

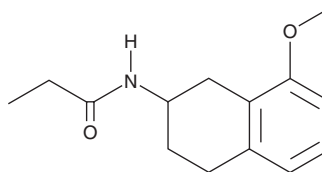
PRODUCT INFORMATION



8-M-PDOT

Item No. 29521

CAS Registry No.: 134865-70-6
Formal Name: N-(1,2,3,4-tetrahydro-8-methoxy-2-naphthalenyl)-propanamide
Synonym: AH 002
MF: C₁₄H₁₉NO₂
FW: 233.3
Purity: ≥98%
Supplied as: A crystalline 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

8-M-PDOT is supplied as a crystalline solid. A stock solution may be made by dissolving the 8-M-PDOT in the solvent of choice, which should be purged with an inert gas. 8-M-PDOT is soluble in the organic solvent DMSO.

Description

8-M-PDOT is a melatonin receptor 2 (MT₂) agonist (K_i = 1.12 nM).¹ It is 5.2-fold selective for MT₂ over MT₁ receptors in a radioligand binding assay. Striatal microinfusion of 8-M-PDOT (10 µg/µl) increases the percentage of time spent in the open arms of the elevated plus maze, indicating anxiolytic-like activity, in a rat model of rotenone-induced Parkinson's disease.²

References

1. Browning, C., Beresford, I., Fraser, N., *et al.* Pharmacological characterization of human recombinant melatonin mt₁ and MT₂ receptors. *Br. J. Pharmacol.* **129**(5), 877-886 (2000).
2. Nosedá, A.C., Targa, A.D., Rodrigues, L.S., *et al.* REM sleep deprivation promotes a dopaminergic influence in the striatal MT2 anxiolytic-like effects. *Sleep Sci.* **9**(1), 47-54 (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, 09/29/2022

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