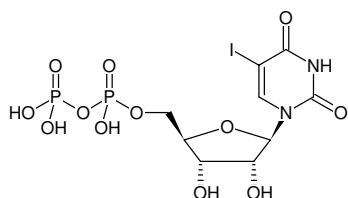


**5-Iodo-UDP**

(5I-UDP)

5-Iodo-uridine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-867S	50 µl (10 mM)
NU-867L	5 x 50 µl (10 mM)



Structural formula of 5-Iodo-UDP

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: 12 months after date of delivery**Molecular Formula:** C₉H₁₃N₂O₁₂P₂I (free acid)**Molecular Weight:** 530.06 g/mol (free acid)**Exact Mass:** 529.90 g/mol (free acid)**CAS#:** 34198-43-1**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 287 nm, ε 7.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Specific Ligands:**Selective agonistic ligand for P2Y₆ receptor^[1,2,3]**Selected References:**

[1] Besada *et al.* (2006) Structure-activity relationships of uridine-5'-diphosphate analogues at the human P2Y₆ receptor. *J. Med. Chem.* **49** (18):5532.

[2] Mamedova *et al.* (2008) Attenuation of apoptosis in vitro and ischemia/reperfusion injury in vivo in mouse skeletal muscle by P2Y₆ receptor activation. *Pharmacol. Res.* **58** (3-4):232.

[3] Jacobson *et al.* (2009) Development of selective agonists and antagonists of P2Y receptors. *Purinergic Signalling* **5**:75.