

# Binding kinetics of GSK690693 on PKN1

PKN1 GSK690693

