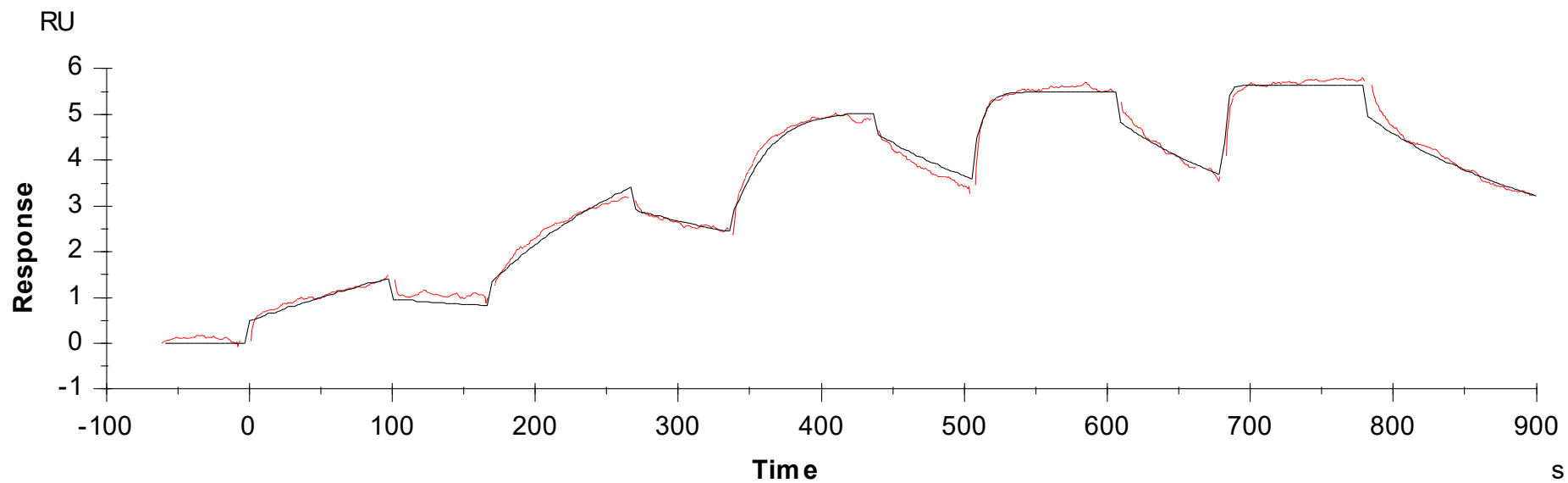


Binding Kinetics of Pictilisib on PIK3CA[C420R]/PIK3R1



Data generated and provided by Carna Biosciences, Inc.