

# 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 »](#)

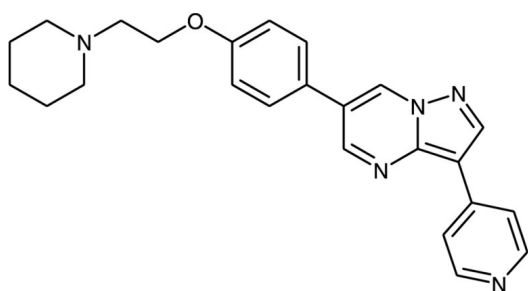
## Ordering Information

[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)



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)