

# Binding kinetics of sunitinib on ITK

