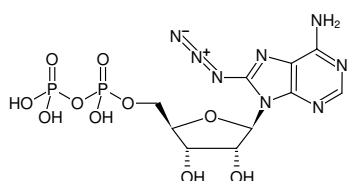


**8-Azido-ADP**

8-Azido-adenosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-159S	500 µl (10 mM)
NU-159L	5 x 500 µl (10 mM)



Structural formula of 8-Azido-ADP

For research use only!**Shipping:** shipped on blue ice**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:** C₁₀H₁₄N₈O₁₀P₂ (free acid)**Molecular Weight:** 468.21 g/mol (free acid)**Exact Mass:** 468.03 g/mol (free acid)**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} 281 nm, ε 13.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Marti *et al.* (1997) 8-Azido-nucleotides as substrates of Torpedo electric organ apyrase. effect of photoactivation on apyrase activity. *Brain Res. Bull.* **44** (6):695.

Sloothaak *et al.* (1985) The use of 8-azido-ATP and 8-azido-ADP as photoaffinity labels of the ATP synthase in submitochondrial particles: evidence for a mechanism of ATP hydrolysis involving two independent catalytic sites? *Biochim. Biophys. Acta.* **809** (1):27.

Wagenvoord *et al.* (1979) Localisation of adenine nucleotide-binding sites on beef-heart mitochondrial ATPase by photolabelling with 8-azido-ADP and 8-azido-ATP. *Biochim. Biophys. Acta.* **548** (1):85.