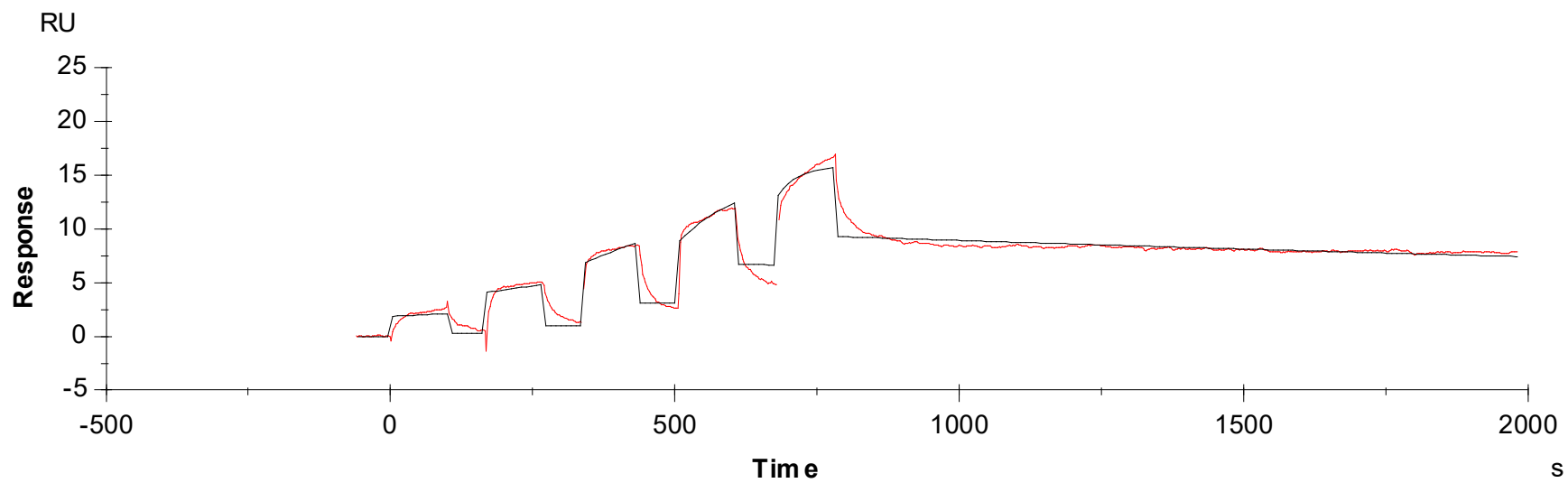


Binding Kinetics of Brigatinib on ALK [G1202R]



Data generated and provided by Carna Biosciences, Inc.

Binding kinetics of lorlatinib on ALK [G1202R]

