

Binding kinetics of GSK690693 on PAK4

