

Binding kinetics of ADW742 on IGF1R

