

Binding kinetics of vandetanib on KDR

KDR-Vandetanib

