

Stick Model of β -Adrenergic Receptor Kinase (GRK2) binding to $G\beta\gamma$

