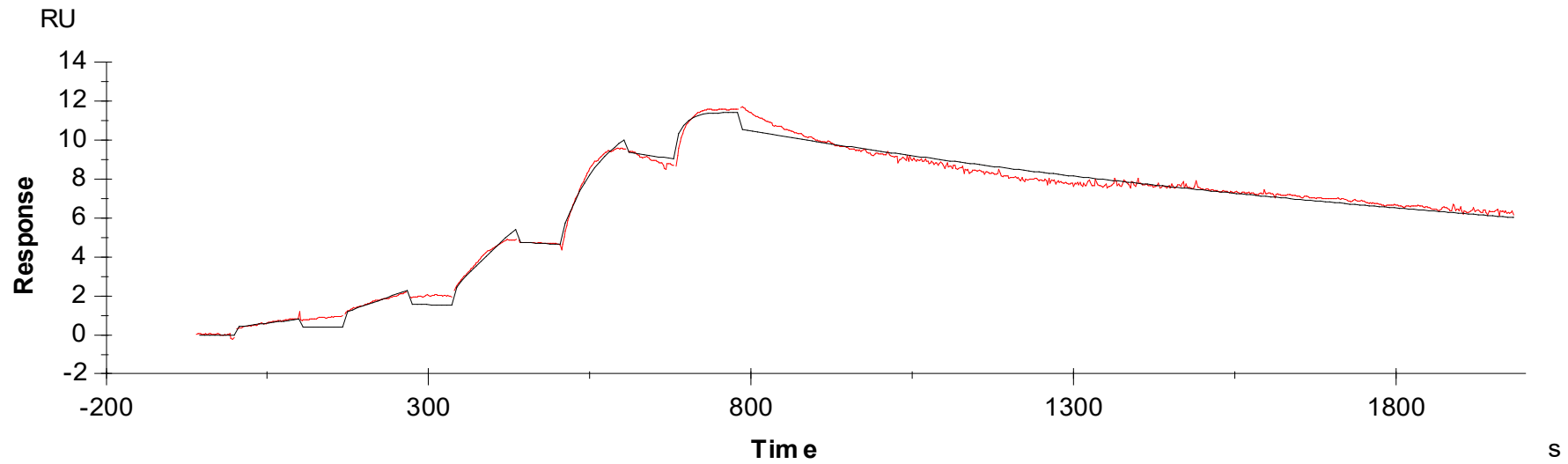


Binding Kinetics of Ceritinib on ALK [I1171S]



Data generated and provided by Carna Biosciences, Inc.