

Binding kinetics of sunitinib on TNIK

TNIK Sunitinib

