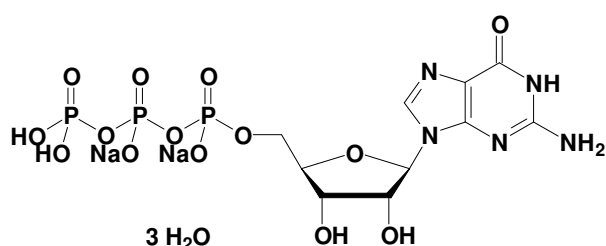


GTP, lyophilized

Guanosine 5'-triphosphate, sodium salt

Cat. No.	Amount
NU-1012-100	100 mg, lyophilized
NU-1012-1G	1 g, lyophilized
NU-1012-10G	10 g, lyophilized
NU-1012-100G	100 g, lyophilized



Molecular Formula: $C_{10}H_{14}N_5Na_2O_{14}P_3 \times 3H_2O$

Molecular Weight: 621.18

Absorbance

absorbance max: 252 nm (pH 7)
 ϵ at absorbance max: $14.2 \text{ mmol}^{-1} \text{ cm}^{-1}$

Purity: >90%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at $< -20^\circ\text{C}$.

For research use only!

Applications:

Assembly of ribosomal units^[1]

Microdomain formation by small GTPases^[2]

Antiviral activity of large GTPases (dynamin superfamily) ^[3]

Regulation of exocytosis by Rho GTPases^[4]

Mechanism of hydrolysis by ADP-ribosylation factors^[5]

Specific Ligands:

Guanylate binding proteins^[6]

Yeast septins^[7]

Selected References:

[1] Blombach *et al.* (2011) Assembling the archeal ribosome: roles for transition factor-related GTPases. *Biochemical Society Transactions* **39**:45.

[2] Stuermer (2011) Microdomain-forming proteins and the role of the reggies/flotillins during axon regeneration in zebrafish. *Biochimica Biophysica Acta, Molecular Basis of Disease* **1812**:415.

[3] Haller *et al.* (2011) Human MxA protein: An Interferon-induced Dynamin-like GTPase with broad antiviral activity. *J. Interferon and Cytokine Research* **31**:79.

[4] Stephane *et al.* (2011) Rho GTPases and exocytosis: what are the molecular links? *Seminars in Cell and Developmental Biology* **22**:27.

[5] East *et al.* (2011) Models for the function of Arf GAPs. *Seminars in Cell and Developmental Biology* **22**:3.

[6] Vestal *et al.* (2011) The guanylate binding proteins: Emerging insights into the biochemical properties and functions of this family of large interferon-induced guanosine triphosphatase. *J. Interferon and Cytokine Research* **31**:89.

[7] Younghoon *et al.* (2011) Septin structure and function in yeast and beyond. *Trends in Cell Biology* **21**:141.