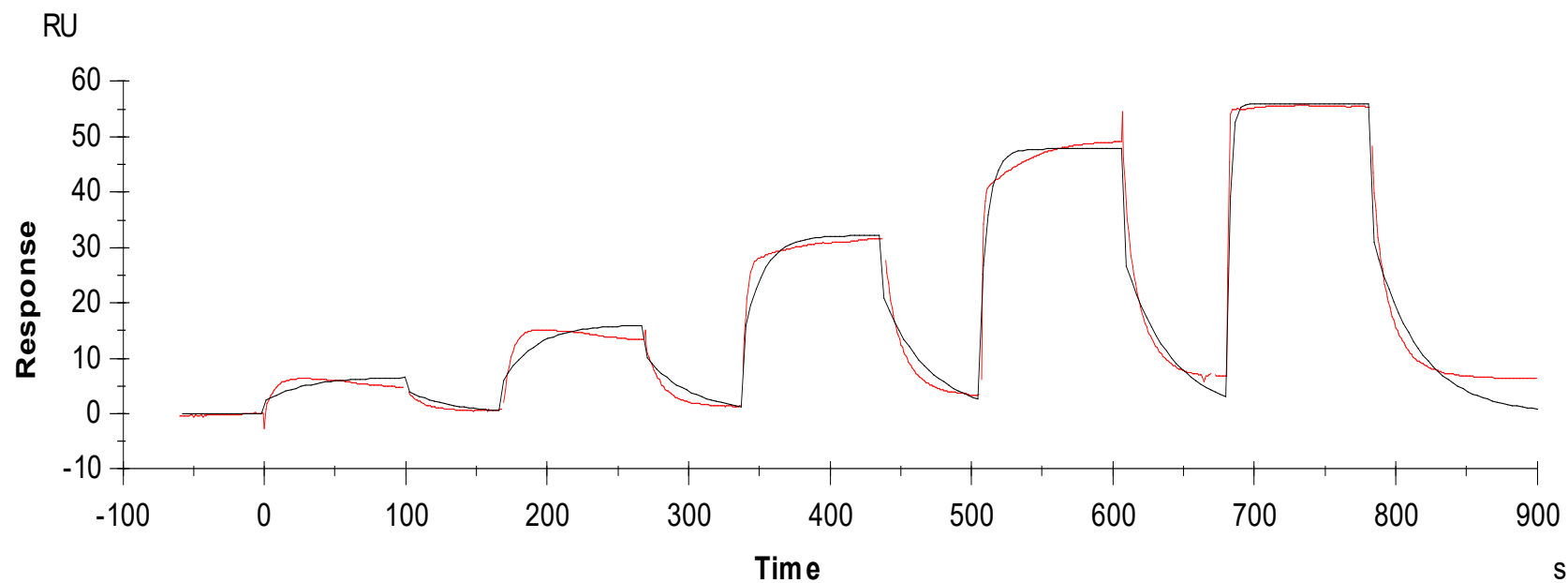


Binding Kinetics of crizotinib on CASK



Data generated and provided by Carna Biosciences, Inc.