

Binding kinetics of sunitinib on IRAK4

IRAK4 Sunitinib

