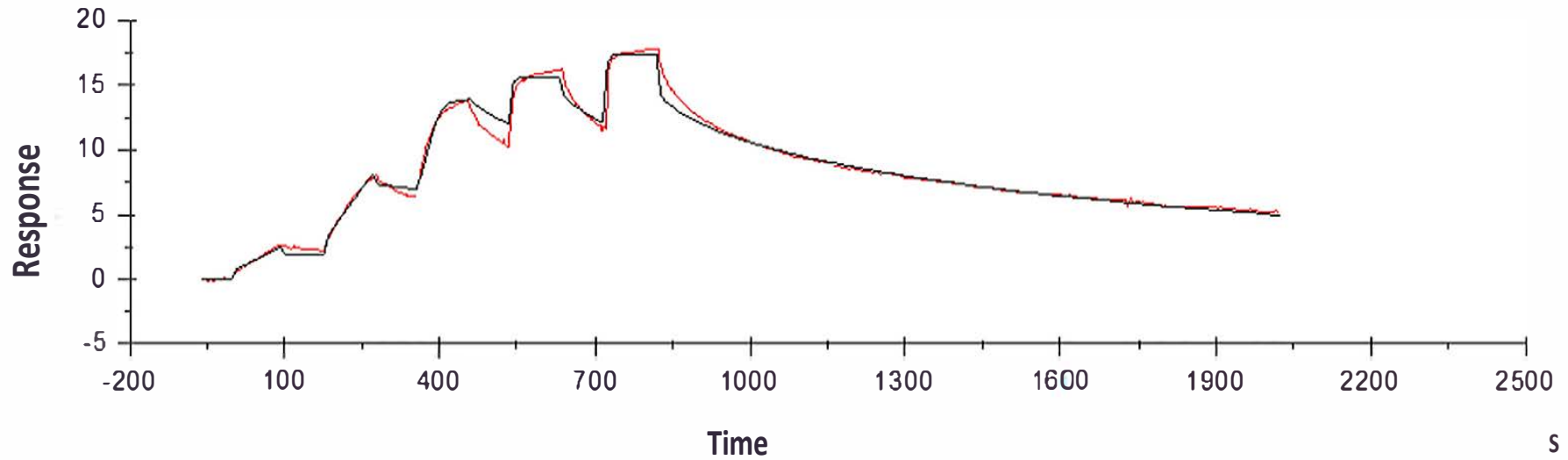




Binding Kinetics of Brigatinib on ALK [I1171T]



Data generated and provided by Carna Biosciences, Inc.