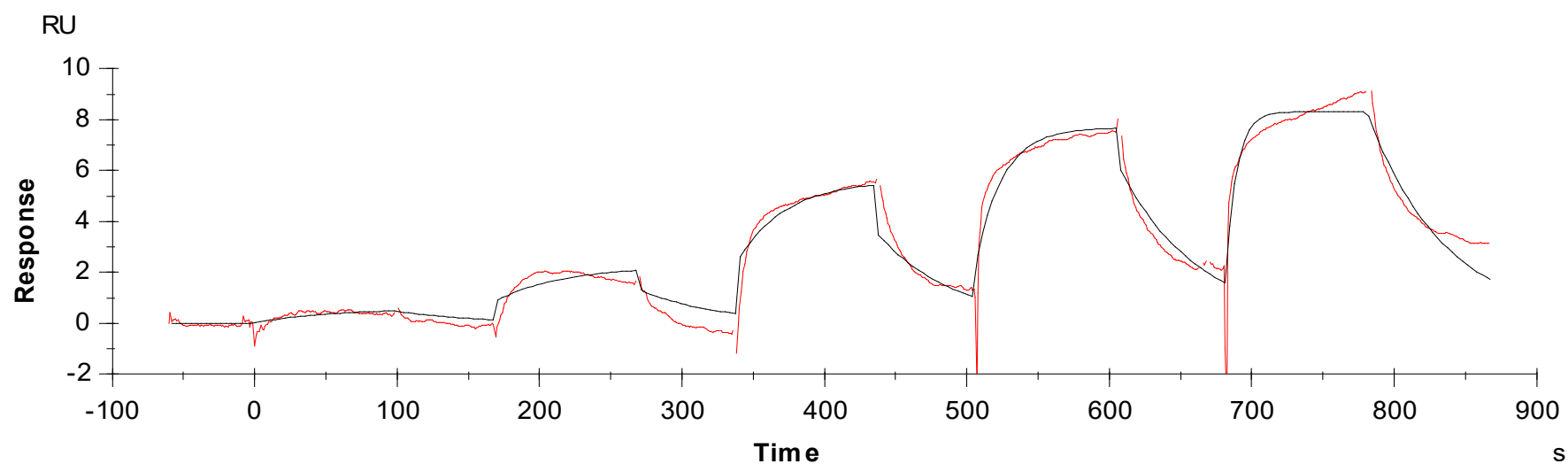




Binding Kinetics of bosutinib on ABL(ABL1)[T315I]



Data generated and provided by Carna Biosciences, Inc.