

PRODUCT INFORMATION



TNP-β-L-GTP (sodium salt)

Item No. 38830

Formal Name: ((3a'S,4'S,6'S,6a'S)-4'-(2-amino-6-oxo-3,6-dihydro-9H-purin-9-yl)-2,4,6-trinitro-3a',4',6',6a'-tetrahydrospiro[cyclohexane-1,2'-furo[3,4-d][1,3]dioxole]-2,5-dien-6'-yl)methyl triphosphate, tetrasodium salt
Synonym: TNP-β-L-Guanosine 5'-triphosphate

MF: C₁₆H₁₃N₈O₂₀P₃ • 4Na

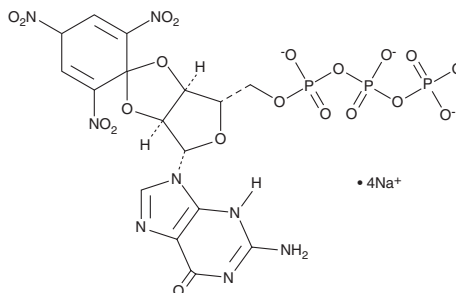
FW: 822.2

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

TNP-β-L-GTP is an enantiomer of the fluorescent GTP derivative TNP-GTP (Item No. 38570). A related derivative, TNP-GTP (Item No. 38570), displays an emission maximum of 552 nm upon excitation at 410 nm.¹

Reference

1. Hiratsuka, T. A chromophoric and fluorescent analog of GTP, 2',3'-O-(2,4,6-trinitrocyclohexadienylidene)-GTP, as a spectroscopic probe for the GTP inhibitory site of liver glutamate dehydrogenase. *J. Biol. Chem.* **260**(8), 4784-4790 (1985).

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, 05/05/2026

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