

# Binding kinetics of erlotinib on EGFR

