

Regulation of Cdc2 kinase in Fission Yeast *S. Pombe*

