

Binding kinetics of vandetanib on TIE2

