

# PRODUCT INFORMATION



## Guanosine-5'-O-triphosphoro-γ-1-(5-sulfonic acid)naphthylamidate (sodium salt)

Item No. 38687

**Formal Name:** P'→P-anhydride with 5-(phosphonoamino)-1-naphthalenesulfonic acid guanosine 5'-(trihydrogen diphosphate), trisodium salt

**Synonym:** GTP-γ-AmNS

**MF:** C<sub>20</sub>H<sub>20</sub>N<sub>6</sub>O<sub>16</sub>P<sub>3</sub>S • 3Na

**FW:** 794.4

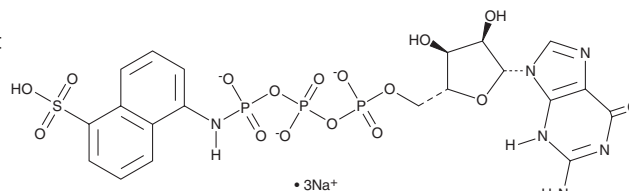
**Purity:** ≥95%

**EX./EM. Max:** 340/460 nm

**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

Guanosine-5'-O-triphosphoro-γ-1-(5-sulfonic acid)naphthylamidate (GTP-γ-AmNS) is a fluorescent derivative of GTP (Item No. 16060).<sup>1,2</sup> It has been used as a substrate for snake venom phosphodiesterases (PDEs) and to monitor tubulin assembly in cell-free assays. GTP-γ-AmNS displays an emission maximum of 444 nm upon excitation at 320 nm, which is quenched by 1.2-1.4-fold upon cleavage of the α-β phosphoryl bond.<sup>1</sup>

### References

1. Pollack, S.E. and Auld, D.S. Fluorescent nucleotide triphosphate substrates for snake venom phosphodiesterase. *Anal. Biochem.* **127**(1), 81-88 (1982).
2. Yarbrough, L.R. and Kirsch, M. Binding of fluorescent analogs of GTP to the exchangeable nucleotide binding site of tubulin. *J. Biol. Chem.* **256**(1), 112-117 (1981).

#### 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, 04/24/2026

#### 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](http://WWW.CAYMANCHEM.COM)