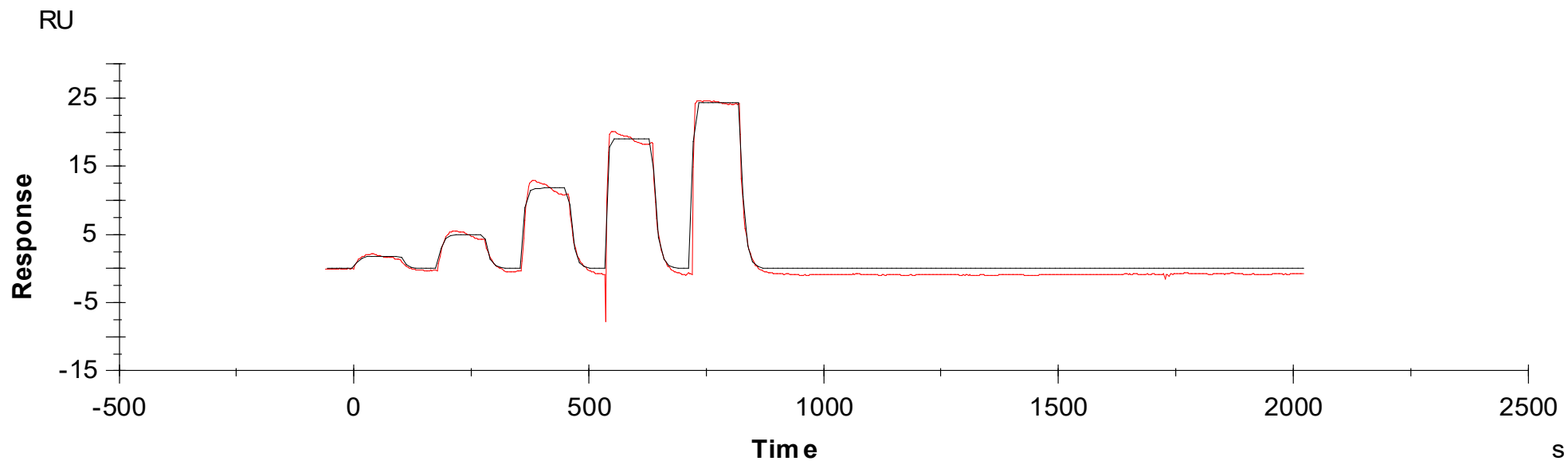


## Binding Kinetics of dovitinib on FGFR2[C491A/V564I]



Data generated and provided by Carna Biosciences, Inc.