

PRODUCT INFORMATION

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

Item No. 38574

Formal Name: 5-[[hydroxy[[hydroxy(phosphonooxy)phosphinyl]oxy]phosphinyl]amino]-1-naphthalenesulfonic acid, P''→5'-ester with uridine, trisodium salt

Synonym: UTP-γ-AmNS

MF: C₁₉H₁₉N₃O₁₇P₃S • 3Na

FW: 755.3

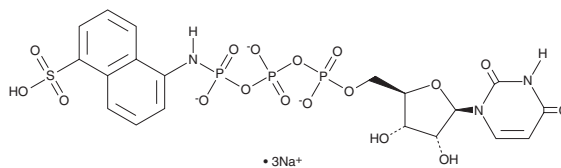
Purity: ≥95%

Ex./Em. Max: 360/465 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

Uridine-5'-O-triphosphoro-γ-1-(5-sulfonic acid)naphthylamidate (UTP-γ-AmNS) is a fluorescent derivative of the nucleotide and dual purinergic P2Y₂ and P2Y₄ receptor agonist UTP (Item No. 9003530).¹ It has been used in high-throughput screening (HTS) assays for inhibitors of bacterial RNA polymerase activity. Upon incorporation of UMP in RNA, γ-AmNS-PPi is released, which has an emission at 465 nm upon excitation at 360 nm and a higher fluorescence intensity than UTP-γ-AmNS.

Reference

1. Bhat, J., Rane, R.A., Solapure, S.M., *et al.* High-throughput screening of RNA polymerase inhibitors using a fluorescent UTP analog. *J. Biomol. Screen.* **11**(8), 968-976 (2006).

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