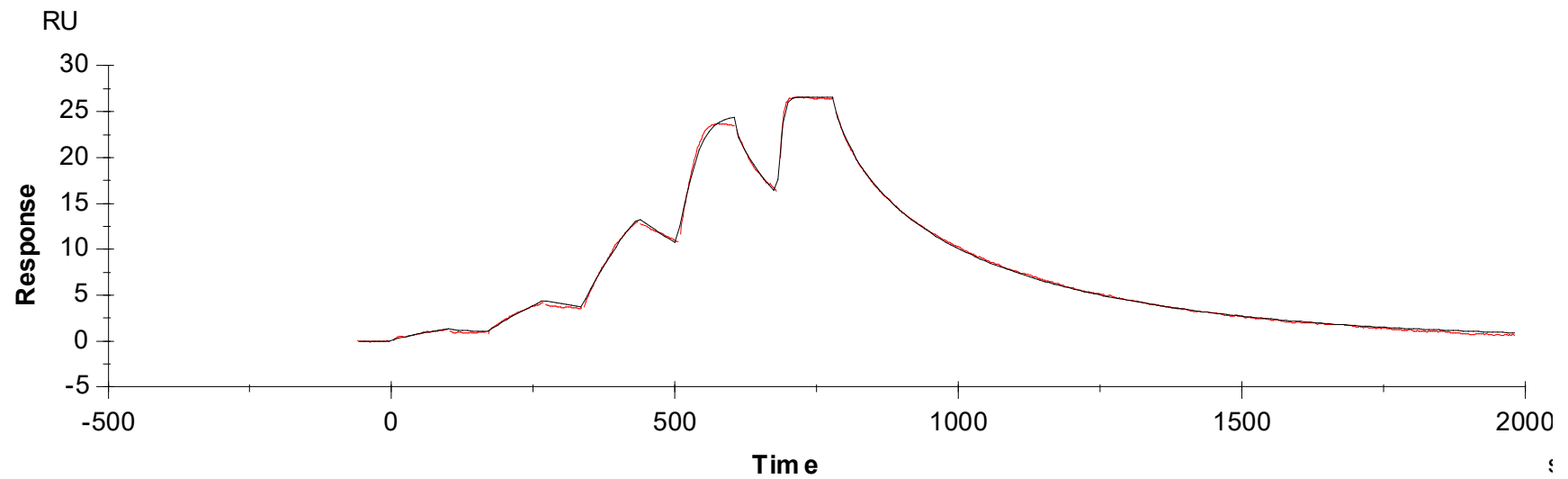


## Binding Kinetics of Cediranib on RET[Y791F] (non-activated)



Data generated and provided by Carna Biosciences, Inc.