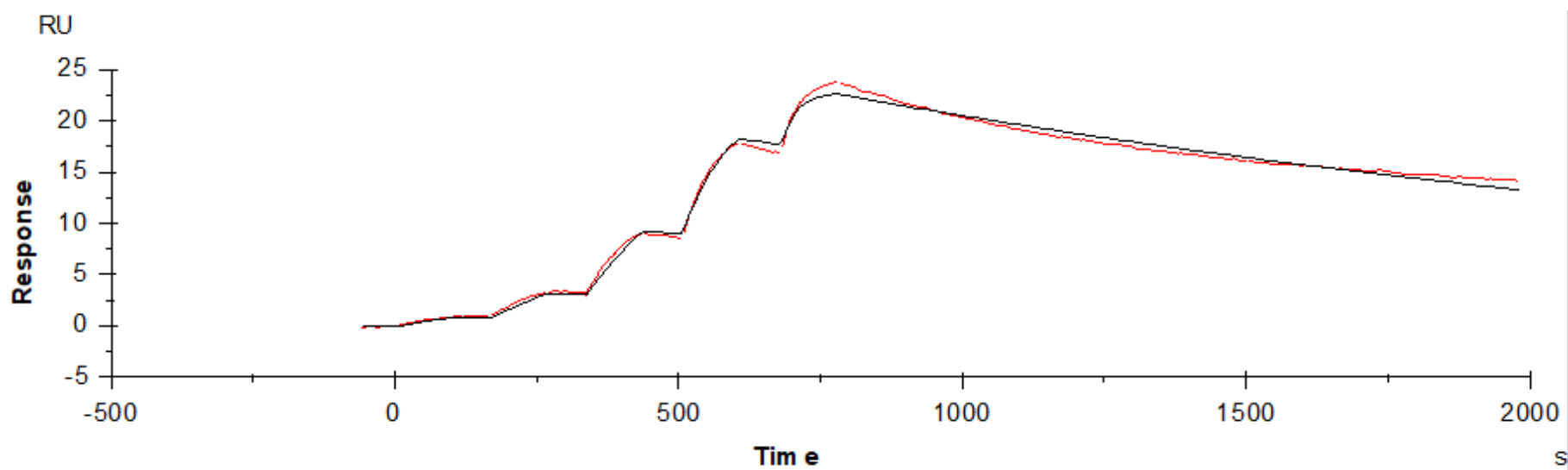
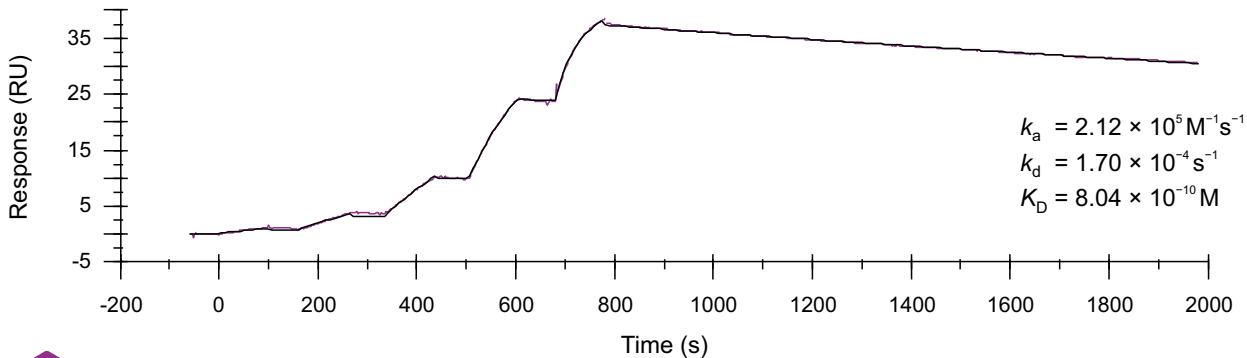


Binding Kinetics of Ponatinib on FGFR2



Data generated and provided by Carna Biosciences, Inc.

Binding kinetics of infigratinib on FGFR2



Oncolines
A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com