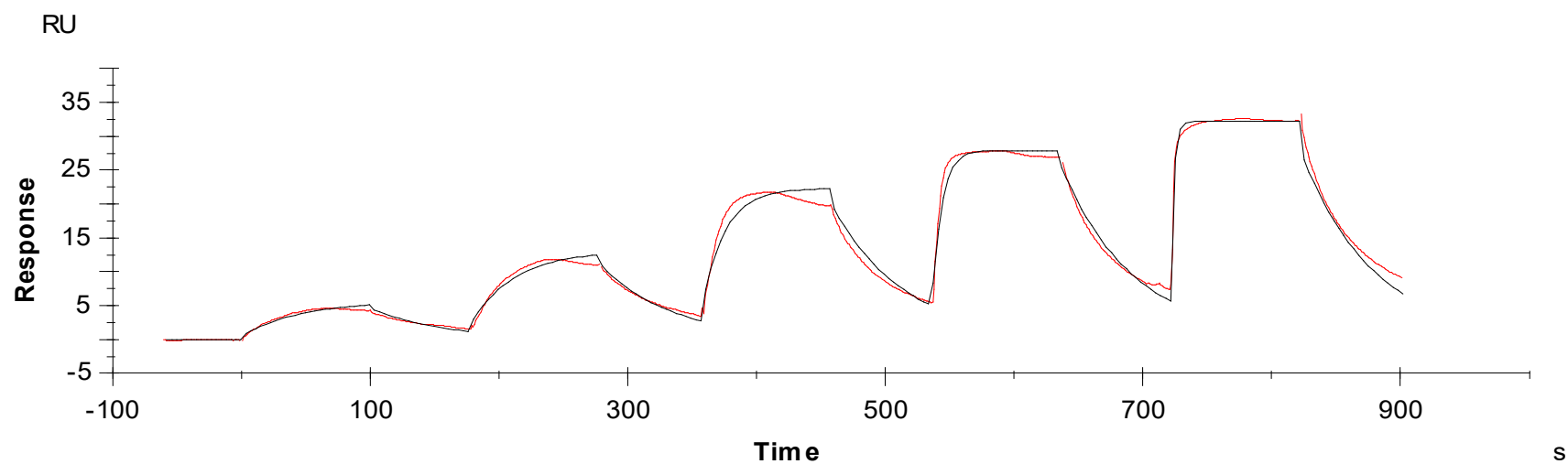




## Binding Kinetics of TAE684 on FGFR2[K659M][non-activated]



Data generated and provided by Carna Biosciences, Inc.