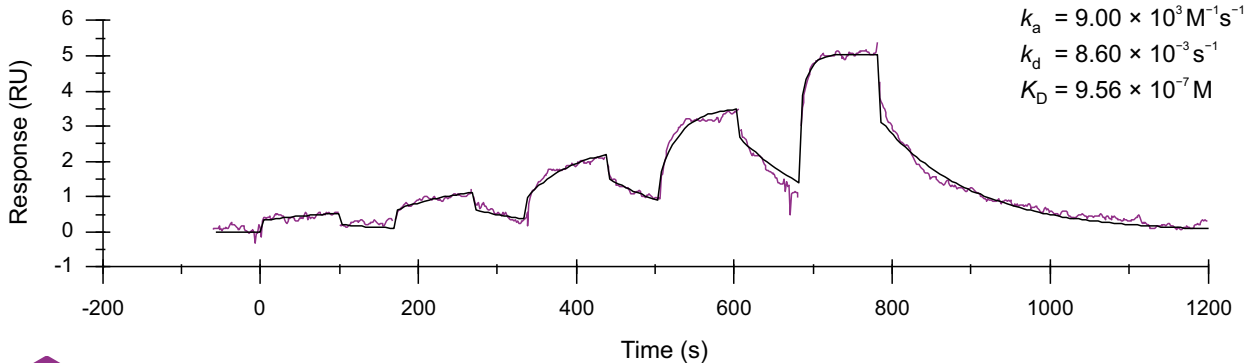


Binding kinetics of ponatinib on BTK non-activated



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com