

# Z (AICAR) Kit

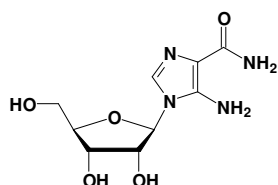
Cat.No.: NK-112

## Kit Contents

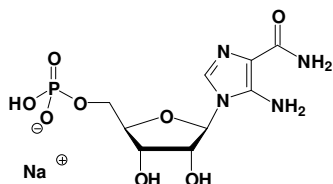
Z (AICAR) Analogs	Cat. No.	Amount
Z (AICAR)	N-1051	1 mg
ZMP (AICAR monophosphate)	NU-1193	50 Units
ZDP (AICAR diphosphate)	NU-1167	50 Units
ZTP (AICAR triphosphate)	NU-1166	50 Units

1 Unit = 1  $\mu$ l of a 10 mM solution

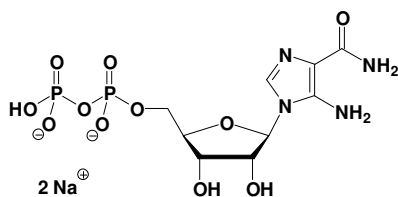
## Structures



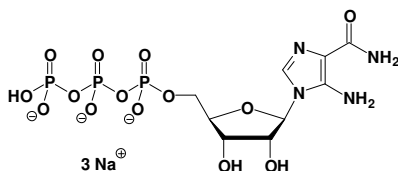
Z  
(AICAR)



ZMP  
(AICAR  
monophosphate)



ZDP  
(AICAR  
diphosphate)



ZTP  
(AICAR  
triphosphate)

## Introduction

Z, (AICAR) analogs are AMP-activated protein kinase activators.

## Kit Description

The Z (AICAR) Kit contains a set of 4 typical Z (AICAR) analogs (Nucleoside, 5'-Mono, Di and Triphosphate).

## 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.

## Storage and Stability

Short term exposure (up to 1 week cumulative) to ambient temperature is possible. Long term storage is recommended at  $< -20^{\circ}\text{C}$ . If properly stored, Jena Bioscience guarantees optimal performance of the compounds for 12 months after date of delivery.