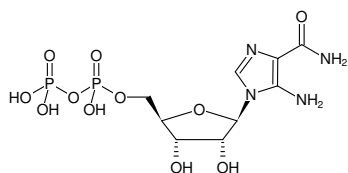


**ZDP**

(AICAR diphosphate)

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

Cat. No.	Amount
NU-1167S	20 µl (10 mM)
NU-1167L	5 x 20 µl (10 mM)



Structural formula of ZDP

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.

For general laboratory use.**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:** C₉H₁₆N₄O₁₁P₂ (free acid)**Molecular Weight:** 418.19 g/mol (free acid)**Exact Mass:** 418.03 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} 265 nm, ε 12.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)