

# Binding kinetics of dasatinib on TGFB $\beta$ 1

TGFB $\beta$ 1 Dasatinib

