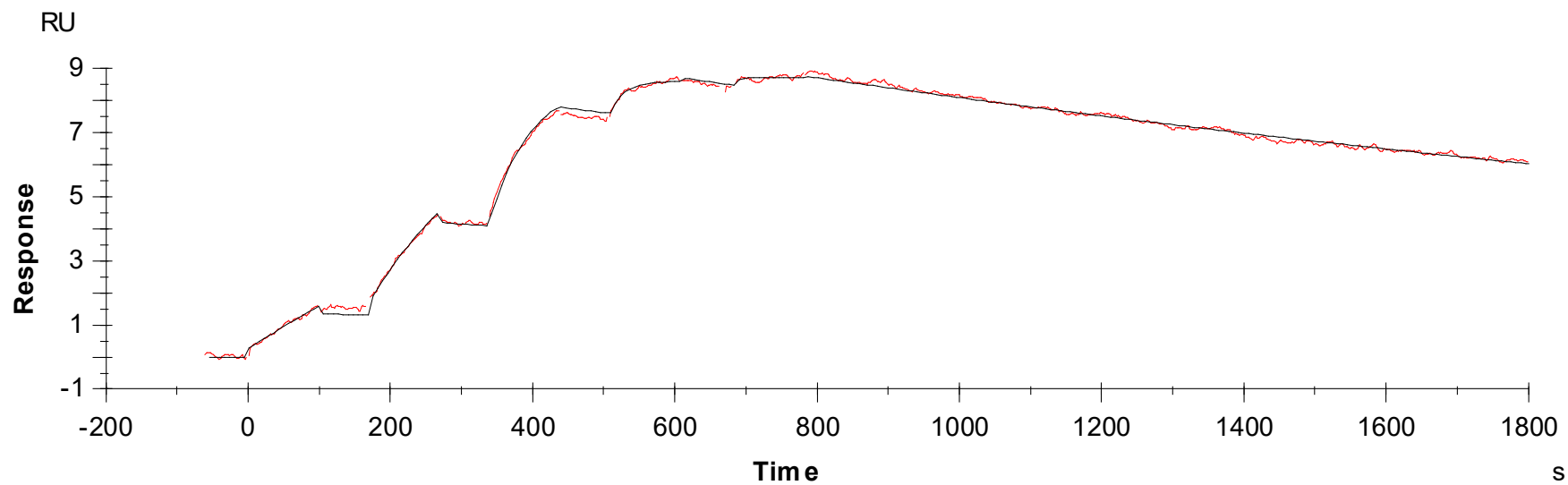


Binding Kinetics of WNK463 on WNK4



Data generated and provided by Carna Biosciences, Inc.