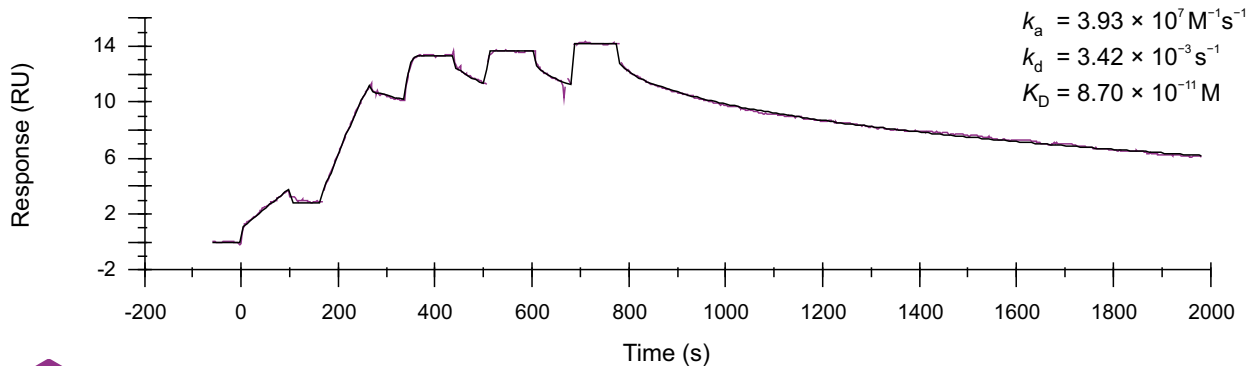


Binding kinetics of gefitinib on EGFR [C797S/L858R]



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com