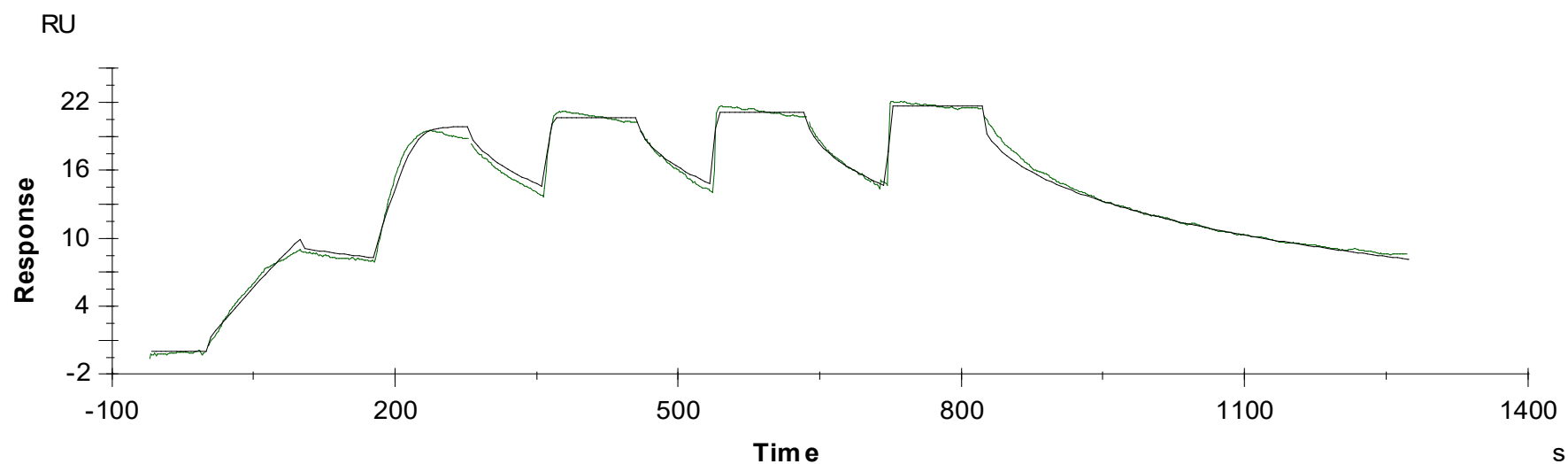


Binding Kinetics of Cediranib on FGFR2[N549K]



Data generated and provided by Carna Biosciences, Inc.