

Binding kinetics of GSK1070916 on AurA

AurA GSK1070916

