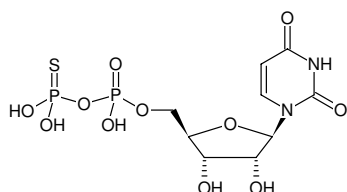


**UDPβS**

Uridine-5'-(β-thio)-diphosphate, Sodium salt

Cat. No.	Amount
NU-442-5	5 mg



Structural formula of UDPβS

**For research use only!****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C<sub>9</sub>H<sub>14</sub>N<sub>2</sub>O<sub>11</sub>P<sub>2</sub>S (free acid)**Molecular Weight:** 420.22 g/mol (free acid)**Exact Mass:** 419.98 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 10.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Specific Ligands:**Ligand for purinergic receptors:Selective agonist at P2Y<sub>6</sub><sup>[1,2]</sup> receptors

**Please note:** For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.

**Selected References:**

[1] Malmjö *et al.* (2003) The stable pyrimidines UDPbetaS and UTPgammaS discriminate between contractile cerebrovascular P2 receptors. *J. Pharmacol.* **458** (3):305.

[2] Malmjö *et al.* (2000) 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.