

Binding kinetics of vandetanib on KDR (VEGFR2)

