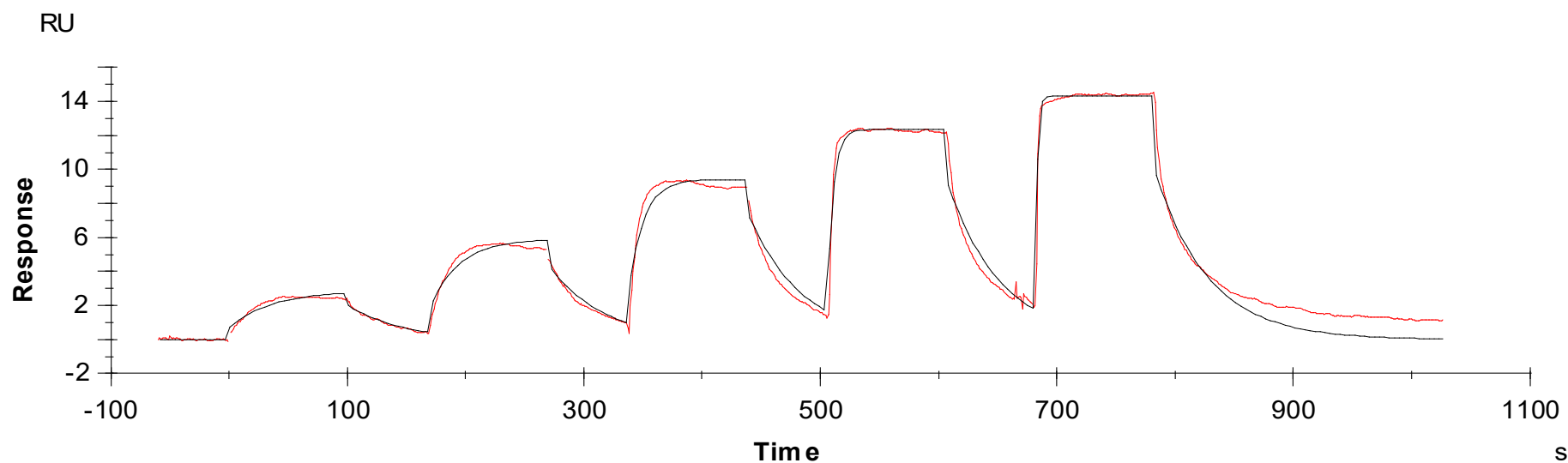




Binding Kinetics of Bosutinib on PKC θ (PRKCQ)



Data generated and provided by Carna Biosciences, Inc.