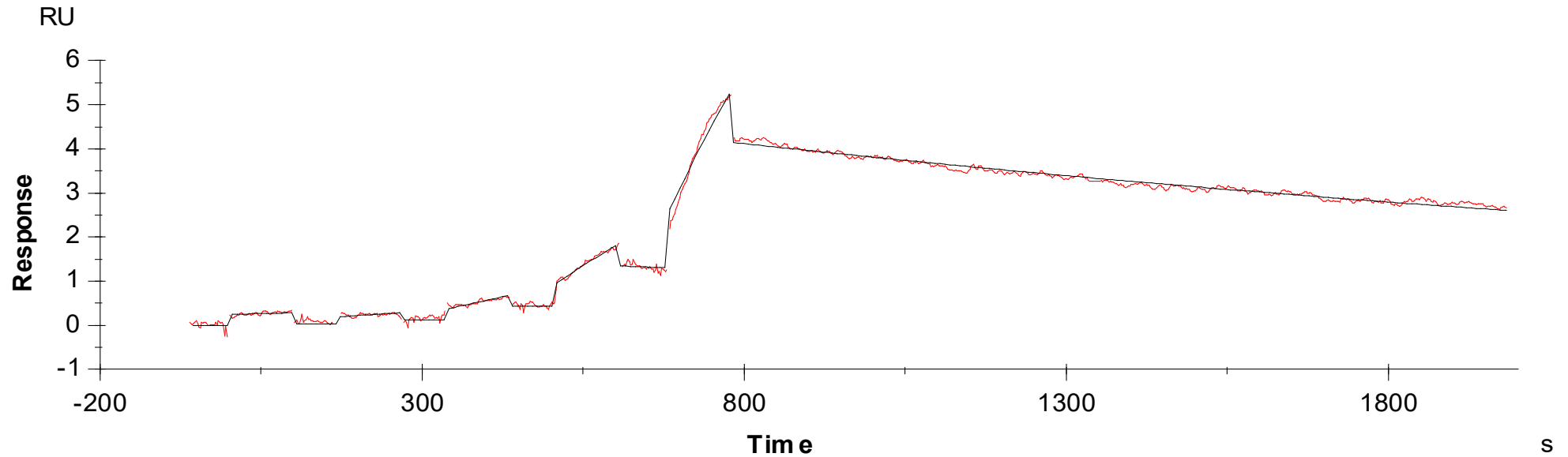


## Binding Kinetics of Compound X (Carna's proprietary) on DGK $\gamma$ (261-791aa)



8 U U [ Y b Y f U h Y X ' U b X ' d f c j ] X Y X ' V m 7 U f b U 6 ] c g V Y b W W g z - b W