

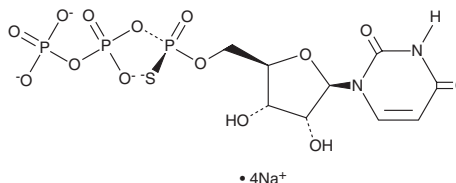
PRODUCT INFORMATION



Sp-Uridine-5'-O-(1-thiotriphosphate) (sodium salt)

Item No. 38573

Synonym: Sp-UTP- α -S
MF: $C_9H_{11}N_2O_{14}P_3S \cdot 4Na$
FW: 588.1
Purity: $\geq 95\%$
Supplied as: A 10 mM solution in water
Storage: $-80^\circ C$
Stability: ≥ 4 years



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

Description

Sp-Uridine-5'-O-(1-thiotriphosphate) (Sp-UTP- α -S) is a nucleotide agonist of the purinergic Y2 (P2Y₂) and P2Y₄ receptors.¹ It induces inositol phosphate accumulation in 1321N1 astrocytoma cells expressing the purinergic P2Y₂ or P2Y₄ receptors (EC₅₀s = 14 and 81 μM , respectively).

Reference

1. Jacobson, K.A., Costanzi, S., Ivanov, A.A., *et al.* Structure activity and molecular modeling analyses of ribose- and base-modified uridine 5'-triphosphate analogues at the human P2Y₂ and P2Y₄ receptors. *Biochem. Pharmacol.* **71**(4), 540-549 (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, 06/25/2026

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