

Binding kinetics of ponatinib on FYN

FYN - Ponatinib

