

Binding kinetics of GSK690693 on PKN1

