

Binding kinetics of GSK 429286A on ROCK

