

Binding kinetics of nintedanib on PDGFR α (PDGFRA)

