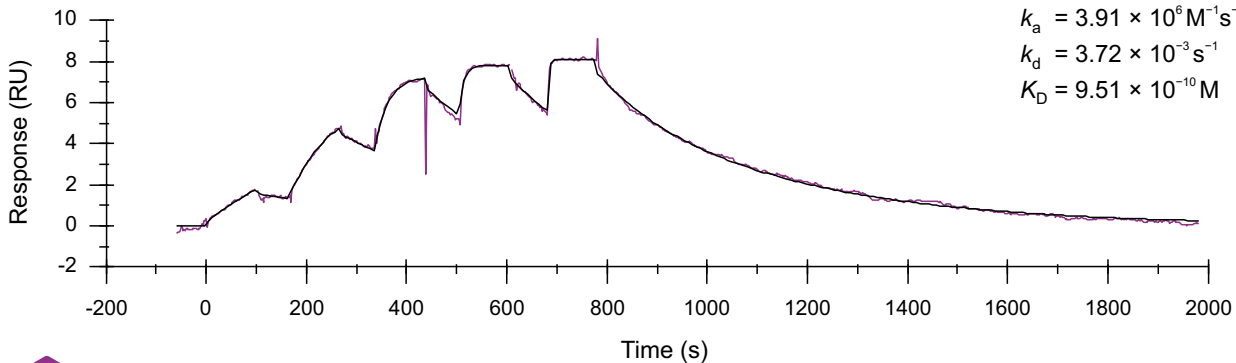


Binding kinetics of pictilisib on PIK3CA/PIK3R1



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com