

Binding kinetics of ponatinib on SRC

ponatinib SRC

