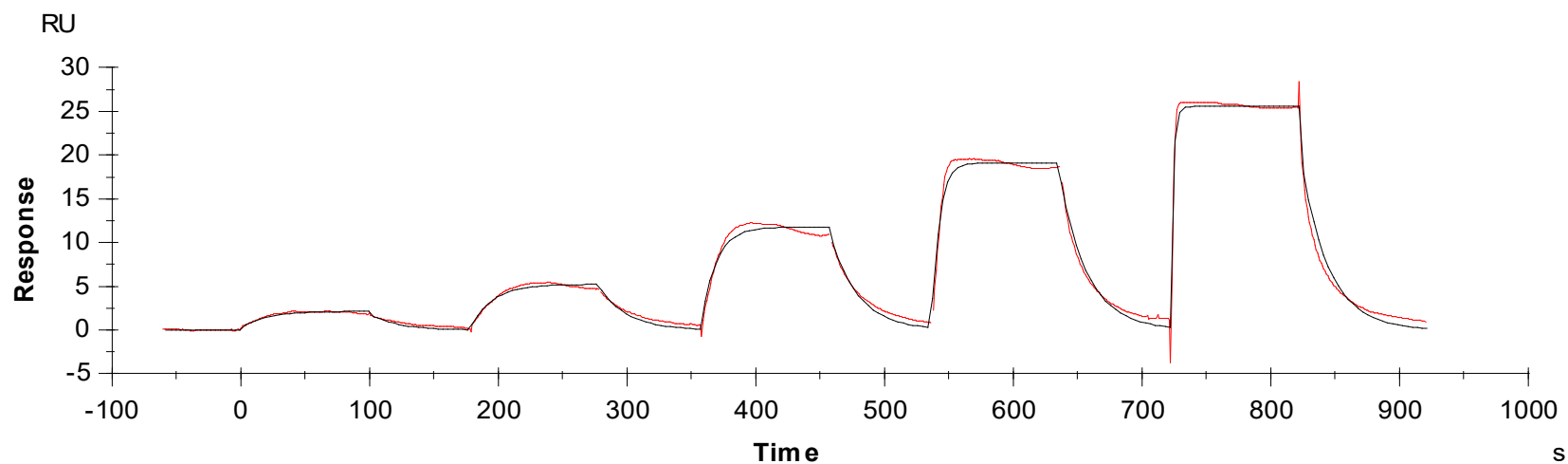




Binding Kinetics of Axitinib on FGFR2[L617V][non-activated]



Data generated and provided by Carna Biosciences, Inc.