

Binding kinetics of sorafenib on p38 α (MAPK14)

