

Binding kinetics of GSK429286A on ROCK1

ROCK1 GSK429286A

