

Binding kinetics of sorafenib on TRKC

TRKC Sorafenib
RU

