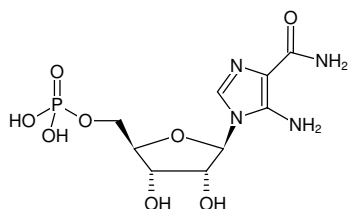


**ZMP**

(AICAR monophosphate)

5'-Aminoimidazole-4-carboxamide-1-β-D-ribofuranosyl-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1193S	50 µl (10 mM)
NU-1193L	5 x 50 µl (10 mM)



Structural formula of ZMP

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:** C₉H₁₅N₄O₈P (free acid)**Molecular Weight:** 338.21 g/mol (free acid)**Exact Mass:** 338.06 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} 265 nm, ε 12.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Corton *et al.* (1995) 5-Aminoimidazole-4-carboxamide ribonucleoside: a specific method for activating protein kinase in intact cells? *Eur J Biochem.* **229**:558.Henin *et al.* (1995) Inhibition of fatty acid and cholesterol synthesis by stimulation of AMP-activated protein kinase. *FASEB J.* **9**:54.Henin *et al.* (1996) Stimulation of rat liver AMP-activated protein kinase by AMP analogues. *Biochim Biophys Acta.* **1290**:197.Sullivan *et al.* (1994) Inhibition of lipolysis and lipogenesis in isolated rat adipocytes with AICAR, a cell-permeable activator of AMP-activated protein kinase. *FEBS Lett.* **353**:33.