

Binding kinetics of BGJ398 on FGFR1 non activated

FGFR1 non activated BGJ398

