

AICAR

AMPK activator

An adenosine analog that is phosphorylated in whole cells to form 5-aminoimidazole-4-carboxamide-1-D-ribofuranosyl-5'-monophosphate (ZMP), which stimulates AMPK activity. AICAR acts by entering nucleoside pools, significantly increasing levels of adenosine during periods of ATP breakdown. Induces the phosphorylation of TORC2 and inhibits its entry into the nucleus. Mimics the effects of insulin on the expression of two gluconeogenic genes PEPCK and glucose-6-phosphatase. Inhibits PPAR α coactivation and adipocyte differentiation. Substrate for the AICAR transformylase activity of ATIC. Inhibits the growth of prostate cancer cells by inhibiting fatty acid synthesis. Suppresses LPS-induced TNF production via the inhibition of Akt activation. Induces apoptosis in Jurkat cells via a non-AMPK pathway.

Citations: 13

[View Online »](#)

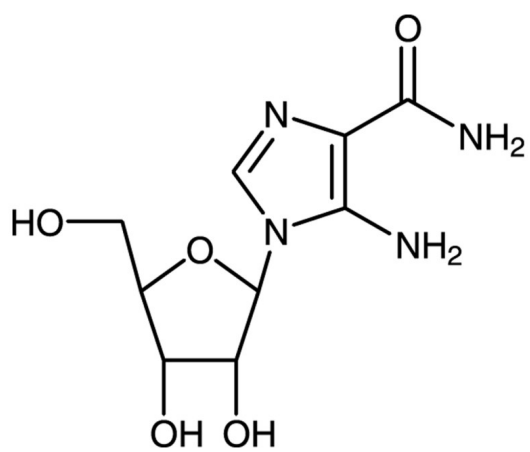
Ordering Information

[Order Online »](#)

BML-EI330-0050	50mg
BML-EI330-0250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

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

Appearance Off-white solid.

CAS 2627-69-2

Couple Target AICAR transformylase, AMPK

Couple Type Activator, Substrate

Formula $C_9H_{14}N_4O_5$

MW 258.2

Purity ≥98% (TLC)

Solubility Soluble in water (9mg/ml).

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)