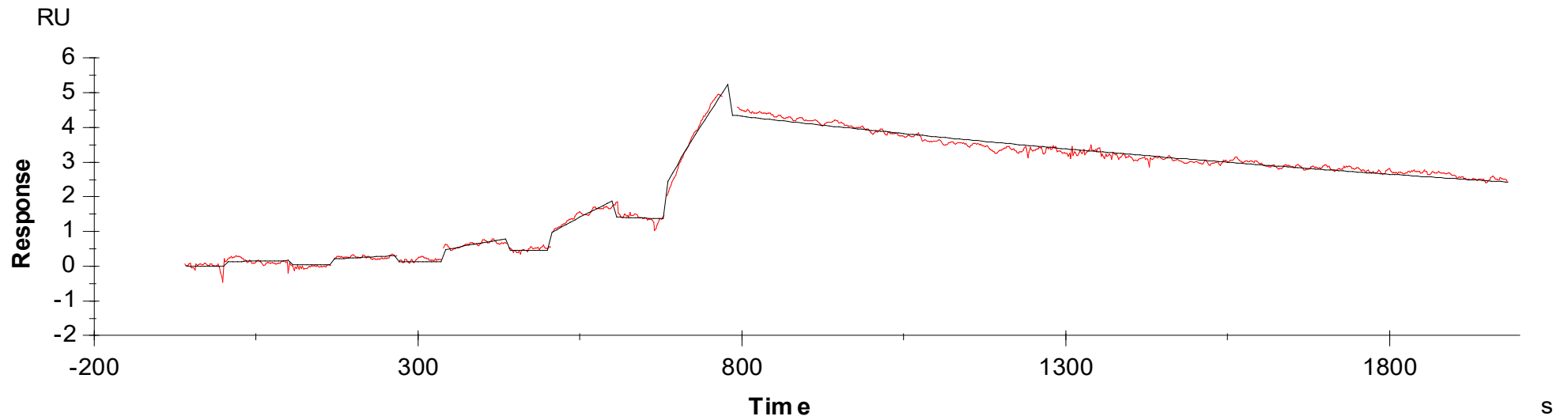


Binding Kinetics of Compound X (Carna's proprietary) on DGK β (235-803aa)



8 UU[YbYfUHX'UbX'dfcj]XYX'Vmi7 UfbU6]cgWYbWgž=bW