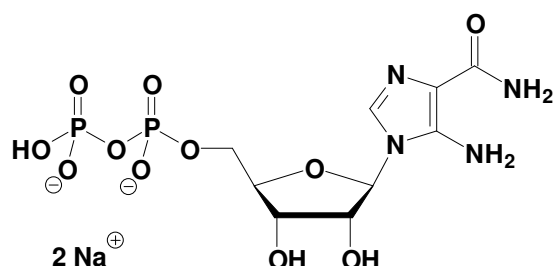


## ZDP

**5'-Aminoimidazole-4-Carboxamide-1- $\beta$ -D-Ribofuranosyl-5'-diphosphate,  
Sodium salt  
(AICAR diphosphate)**

Cat. No.	Amount
NU-1167S	20 Units
NU-1167L	100 Units



### 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

**Cat. No.:** NU-1167

**Molecular Formula:** C<sub>9</sub>H<sub>14</sub>N<sub>4</sub>O<sub>11</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 416.17 (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  $\mu$ l of a 10 mM solution

**For research use only!**