

Binding kinetics of BI2536 on PLK1

PLK1 BI2536

