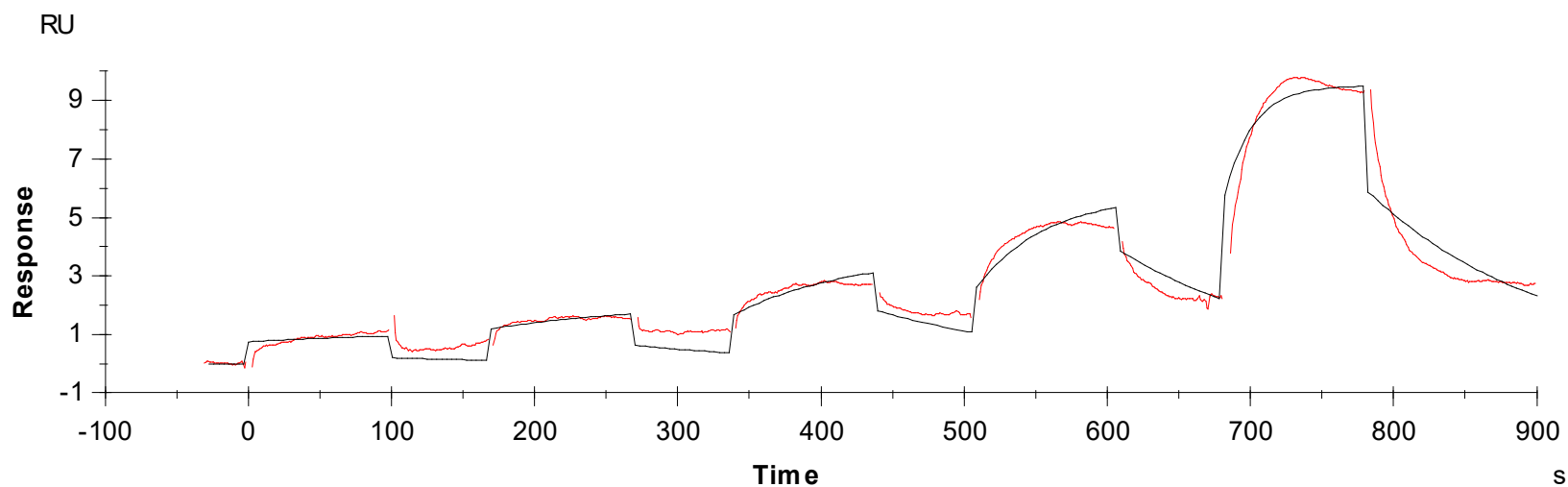


Binding Kinetics of ponatinib on ROR2



Data generated and provided by Carna Biosciences, Inc.