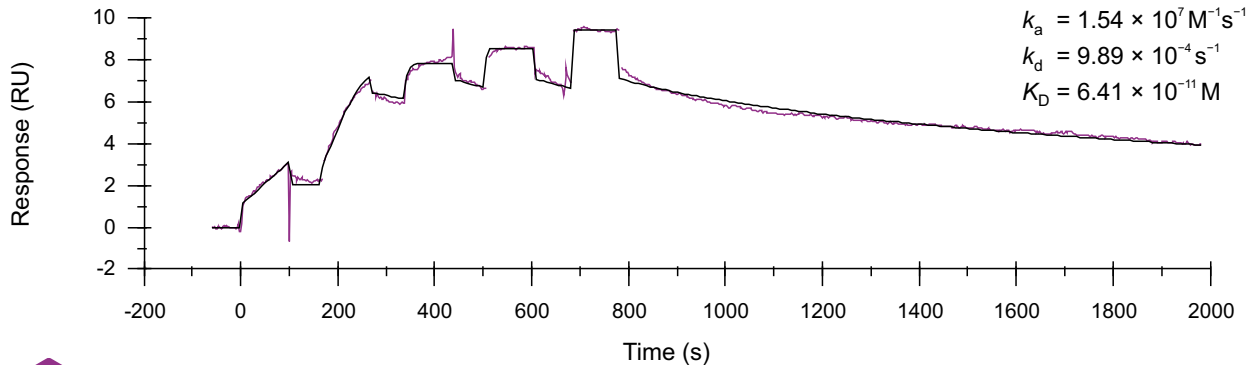


Binding kinetics of gefitinib on EGFR [d746-750]



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com