

Stick Model of Trimeric G Protein Complex - G_t

