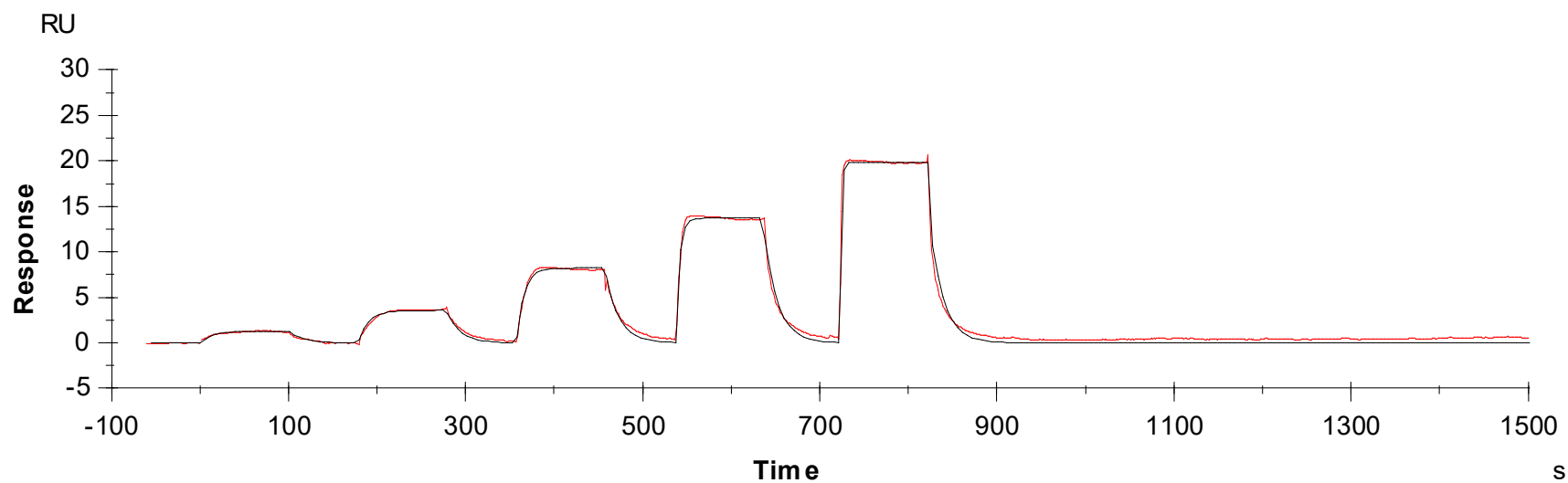


Binding Kinetics of Axitinib on FGFR2[V564I]



Data generated and provided by Carna Biosciences, Inc.