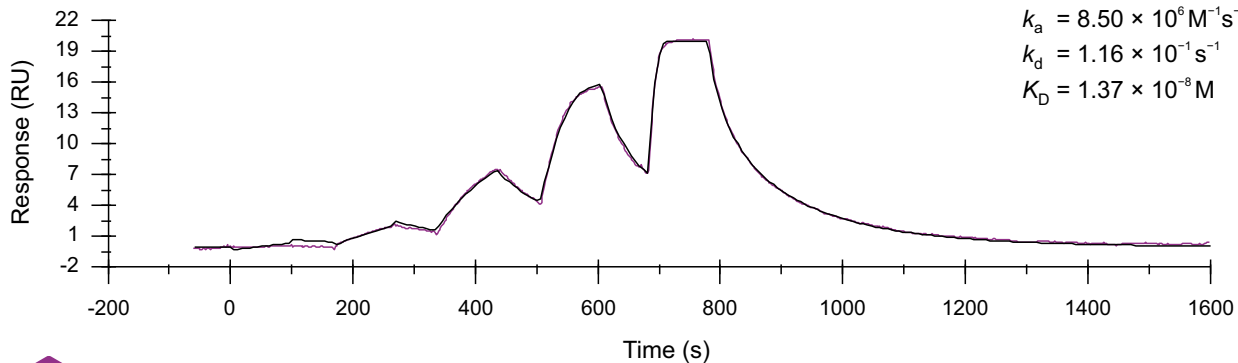


Binding kinetics of crizotinib on FAK (PTK2)



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com