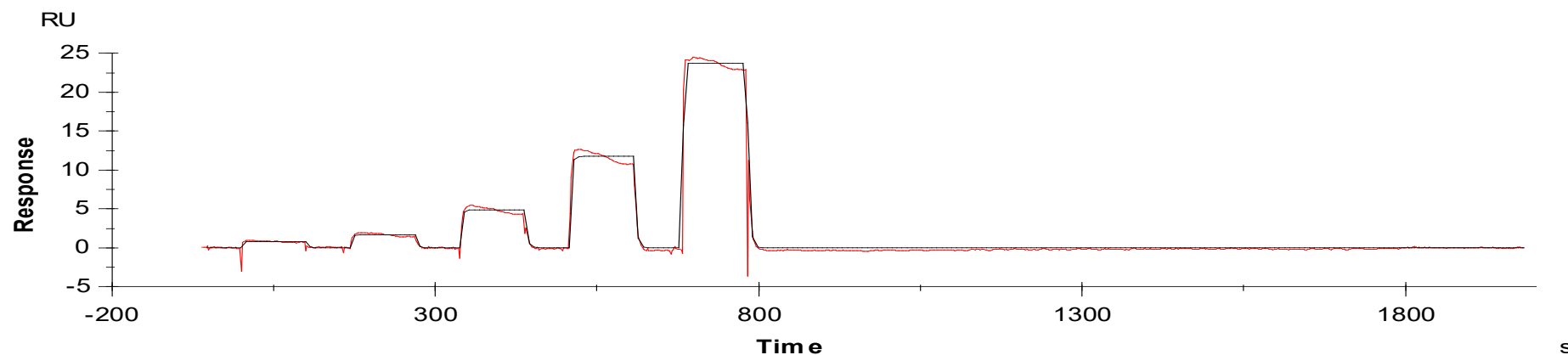


Binding Kinetics of Sunitinib on CaMK2D



Data generated and provided by Carna Biosciences, Inc.