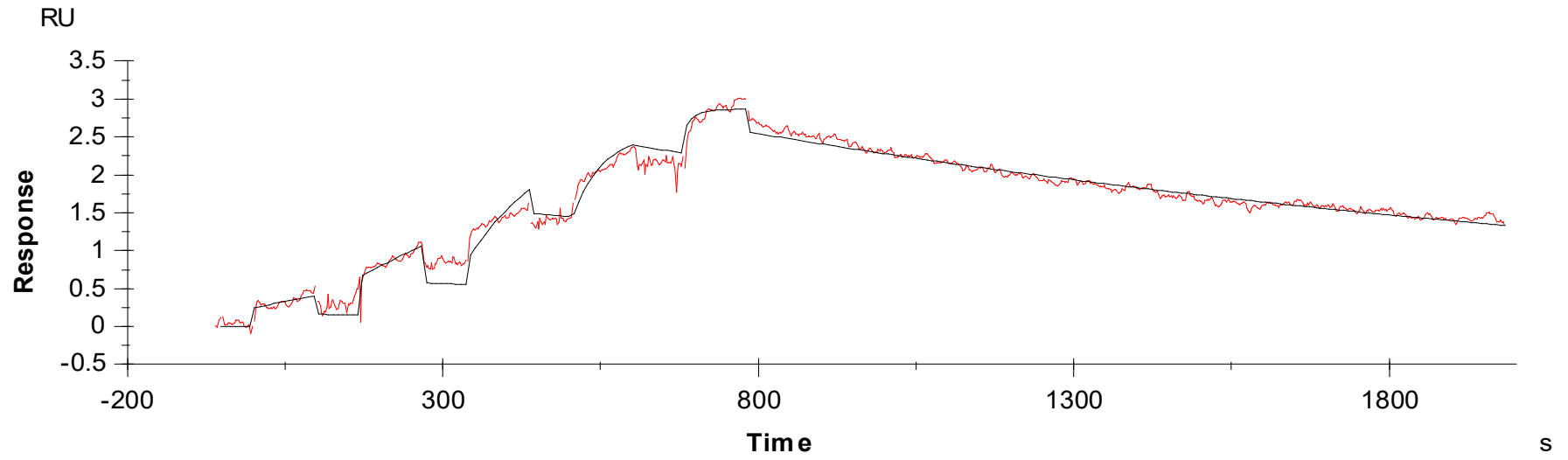
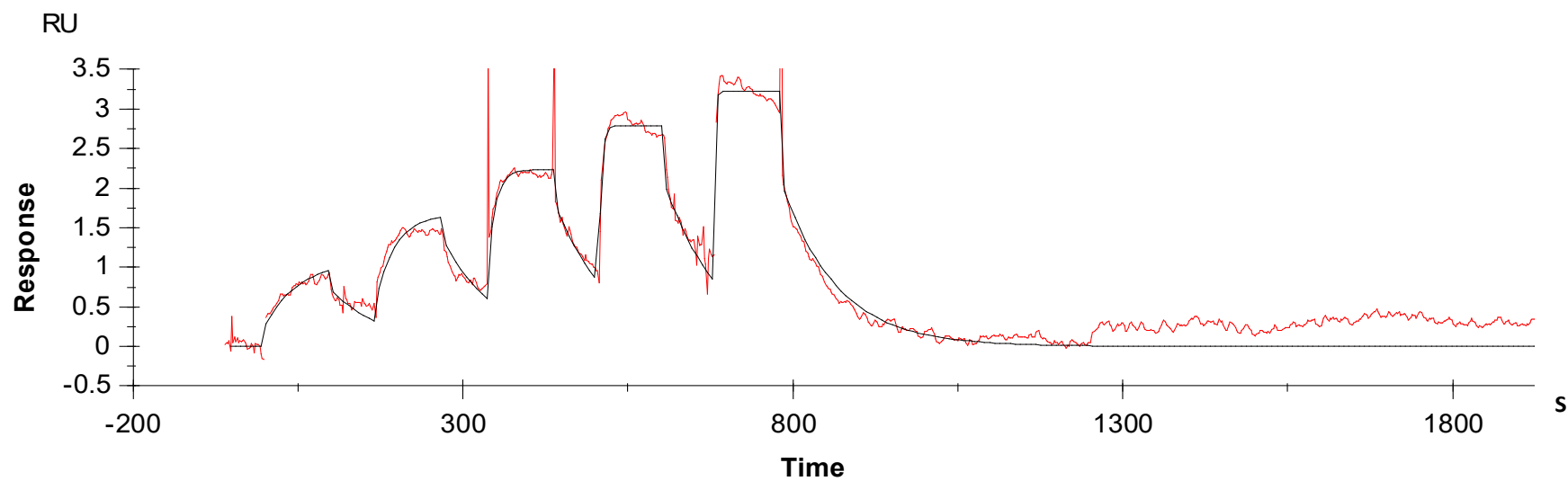


Binding Kinetics of Dasatinib on FMS



Data generated and provided by Carna Biosciences, Inc.

Binding Kinetics of Staurosporine on FMS



Data generated and provided by Carna Biosciences, Inc.