

## Binding kinetics of dabrafenib on RAF1

