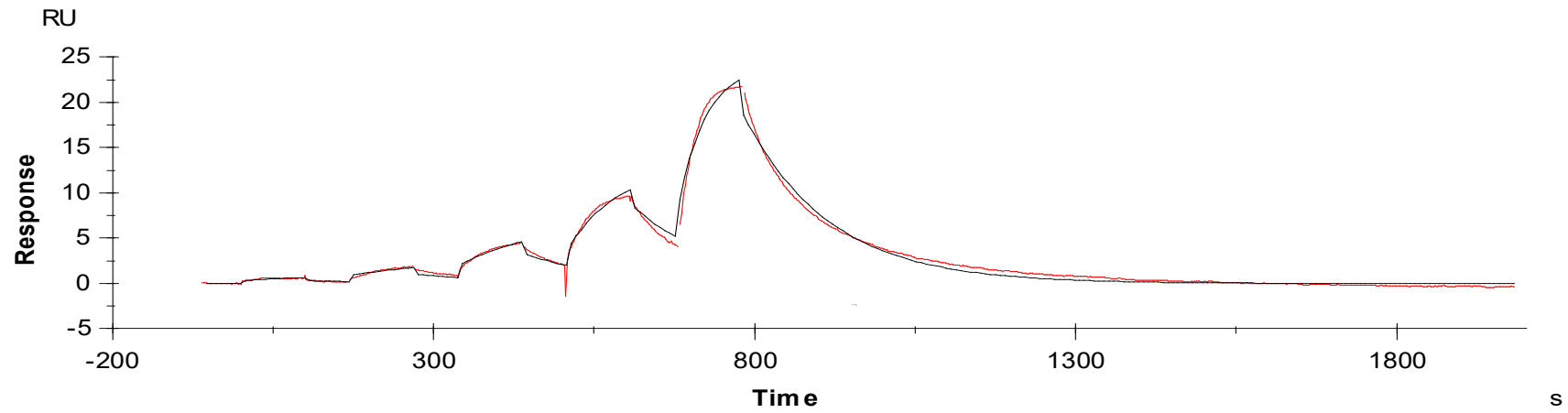


Binding Kinetics of Midostaurin on CaMK2B



Data generated and provided by Carna Biosciences, Inc.