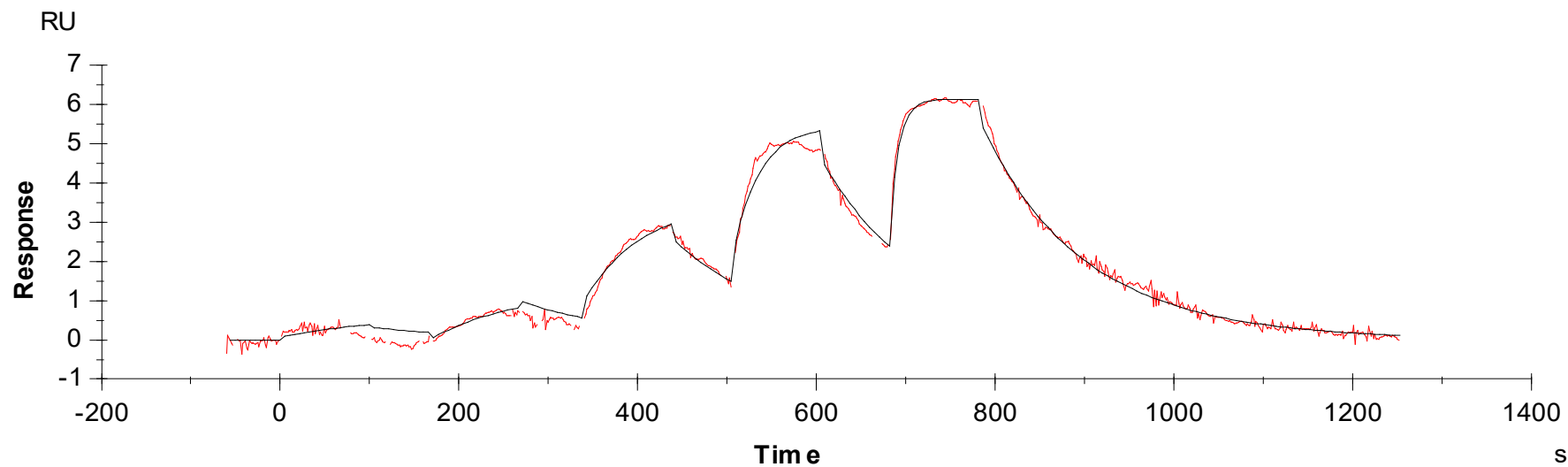


Binding Kinetics of Crizotinib on ALK[F1174L]



Data generated and provided by Carna Biosciences, Inc.