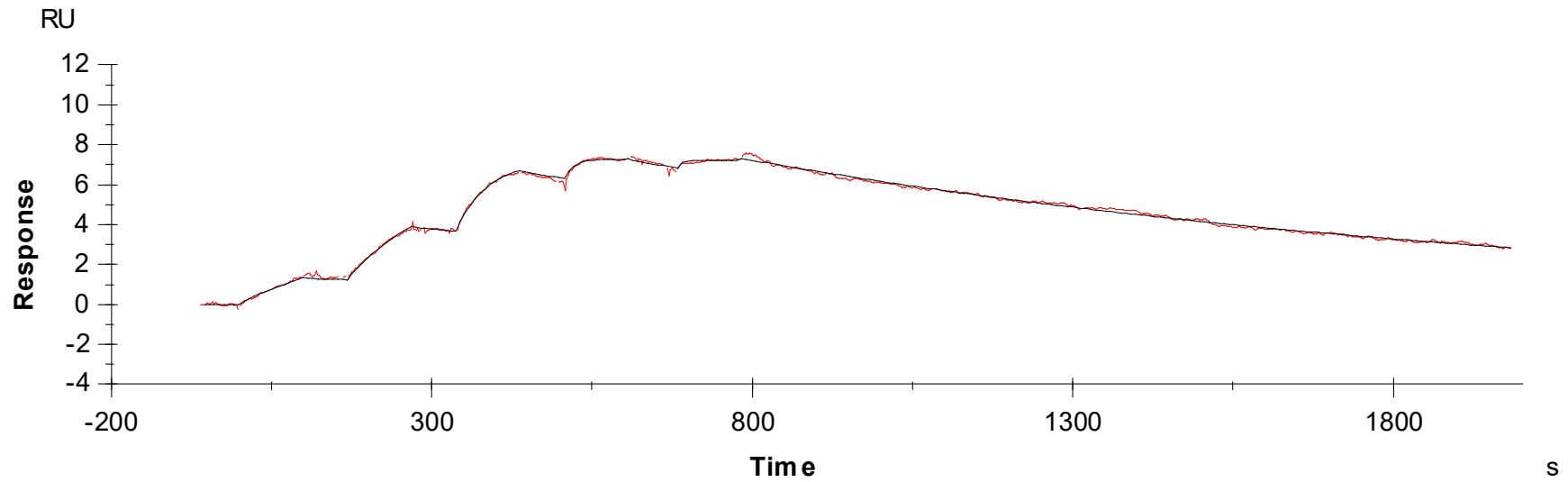


Binding Kinetics of WNK463 on WNK1



Data generated and provided by Carna Biosciences, Inc.