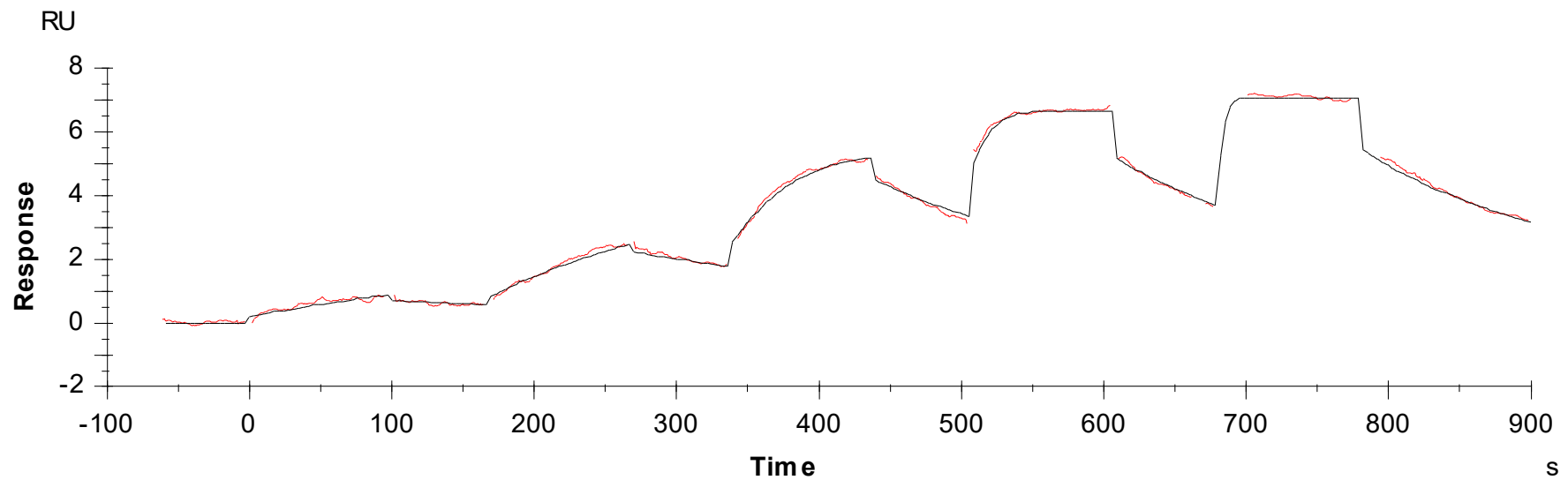


## Binding Kinetics of Pictilisib on PIK3CA[H1047L]/PIK3R1



Data generated and provided by Carna Biosciences, Inc.