

Binding kinetics of BGJ398 on FGFR4

FGFR4 BGJ398

