

$G_{s\alpha L}$ -Leu²²⁷

stimulatory heterotrimeric G-protein, long splice variant of the α -subunit
rat, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-504	1 ml

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Membrane suspension. Supplied in 75 mM Tris-HCl, pH 7.4, 12.5 mM MgCl₂, and 1 mM EDTA.

Protein concentration

1.3 mg/ml

Description

$G_{s\alpha L}$ is the long splice variant of the α -subunit of stimulatory heterotrimeric G_s-proteins. It differs from the short splice variant ($G_{s\alpha S}$) by 15-amino acid insert between the Ras-like domain and the α -helical domain.

$G_{s\alpha L}$ activates adenylate cyclase (AC) and possesses a lower GDP-affinity than $G_{s\alpha S}$ (cat.# PR-505).

These differences in GDP-affinity have important consequences for receptor/G-protein coupling and AC activation.

The exchange of Gln²²⁷ to Leu²²⁷ inhibits the intrinsic GTPase activity, resulting in a constitutively active G_{α} , and increases GDP-affinity of G_{α} .

Selected References:

Graziano et al. (1989) Expression of $G_{s\alpha}$ in Escherichia coli. Purification and properties of two forms of the protein. *J. Biol. Chem.* **264**:409.

Yu, B. and Simon, M. (1998) Interaction of the Xanthine Nucleotide Binding $G_{0\alpha}$ Mutant with G Protein-coupled Receptors. *J. Biol. Chem.* **273**:30183.

Gille, A. and Seifert, R. (2003) 2'-(3')-O-(N-Methylantraniloyl)-substituted GTP Analogs: A Novel Class of Potent Competitive Adenylyl Cyclase Inhibitors. *J. Biol. Chem.* **278**:12672.

Gille et al. (2003) GDP Affinity and Order State of the catalytic Site Are Critical for Function of Xanthine Nucleotide-selective $G_{\alpha S}$ Proteins. *J. Biol. Chem.* **278**:7822.

Associated products available from Jena Bioscience

Non-hydrolyzable GTP-analogs (such as GTP γ S, GppCp, GppNHp, NPE-caged-GTP...)

Fluorescent GTPs (such as MANT, ANT, TNP, or GTP labeled with a dye of your choice)

Recombinant G-protein signaling reagents including GAPs, GEFs, small GTPases, and many more...

For detailed information please view the sections on

G_{sαL}-Leu²²⁷

stimulatory heterotrimeric G-protein, long splice variant of the α-subunit
rat, recombinant, Sf9 insect cells

<http://www.jenabioscience.com>