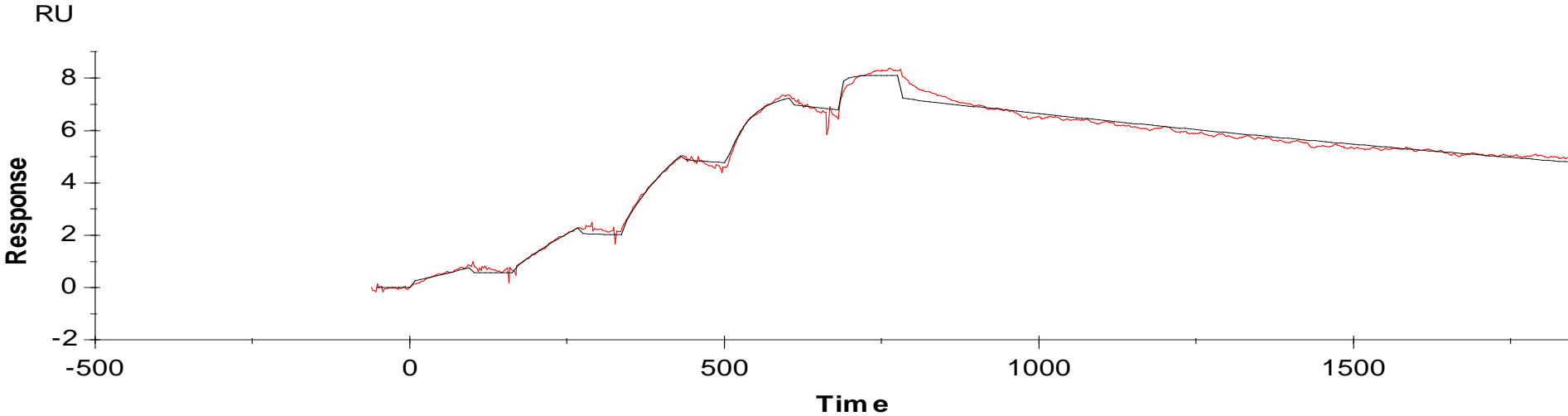
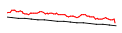


Binding kinetics of dabrafenib on RAF1





—|
2000
s