

Binding kinetics of ponatinib on RET

