

AICAR is a cell-permeable activator of AMP-activated protein kinase (AMPK).

A cell-permeable nucleoside compound whose phosphorylated metabolite activates adenosine monophosphate-activated protein kinase (AMPK).

Acts as a regulator of *de novo* purine synthesis.

Stimulates glucose uptake in perfused and isolated muscle.

Offers protection against cell death induced by glucose deprivation, chemical hypoxia, and exposure to glutamate and Ab peptide.

*BLG-A103-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

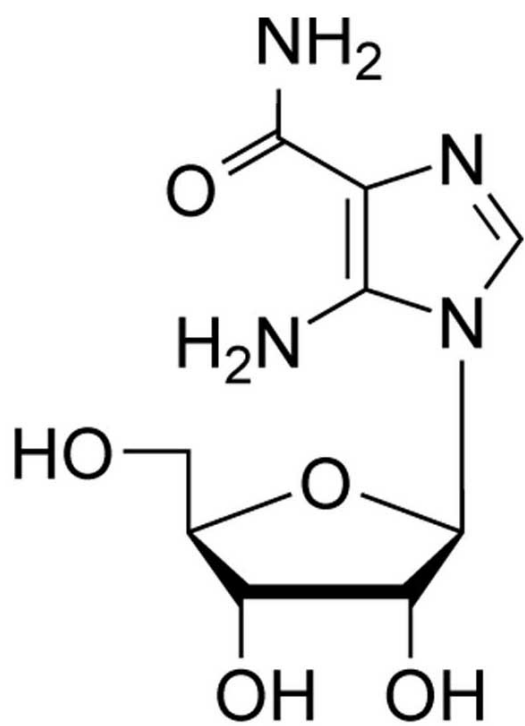
## Ordering Information

[Order Online »](#)

BLG-A103-50	5x10 $\mu$ mol
-------------	----------------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 5-Aminoimidazole-4-carboxamide-1-β-D-ribofuranoside, Acadesine, Z-riboside

**CAS** 2627-69-2

**Couple Type** Modified nucleotides

**Formula**  $C_9H_{14}N_4O_5$

**MW** 258.2

**Purity** > 98% HPLC

**Quantity** 10μmol≈2.6mg

**Solubility** Soluble in water.

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)