

Binding kinetics of nintedanib on PDGFR α (PDGFRA) non-activated

