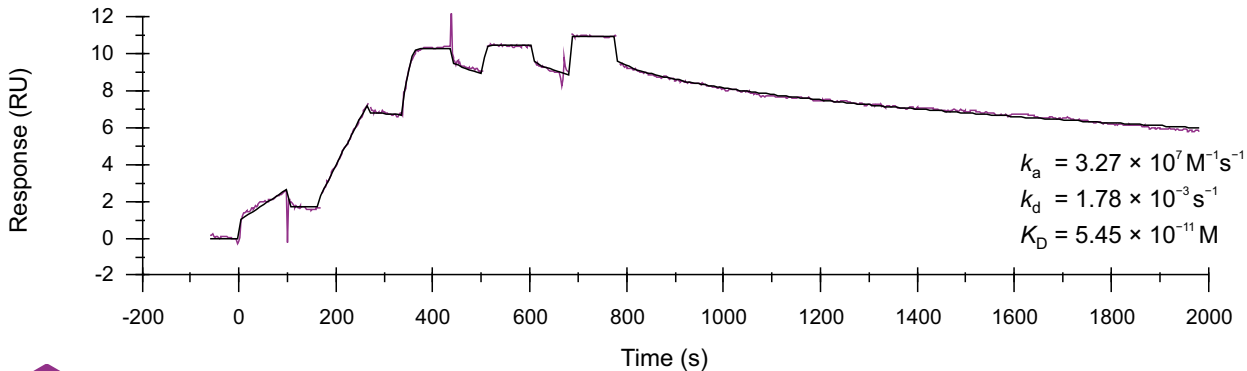


Binding kinetics of gefitinib on EGFR [d746-750/C797S]



Oncolines
A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com