

Binding kinetics of nintedanib on PDGFR β (PDGFRB)

