

Binding kinetics of brigatinib on ALK [L1196M]

