

Binding kinetics of GSK690693 on AKT2

