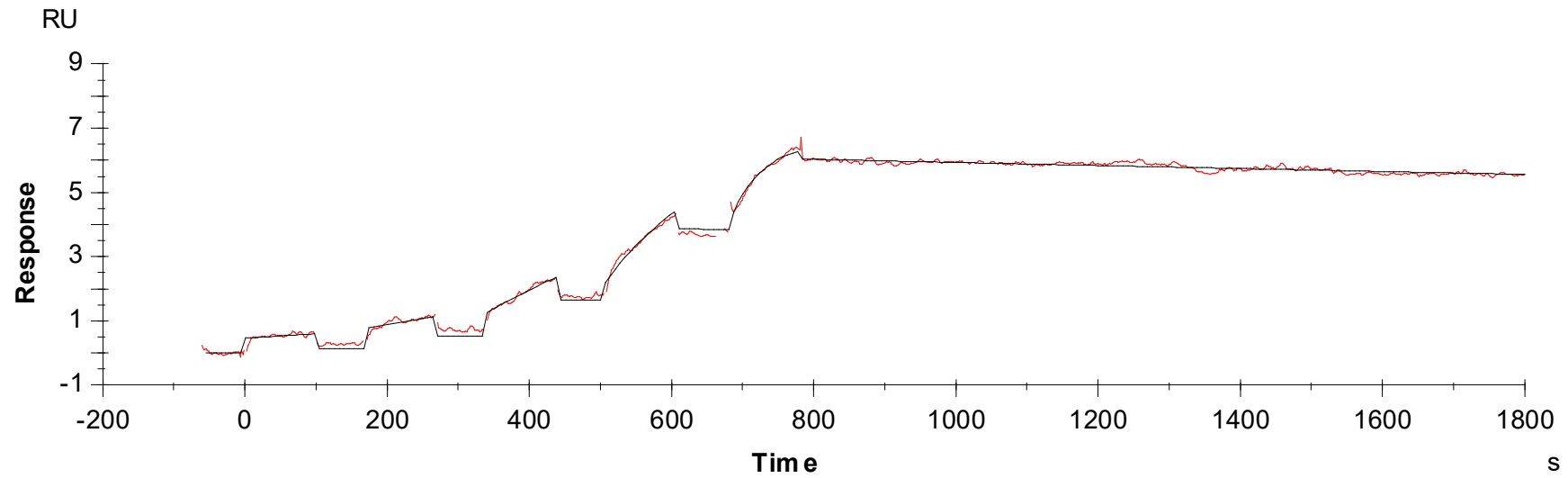




Binding Kinetics of Quizartinib on RET[S891A](non-activated)



Data generated and provided by Carna Biosciences, Inc.