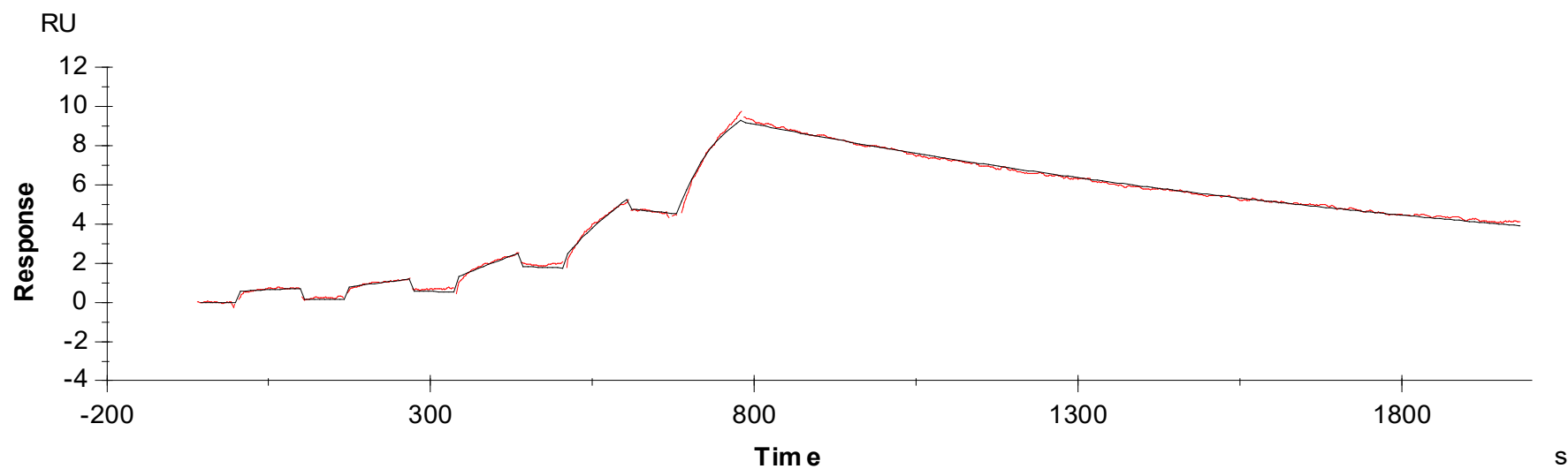
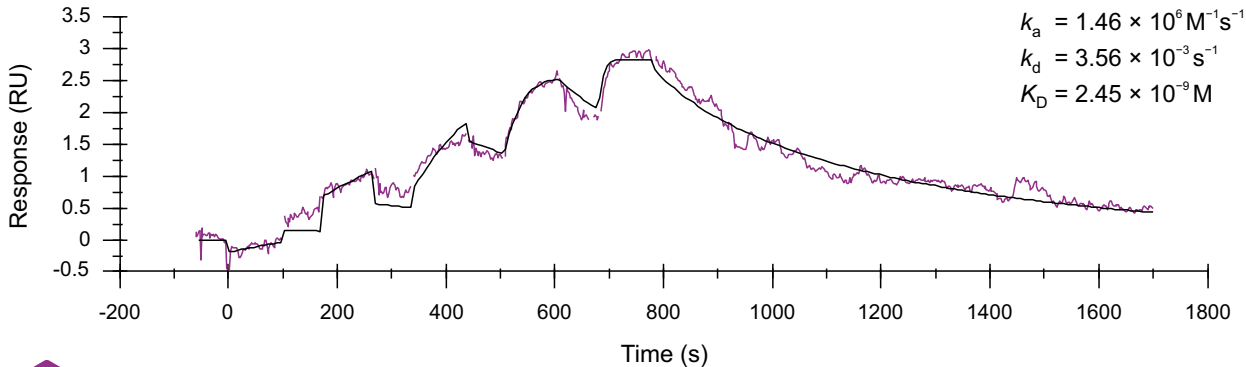


Binding Kinetics of Quizantinib on RET[M918T]



Data generated and provided by Carna Biosciences, Inc.

Binding kinetics of staurosporine on RET [M918T]



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com