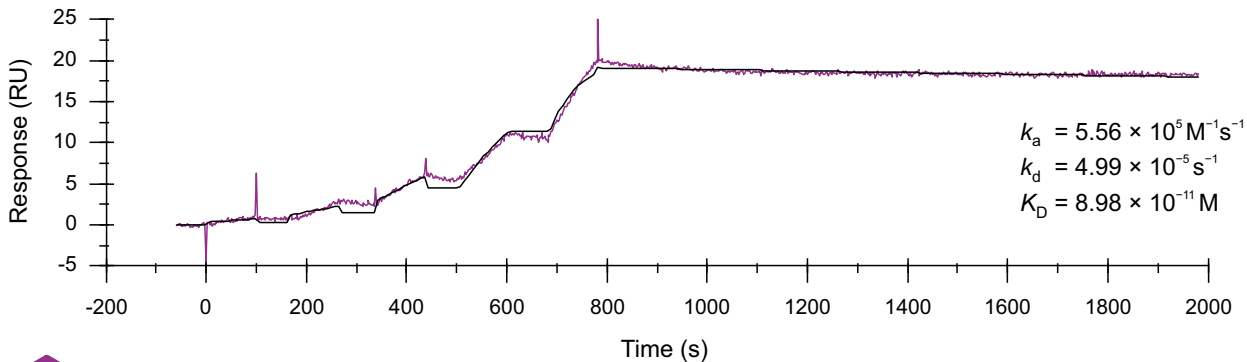


Binding kinetics of ponatinib on LCK



Oncolines
A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com