

## Binding kinetics of dabrafenib on BRAF

