

Binding kinetics of osimertinib on EGFR[T790M/L858R]

