

Binding kinetics of GSK1838705 on IGF1R

