

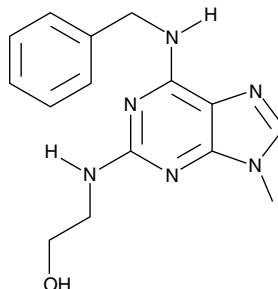
Product Information



Olomoucine

Item No. 10010240

CAS Registry No.: 101622-51-9
Formal Name: 2-[[9-methyl-6-[(phenylmethyl)amino]-9H-purin-2-yl]amino]-ethanol
MF: C₁₅H₁₈N₆O
FW: 298.3
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 231, 290 nm



Laboratory Procedures

For long term storage, we suggest that olomoucine be stored as supplied at -20°C. It should be stable for at least two years.

Olomoucine is supplied as a crystalline solid. A stock solution may be made by dissolving the olomoucine in an organic solvent purged with an inert gas. Olomoucine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of olomoucine in these solvents is approximately 10, 20, and 30 mg/ml, respectively.

Olomoucine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, olomoucine should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Olomoucine 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.

Cyclin-dependent kinases (CDKs) are key regulators of cell cycle progression and are therefore promising targets for cancer therapy. Olomoucine is a CDK inhibitor that acts by competing for the ATP binding site of the kinase. This purine derivative is an inhibitor for CDC2/cyclin B (IC₅₀ = 7 μM), Cdk2/cyclin A (IC₅₀ = 7 μM), Cdk2/cyclin E (IC₅₀ = 7 μM), CDK/p35 kinase (IC₅₀ = 3 μM), and ERK1/p44 MAP kinase (IC₅₀ = 25 μM).¹ Olomoucine inhibits DNA synthesis in interleukin-2 stimulated T lymphocytes, triggers G₁ arrest, and is also used to synchronize cells in G₁.¹ This inhibitor can attenuate astroglial proliferation and glial scar formation, decrease lesion cavity and mitigate functional deficits after spinal cord injury (SCI) in rats.² Administration of olomoucine (3 mg/kg/day) to rats after SCI significantly suppressed microglial proliferation and proinflammatory cytokine expression, reduced tissue edema formation, and attenuated the number of apoptotic neurons.²

References

1. Abraham, R.T., Acquarone, M., Andersen, A., *et al.* Cellular effects of olomoucine, an inhibitor of cyclin-dependent kinases. *Biology of the Cell* **83**(2), 105-120 (1995).
2. Tian, D.-S., Xie, M.-J., Yu, Z.-Y., *et al.* Cell cycle inhibition attenuates microglia induced inflammatory response and alleviates neuronal cell death after spinal cord injury in rats. *Brain Res.* **1135**, 177-185 (2007).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10010240

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

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 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, 10/11/2013

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