

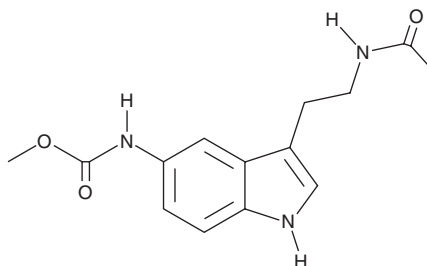
PRODUCT INFORMATION



5-Methoxycarbonylamino-N-acetyltryptamine

Item No. 39553

CAS Registry No.: 190277-13-5
Formal Name: N-[3-[2-(acetylamino)ethyl]-1H-indol-5-yl]-carbamic acid, methyl ester
Synonyms: GR135531, 5-MCA-NAT, MCA-NAT
MF: C₁₄H₁₇N₃O₃
FW: 275.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

5-Methoxycarbonylamino-N-acetyltryptamine (5-MCA-NAT) is supplied as a solid. A stock solution may be made by dissolving the 5-MCA-NAT in the solvent of choice, which should be purged with an inert gas. 5-MCA-NAT is slightly soluble in methanol.

Description

5-MCA-NAT is a melatonin (Item No. 14427) derivative and a partial agonist of melatonin receptor 1 (MT₁) and MT₂.¹ It inhibits forskolin-induced production of cAMP in CHO cells expressing human MT₁ or MT₂ (EC₅₀s = 440 and 98 nM, respectively). *In vivo*, ocular administration of 5-MCA-NAT reduces intraocular pressure in control and DBA/2J glaucomatous mice.²

References

1. Vincent, L., Cohen, W., Delagrang, P., *et al.* Molecular and cellular pharmacological properties of 5-methoxycarbonylamino-N-acetyltryptamine (MCA-NAT): A nonspecific MT₃ ligand. *J. Pineal Res.* **48(3)**, 222-229 (2010).
2. Martínez-Águila, A., Fonseca, B., Pérez de Lara, M.J., *et al.* Effect of melatonin and 5-methoxycarbonylamino-N-acetyltryptamine on the intraocular pressure of normal and glaucomatous mice. *J. Pharmacol. Exp. Ther.* **357(2)**, 293-299 (2016).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/11/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM