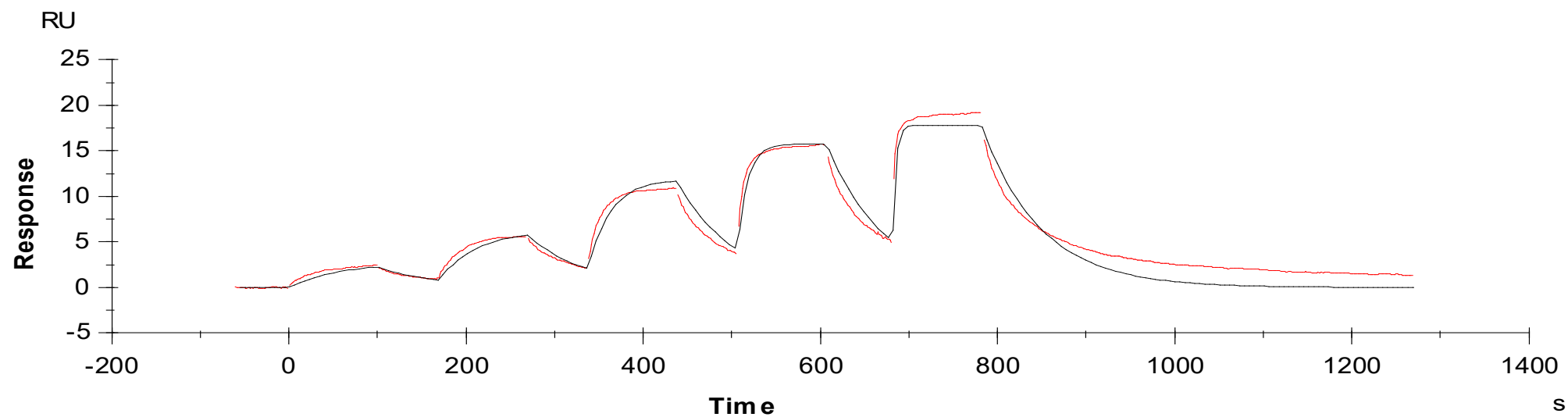


Binding Kinetics of Dasatinib on LIMK2



Data generated and provided by Carna Biosciences, Inc.