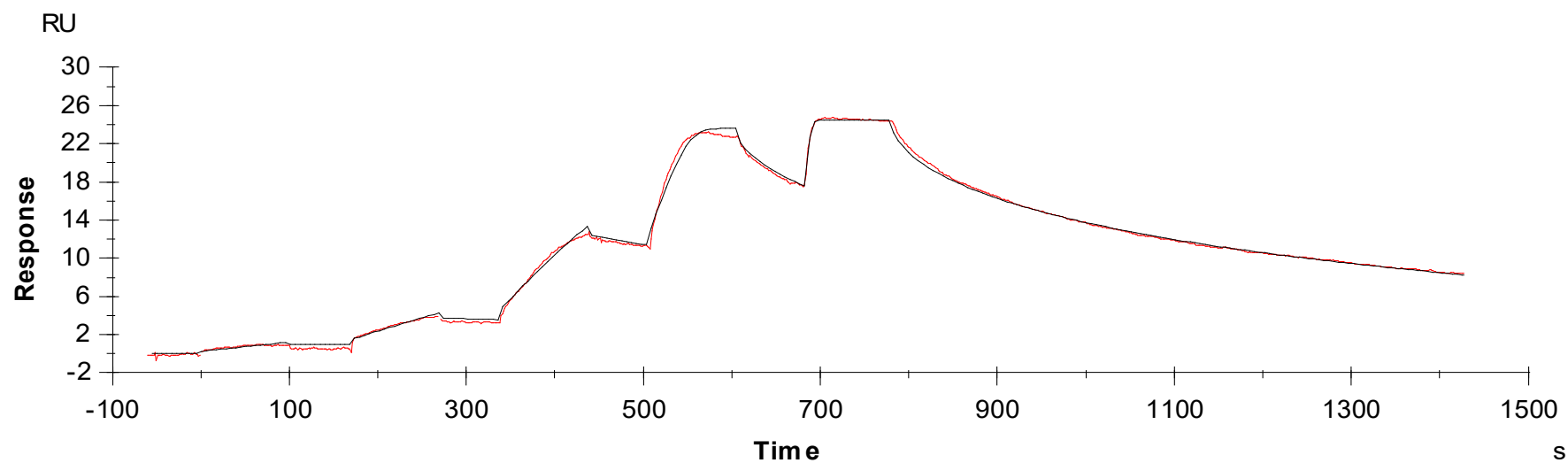




Binding Kinetics of Cediranib on FGFR2[N549D]



Data generated and provided by Carna Biosciences, Inc.