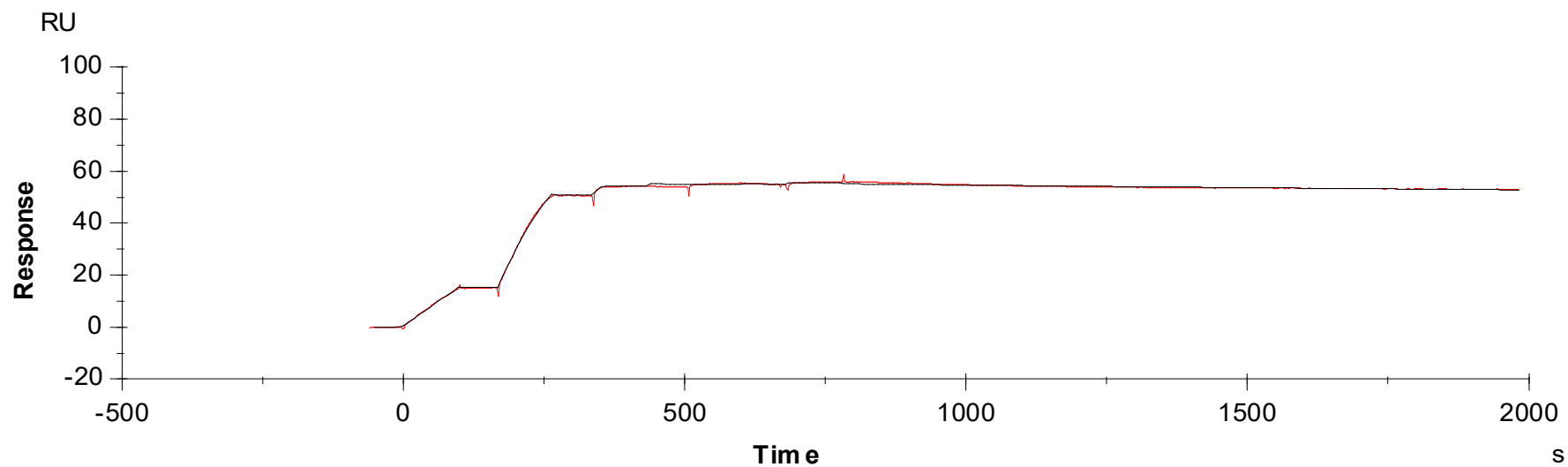


Binding Kinetics of Dasatinib on EphA2 [non-activated]



Data generated and provided by Carna Biosciences, Inc.