

Binding kinetics of sunitinib on FGFR3

FGFR3 - Suninitib

