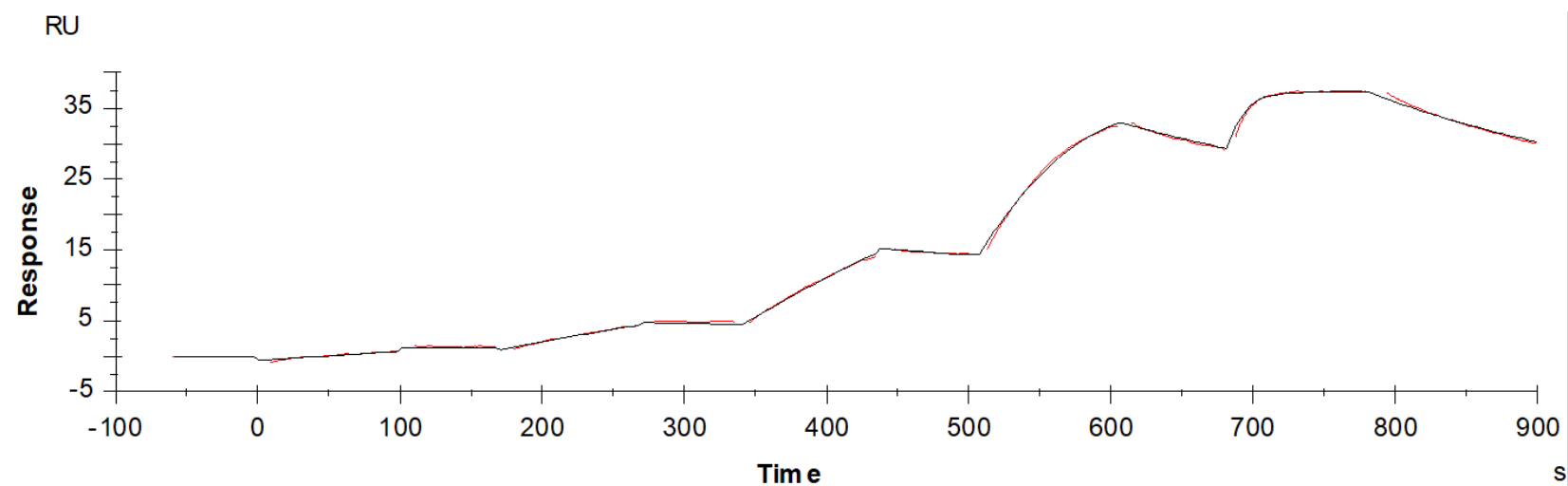




Binding Kinetics of Bosutinib on TNIK [non-activated]



Data generated and provided by Carna Biosciences, Inc.