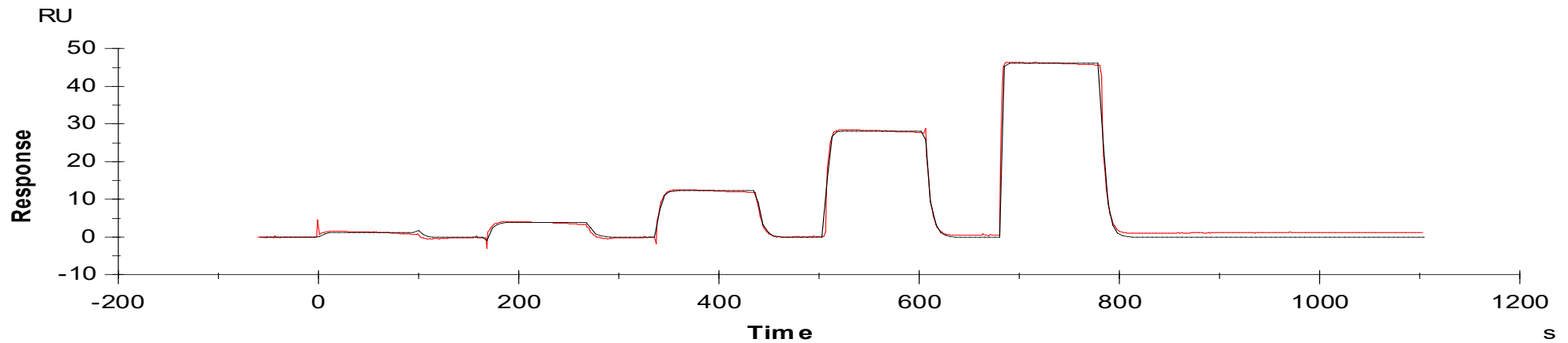


## Binding Kinetics of IGF1R[non-activated] on Brigatinib



Data generated and provided by Carna Biosciences, Inc.