

PRODUCT INFORMATION



Uridine-5'-O-(3-thiotriphosphate) (sodium salt)

Item No. 38692

Formal Name: uridine 5'-(trihydrogen diphosphate), P'-anhydride with phosphorothioic acid, tetrasodium salt

Synonym: UTP- γ -S

MF: C₉H₁₁N₂O₁₄P₃S • 4Na

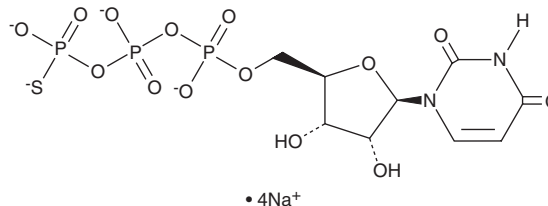
FW: 588.1

Purity: \geq 95%

Supplied as: A solution in water

Storage: -80°C

Stability: \geq 2 years



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

Description

Uridine-5'-O-(3-thiotriphosphate) (UTP- γ -S) is a hydrolysis-resistant derivative of UTP (Item No. 9003530) an agonist of the purinergic P2Y₂ receptor.¹ It induces inositol phosphate formation in 1321N1 astrocytoma cells expressing the human P2Y₂ receptor (EC₅₀ = 240 nM) and chloride secretion in primary nasal epithelial cells from patients with cystic fibrosis in a concentration-dependent manner. UTP- γ -S induces vasoconstriction in isolated human coronary arteries (EC₅₀ = 25.1 μ M).²

References

1. Lazarowski, E.R., Watt, W.C., Stutts, M.J., *et al.* Enzymatic synthesis of UTP γ S, a potent hydrolysis resistant agonist of P2U-purinoceptors. *Br. J. Pharmacol.* **117(1)**, 203-209 (1996).
2. Malmsjö, M., Hou, M., Harden, T.K., *et al.* Characterization of contractile P2 receptors in human coronary arteries by use of the stable pyrimidines uridine 5'-O-thiodiphosphate and uridine 5'-O-3-thiotriphosphate. *J. Pharmacol. Exp. Ther.* **293(3)**, 755-760 (2000).

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, 08/22/2023

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