

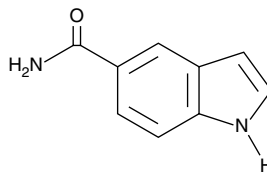
Product Information



SD 169

Item No. 14156

CAS Registry No.: 1670-87-7
Formal Name: 1H-indole-5-carboxamide
Synonym: 5-Carbamoylindole
MF: C₉H₈N₂O
FW: 160.2
Purity: ≥97%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 237, 276 nm



Laboratory Procedures

For long term storage, we suggest that SD 169 be stored as supplied at -20°C. It should be stable for at least two years. SD 169 is supplied as a crystalline solid. A stock solution may be made by dissolving the SD 169 in the solvent of choice. SD 169 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of SD 169 in these solvents is approximately 1.4, 5, and 16 mg/ml, respectively.

SD 169 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SD 169 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SD 169 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

SD 169 is a selective ATP competitive inhibitor of the MAP kinases p38α (IC₅₀ = 3.2 nM) and p38β (IC₅₀ = 122 nM).¹ It has no inhibitory effect against a panel of other kinases, including p38γ MAP kinase, ERK2, JNK-1, and MAPKAPK-2, when tested *in vitro* at 50 μM.¹ SD 169 is orally active, significantly reducing p38 MAP kinase expression in T cells of nonobese diabetic mice when delivered in chow at 600 mg/kg.¹ In this model, SD 169 also reduced the incidence of diabetes, lowered blood glucose, and improved glucose homeostasis.¹ SD 169 also enhances axonal regeneration after sciatic nerve crush injury in rats, when given by gavage (30 mg/kg) before and after nerve damage.²

References

1. Medicherla, S., Protter, A.A., Ma, J.Y., *et al.* Preventive and therapeutic potential of p38α-selective mitogen-activated protein kinase inhibitor in nonobese diabetic mice with type 1 diabetes. *J. Pharmacol. Exp. Ther.* **318(1)**, 99-107 (2006).
2. Myers, R.R., Sekiguchi, Y., Kikuchi, S., *et al.* Inhibition of p38 MAP kinase activity enhances axonal regeneration. *Exp. Neurol.* **184(2)**, 606-614 (2003).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/14156

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 02/15/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com