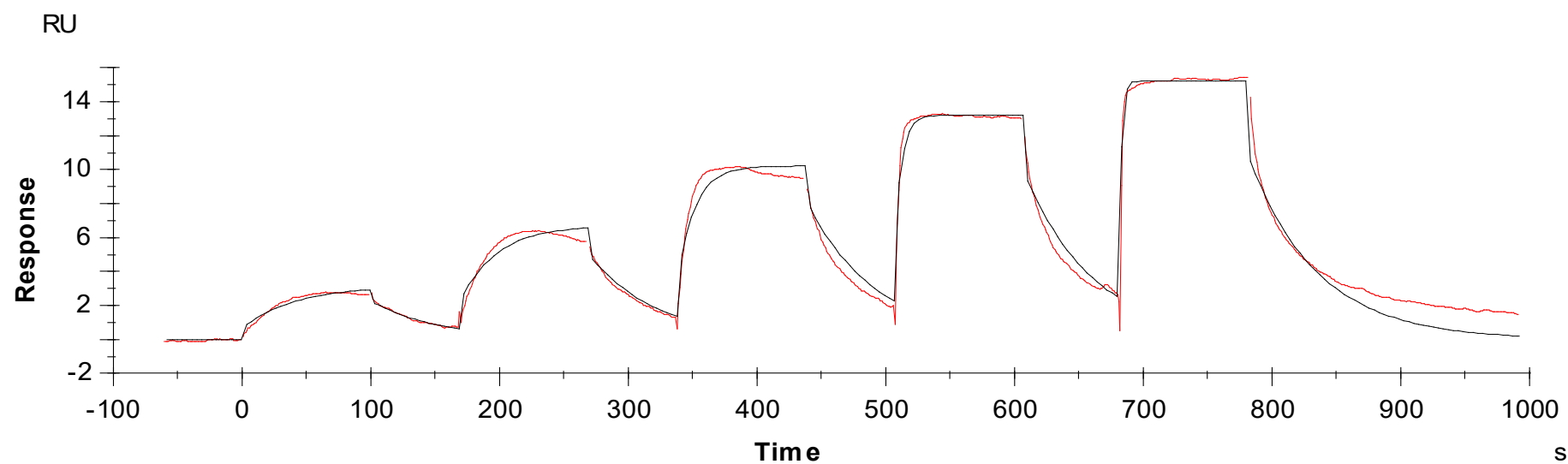




Binding Kinetics of Bosutinib on PKC θ (PRKCQ)[non-activated]



Data generated and provided by Carna Biosciences, Inc.