

Binding kinetics of BGJ398 on FGFR1

FGFR1 - BGJ398

