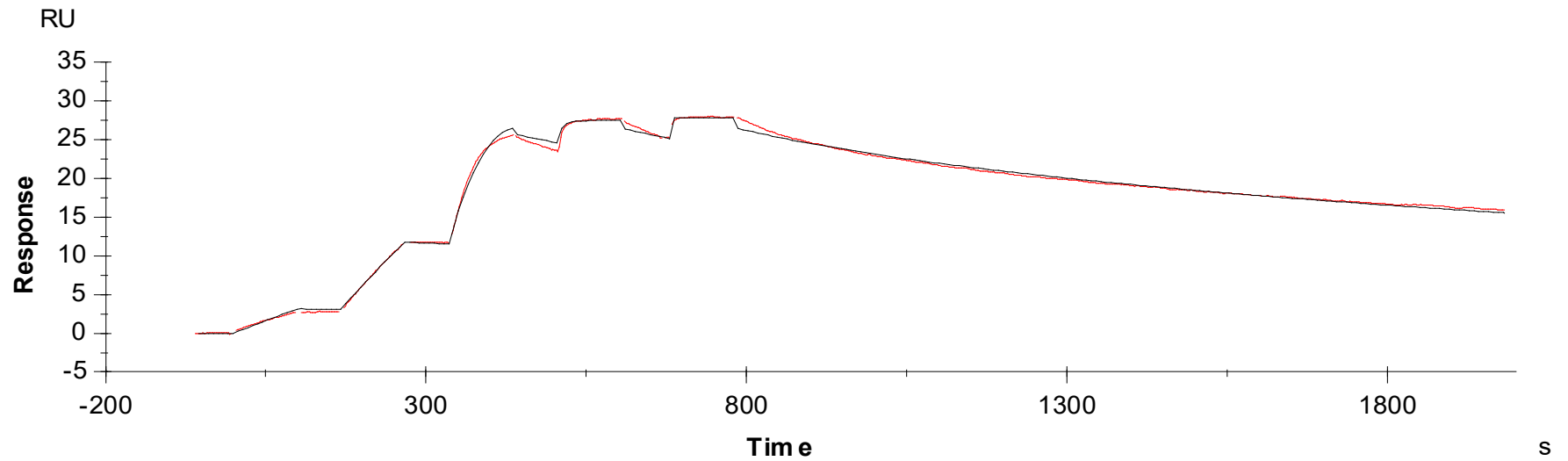


Binding Kinetics of Gilteritinib on ALK [L1198F]



Data generated and provided by Carna Biosciences, Inc.