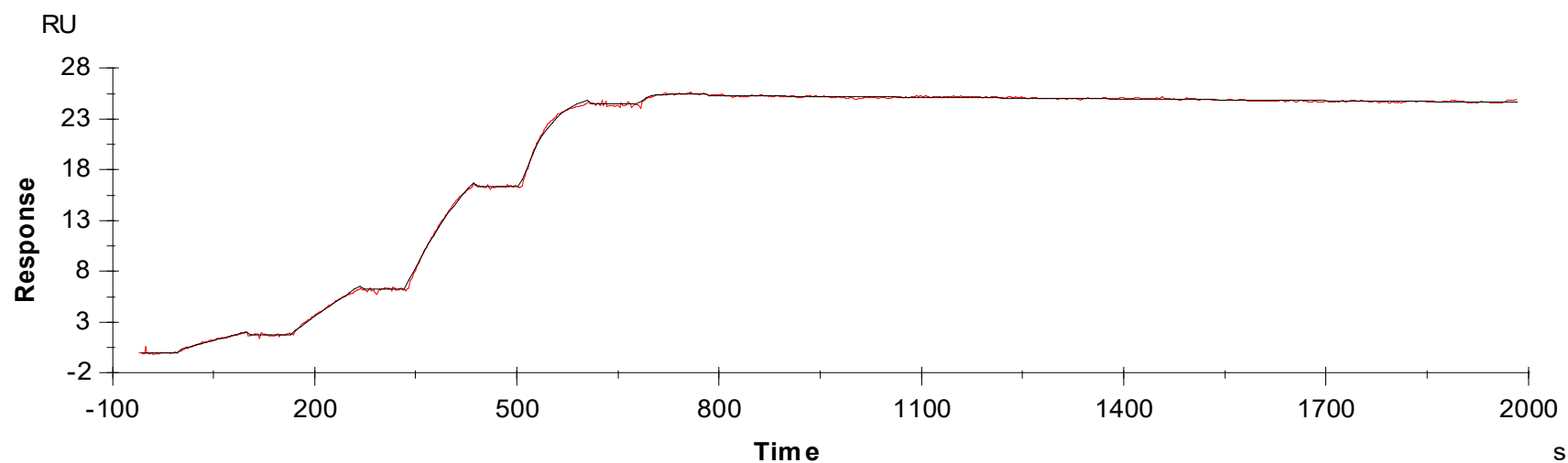




Binding Kinetics of Ponatinib on RET[non-activated]



Data generated and provided by Carna Biosciences, Inc.