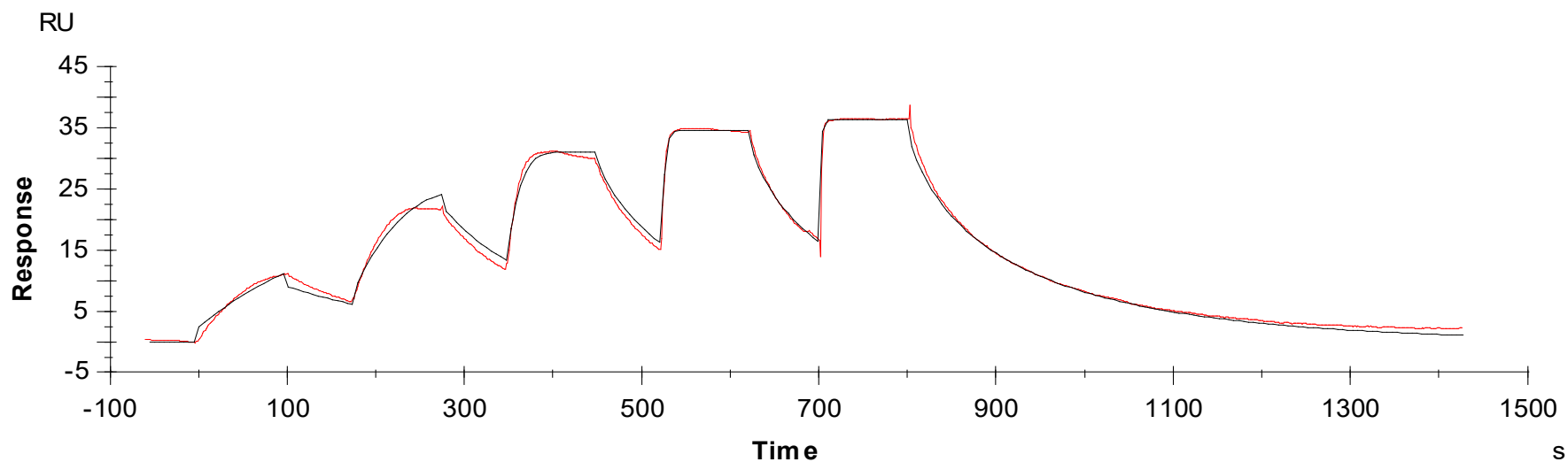




Binding Kinetics of TAE684 on FGFR2[V564F][non-activated]



Data generated and provided by Carna Biosciences, Inc.