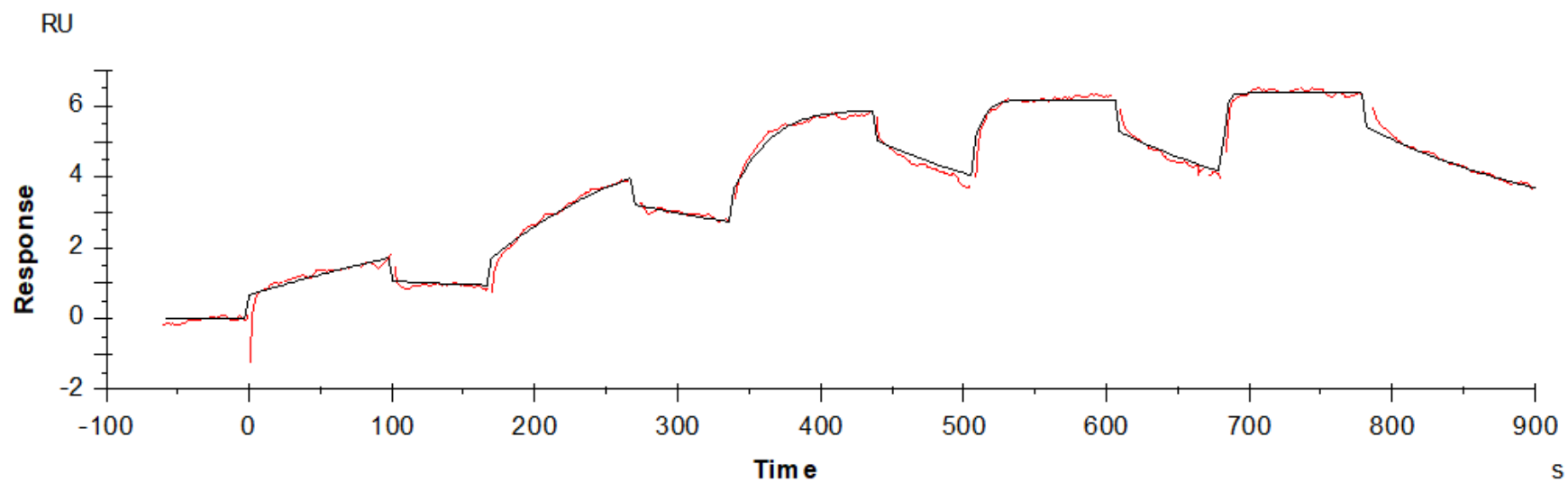


Binding Kinetics of Pictilisib on PIK3CA[N345K]/PIK3R1



Data generated and provided by Carna Biosciences, Inc.