

Binding kinetics of A-92 on GCN2

