

# Binding kinetics of vemurafenib on BRAF

