838705 on ALK[C1156Y]



Data generated and provided by Carna Biosciences, Inc.

Binding Kinetics of TAE684 on ALK[C1156Y]



Data generated and provided by Carna Biosciences, Inc.