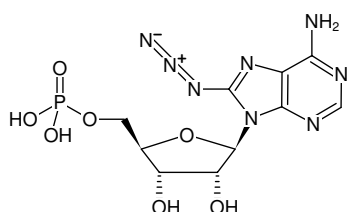


**8-Azido-AMP**

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

Cat. No.	Amount
NU-158S	500 µl (10 mM)
NU-158L	5 x 500 µl (10 mM)



Structural formula of 8-Azido-AMP

Selected References:

Cheunget *al.* (2000) Characterization of AMP-activated protein kinase gamma-subunit isoforms and their role in AMP binding. *Biochem J.* **346**:659.

Marcus *et al.* (1979) Inhibition of fructose-1,6-biphosphatase by the photoaffinity AMP analog, 8-azidoadenosine 5'-monophosphate. *J. Biol. Chem.* **254** (2):259.

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:** C₁₀H₁₃N₈O₇P (free acid)**Molecular Weight:** 388.23 g/mol (free acid)**Exact Mass:** 388.06 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**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)