

Binding kinetics of dabrafenib on BRAF

