

PRODUCT INFORMATION



8-Bromoguanosine-5'-O-triphosphate (sodium salt)

Item No. 38551

Formal Name: ((2R,3S,4R,5R)-5-(2-amino-8-bromo-6-oxo-3,6-dihydro-9H-purin-9-yl)-3,4-dihydroxytetrahydrofuran-2-yl)methyl triphosphate, tetrasodium salt

Synonym: 8-bromo GTP

MF: C₁₀H₁₁BrN₅O₁₄P₃ • 4Na

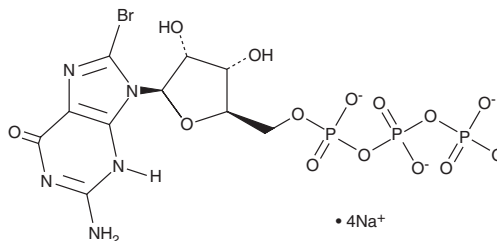
FW: 690.0

Purity: ≥95%

Supplied as: A solution in water

Storage: -80°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

8-Bromoguanosine-5'-O-triphosphate is a derivative of the energy substrate for protein synthesis and gluconeogenesis, guanosine 5'-triphosphate (GTP; Item No. 16060).^{1,2} It inhibits the *E. coli* GTPase FtsZ ($K_i = 31.8 \mu\text{M}$) but promotes the assembly of porcine brain microtubules in a cell-free assay when used at a concentration of 500 μM .

References

1. Lappchen, T., Hartg, A.F., Pinas, V.A., *et al.* GTP analogue inhibits polymerization and GTPase activity of the bacterial protein FtsZ without affecting its eukaryotic homologue tubulin. *Biochemistry* **44**(21), 7879-7884 (2005).
2. Muraoka, M., Fukuzawa, H., Nishida, A., *et al.* The effects of various GTP analogues on microtubule assembly. *Cell Struct. Funct.* **24**(2), 101-109 (1999).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/23/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM