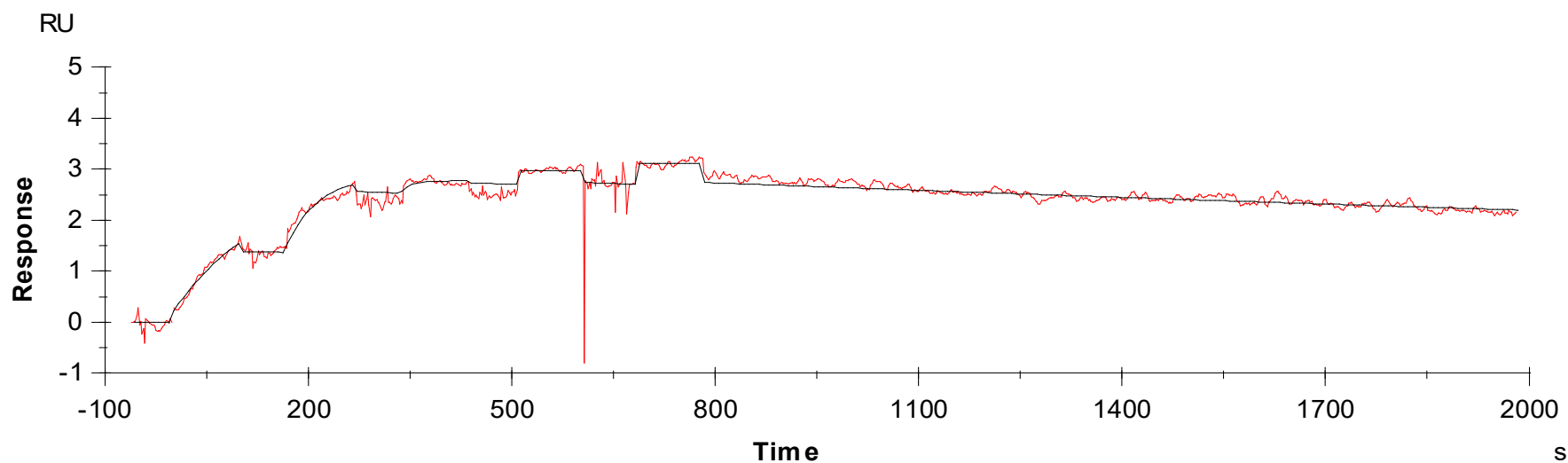




Binding Kinetics of Staurosporine on EGFR[T790M/C797S/L858R]



Data generated and provided by Carna Biosciences, Inc.