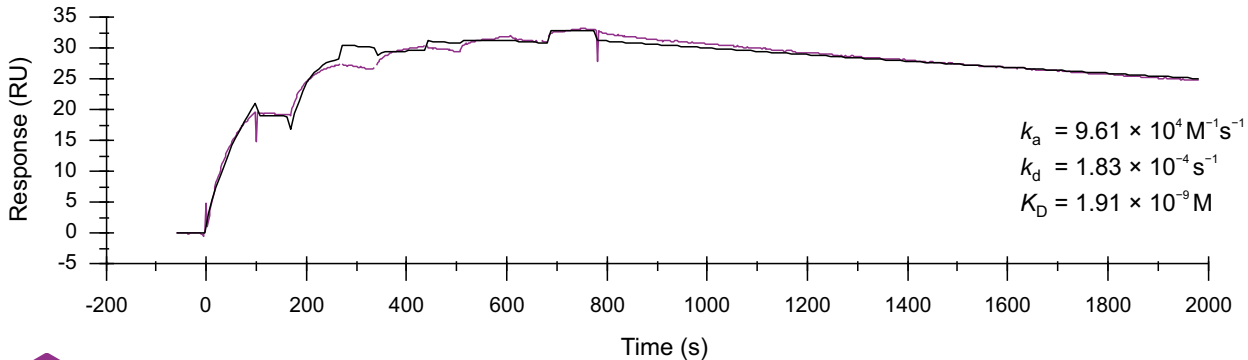


Binding kinetics of BMS-777607 on TYRO3



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com