

Binding kinetics of sorafenib on RET

