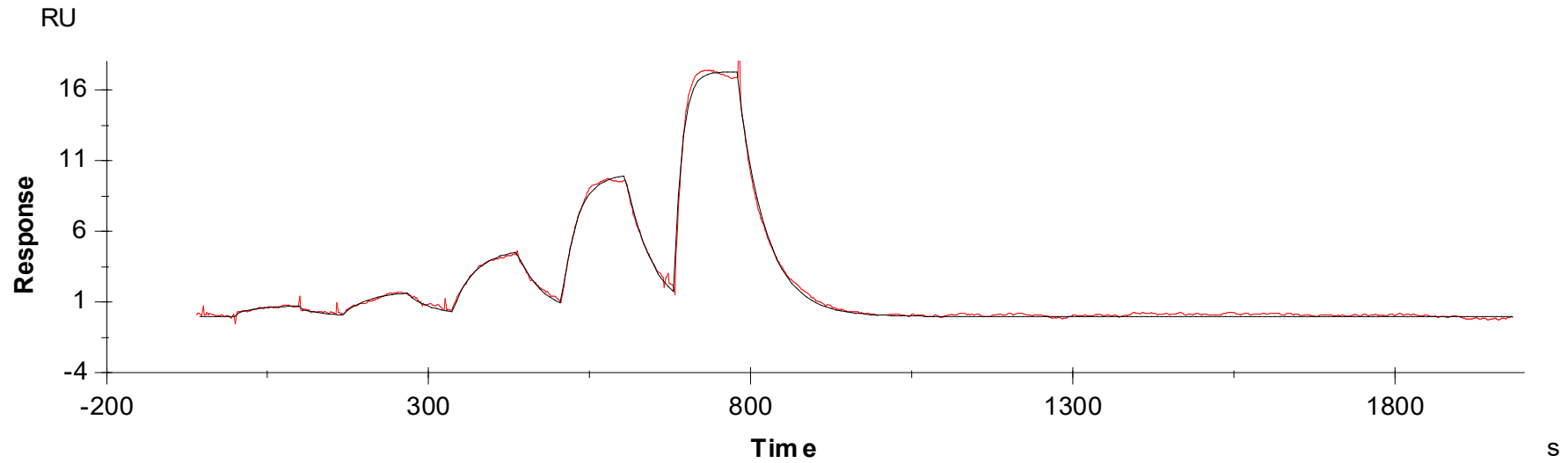


Binding Kinetics of Dovitinib on MYLK4



Data generated and provided by Carna Biosciences, Inc.