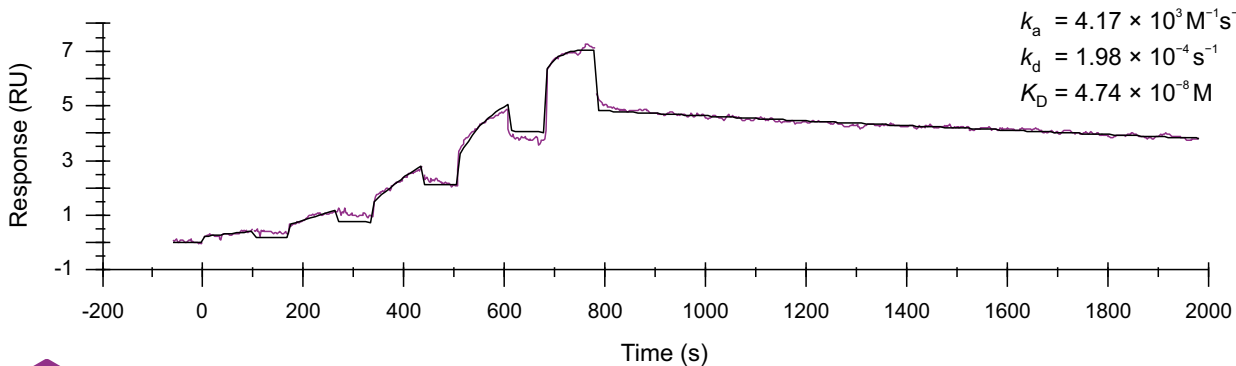


Binding kinetics of pazopanib on KIT



Oncolines
A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com