

Binding kinetics of cabozantinib on ALK[L1196M]

ALK[L1196M] Cabozantinib

