

BML-275

AMPK inhibitor

Cell-permeable pyrazolopyrimidine derivative that inhibits AMP kinase (K_i =109nM in the absence of AMP) in an ATP-competitive manner. It displays no significant inhibition of ZAPK, SYK, PKCT, PKA and JAK3. Decreases food intake in mice and inhibits the effects of AICAR and metformin. It has been shown to inhibit BMP type I receptors ALK2, ALK3 and ALK6. Promotes cardiomyogenesis in mouse embryonic stem cells. Induces protective autophagy in cancer cell lines.

Citations: 16

[View Online »](#)

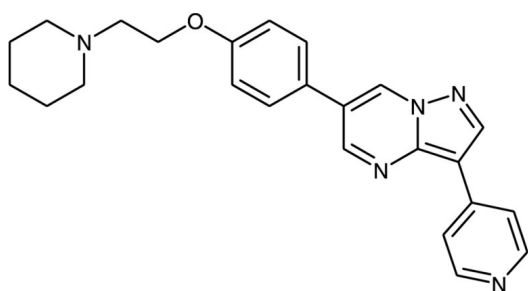
Ordering Information

[Order Online »](#)

BML-EI369-0005	5mg
BML-EI369-0025	25mg

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Compound C, Dorsomorphin

Appearance Yellow solid.

CAS 866405-64-3

Couple Target Activin receptor, AMPK, BMP receptor

Couple Type Inhibitor

Formula $C_{24}H_{25}N_5O$

MW 399.5

Purity ≥98% (TLC)

Solubility Soluble in DMSO (3mg/ml, warm)

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)