

# Binding kinetics of WZ4003 on AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1

