

Binding kinetics of GSK 1070916 on AurB

GSK 1070916 on AURB

