

Binding kinetics of A-674563 on PKACa (PRKACA)

