

# Binding kinetics of AZD1775 on WEE1

