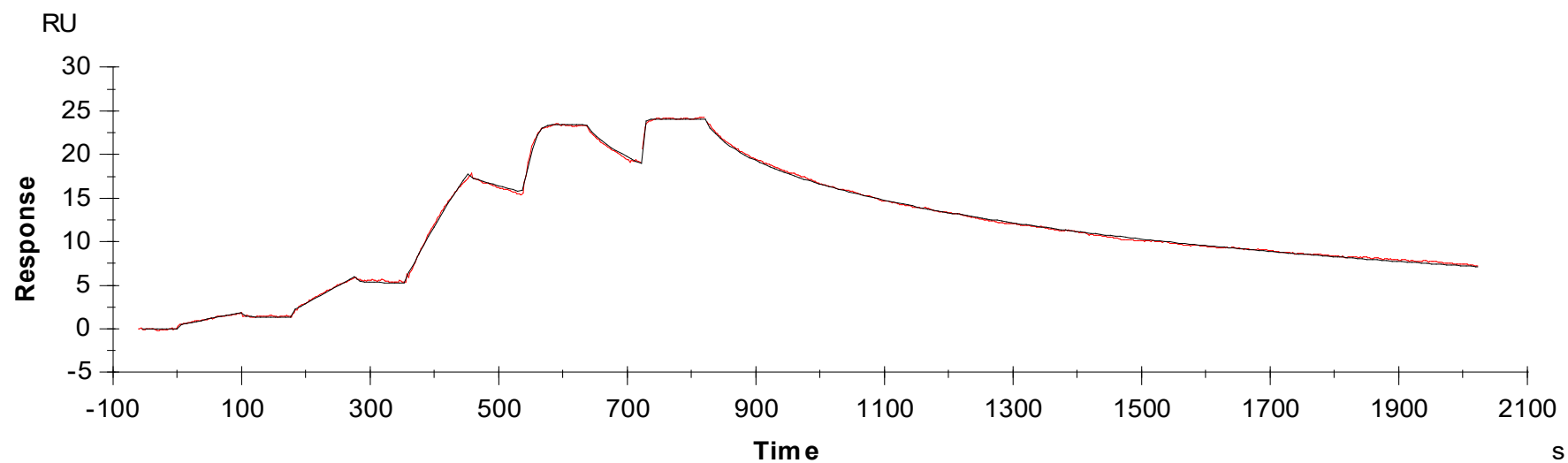




Binding Kinetics of Cediranib on FGFR2[E565A][non-activated]



Data generated and provided by Carna Biosciences, Inc.