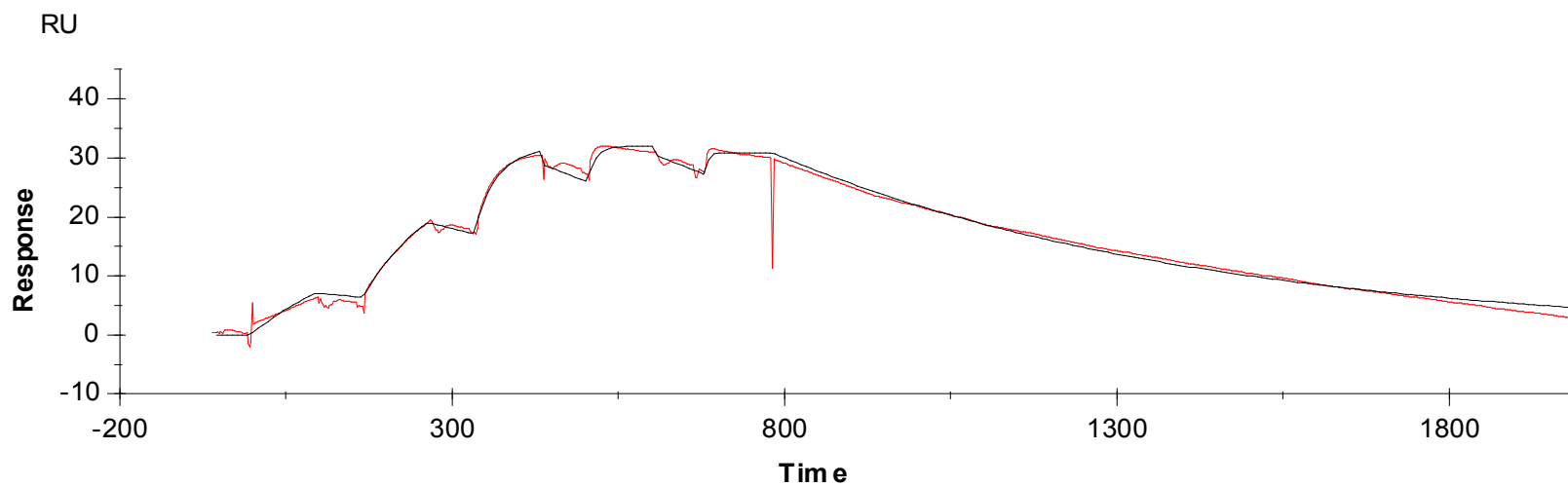


Binding Kinetics of CGI-1746 on BTK(non-activated)



Data generated and provided by Carna Biosciences