

SD-169

p38 inhibitor

A selective inhibitor of p38 α (IC_{50} = 3.2nM) and p38 β (IC_{50} = 3.2nM), but not p38 γ (IC_{50} > 50 μ M), ERK2, JNK1 or MAPKAP2. Lowers blood glucose and improves glucose homeostasis in nonobese diabetic mice. Enhances axonal regeneration in animals with nerve transection.

Citations: 2

[View Online »](#)

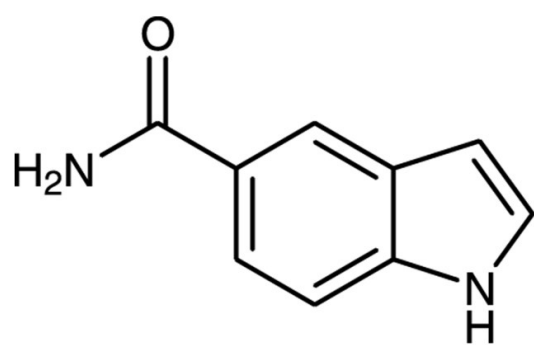
Ordering Information

[Order Online »](#)

BML-EI371-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1H-Indole-5-carboxamide

Appearance White to off-white crystalline powder.

CAS 1670-87-7

Couple Target p38

Couple Type Inhibitor

Formula $C_9H_8N_2O$

Identity Determined by IR.

MW 160.2

Purity $\geq 96.5\%$ (HPLC)

Solubility Soluble in water (1.5mg/ml); DMSO (15mg/ml) or Ethanol (15mg/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-
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)