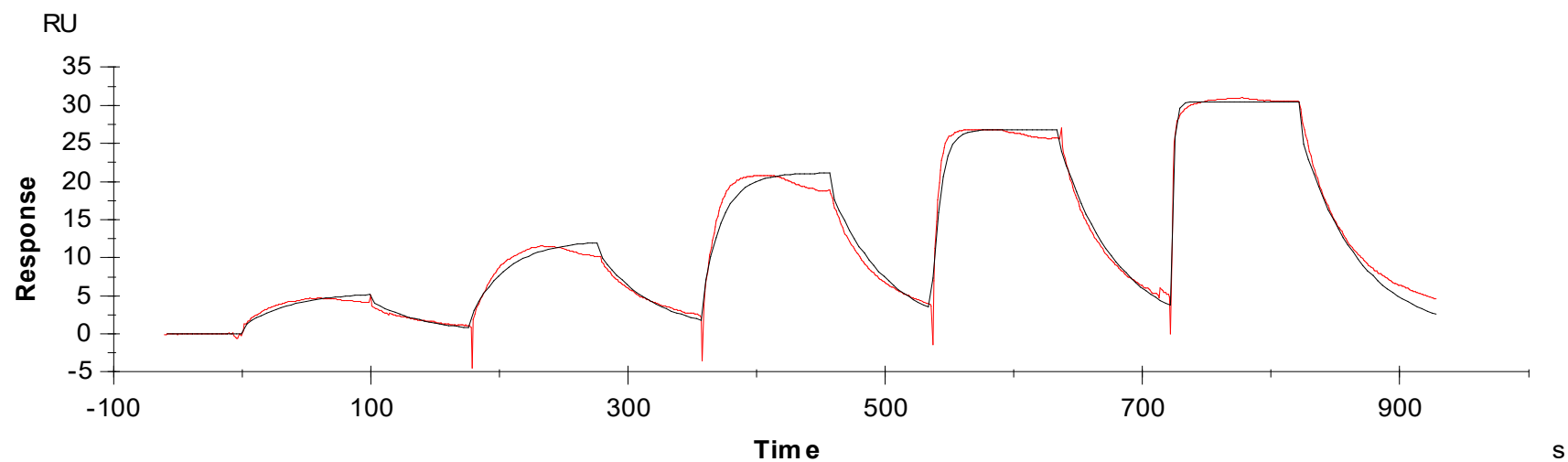




Binding Kinetics of TAE684 on FGFR2[K659M]



Data generated and provided by Carna Biosciences, Inc.