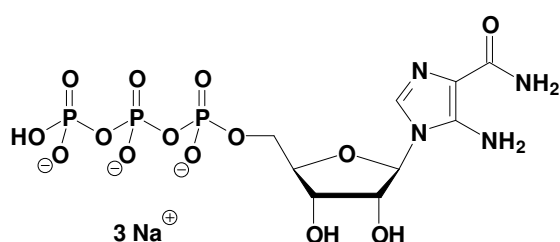


ZTP

**5'-Aminoimidazole-4-Carboxamide-1- β -D-Ribofuranosyl-5'-triphosphate,
Sodium salt
(AICAR triphosphate)**

Cat. No.	Amount
NU-1166S	50 Units
NU-1166L	250 Units



Cat. No.: NU-1166

Molecular Formula: C₉H₁₄N₄O₁₄P₃ (Anion)

Molecular Weight: 495.14 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

* 1 unit = 1 μ l of a 10 mM solution

For research use only!

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