

G α Protein Couples to Adenylate Cyclase

