

Binding kinetics of GSK461364 on PLK1

