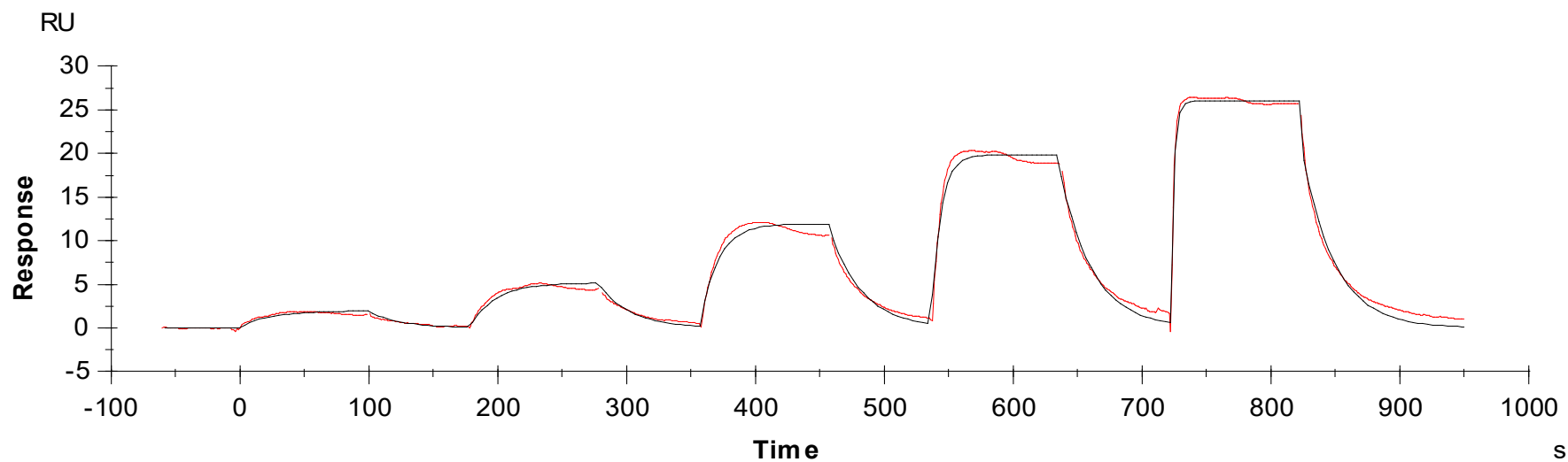




Binding Kinetics of Pazopanib on FGFR2[M538L][non-activated]



Data generated and provided by Carna Biosciences, Inc.