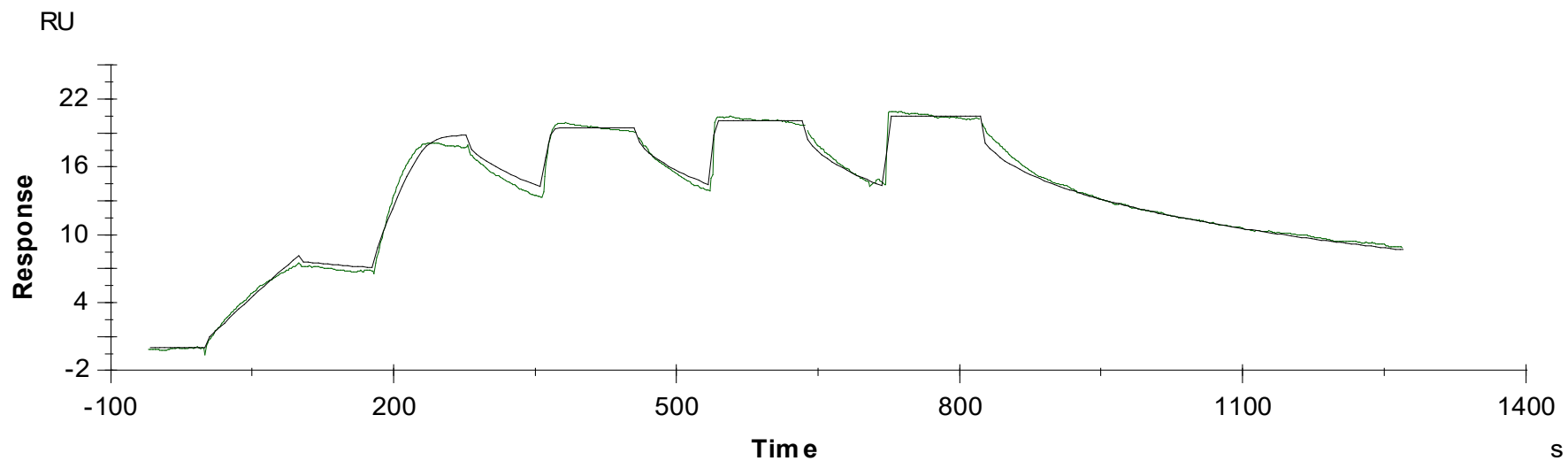


Binding Kinetics of Cediranib on FGFR2[N549K](non-activated)



Data generated and provided by Carna Biosciences, Inc.