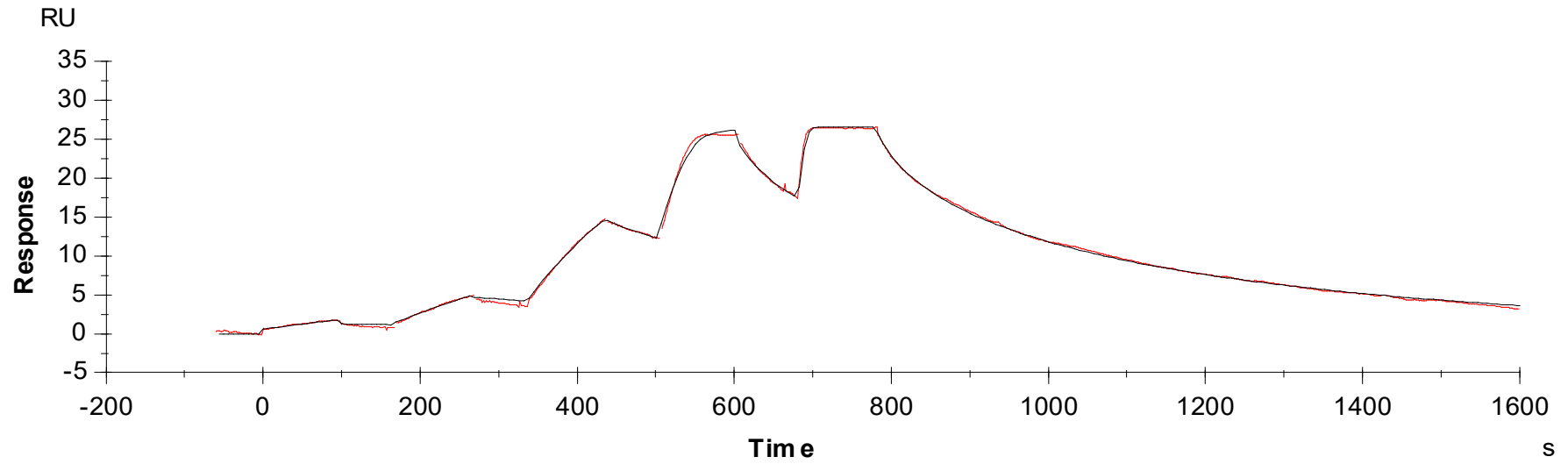


Binding Kinetics of Dovitinib on RET[S891A]



Data generated and provided by Carna Biosciences, Inc.

Binding kinetics of of staurosporine on RET [S891A]

