

## Binding kinetics of ATP on DGK $\alpha$

