

Binding kinetics of sorafenib on p38 α inactive (MAPK14)

