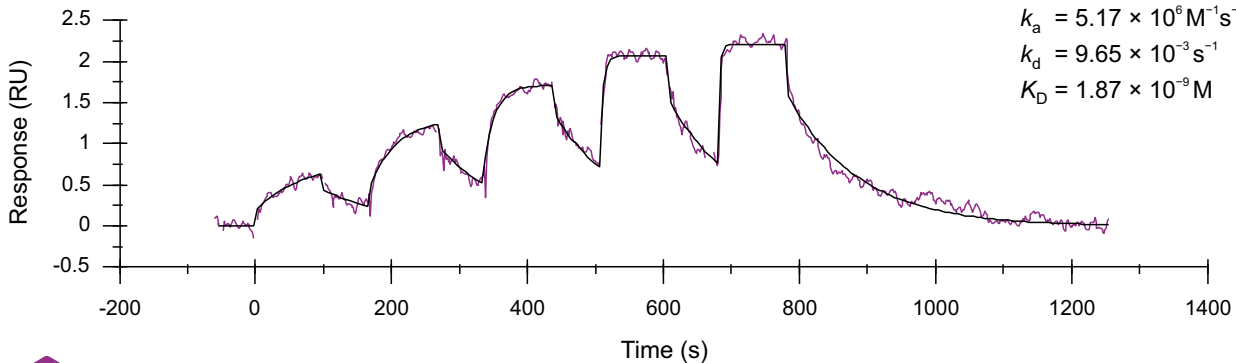


Binding kinetics of afatinib on EGFR [T790M]



Oncolines
A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com