

# Binding kinetics of pazopanib on KIT

