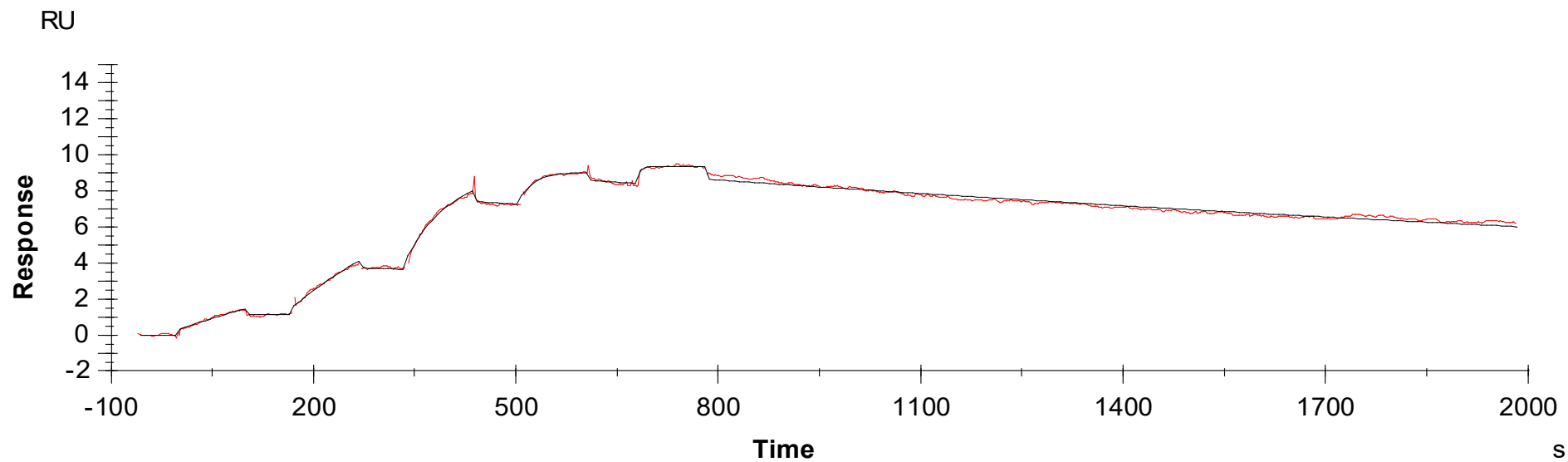


Binding Kinetics of Brigatinib on EGFR[d746-750/T790M/C797S/L858R]



Data generated and provided by Carna Biosciences, Inc.