

Binding kinetics of nintedanib on PDGFR β (PDGFRB) non-activated

