

Binding kinetics of tamatinib on LYNa

