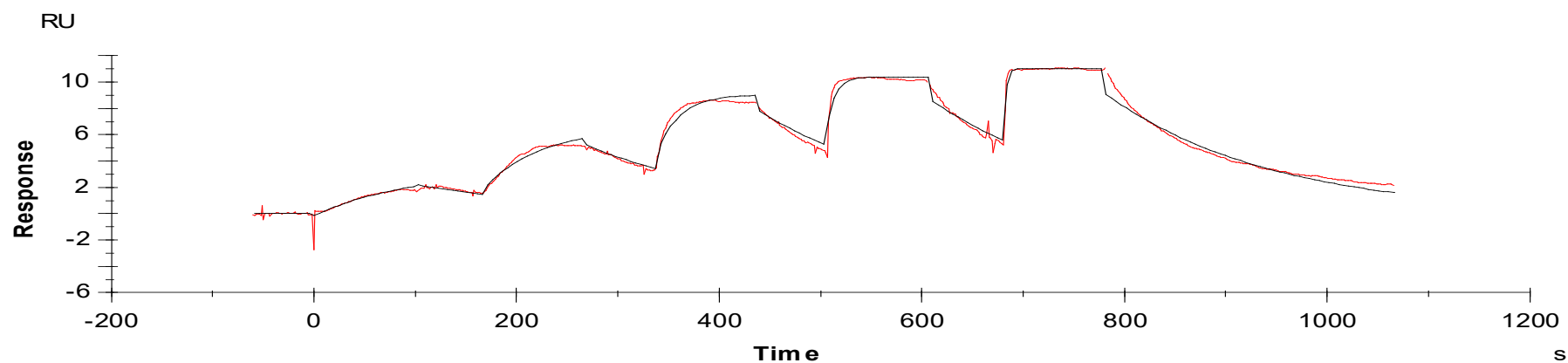


Binding Kinetics of TAE684 on AMPK α 1 β 2 γ 1



Data generated and provided by Carna Biosciences, Inc.