

Binding kinetics of gefitinib on EGFR

