

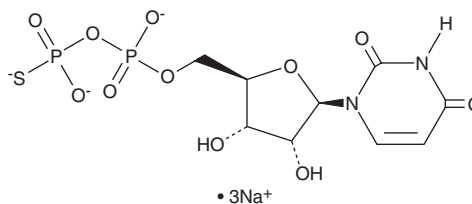
PRODUCT INFORMATION



UDP-β-S (sodium salt)

Item No. 38493

Formal Name:	5'-uridylic acid, monoanhydride with phosphorothioic acid, trisodium salt
Synonym:	Uridine-5' O-(2-thiodiphosphate)
MF:	C ₉ H ₁₁ N ₂ O ₁₁ P ₂ S • 3Na
FW:	486.2
Purity:	≥95%
Supplied as:	A solution in water
Storage:	-80°C
Stability:	≥2 years



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

Description

UDP-β-S is an agonist of the purinergic P2Y₆ receptor.¹ It selectively induces inositol phosphate accumulation in 1321N1 cells expressing P2Y₆ receptors (EC₅₀ = 28 nM) over those expressing P2Y₁, P2Y₂, or P2Y₄ receptors at 10 μM. UDP-β-S (1 μM) induces DNA synthesis in isolated rat aorta smooth muscle cells.

Reference

1. Hou, M., Harden, T.K., Kuhn, C.M., *et al.* UDP acts as a growth factor for vascular smooth muscle cells by activation of P2Y₆ receptors. *Am. J. Physiol. Heart Circ. Physiol.* **282**(2), H784-H792 (2002).

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, 06/08/2023

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