

G Protein Stimulation of Phospholipase C Isoforms

PLC subtype	Tissue distribution	G protein subunit
PLC- β 1	Most in brain	G α q, G $\beta\gamma$, G α 11 G α 14, G α 16
PLC- β 2	Hemopoietic cells	G α 16, G $\beta\gamma$, G α q
PLC- β 3	Wide	G α q, G $\beta\gamma$, G α 16
PLC- β 4	Retina	G α q
PLC-X β	Frog oocytes	G α o

