

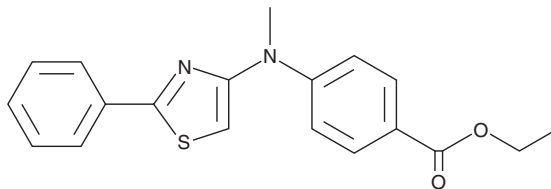
PRODUCT INFORMATION



Neuropathiazol

Item No. 43953

CAS Registry No.: 880090-88-0
Formal Name: 4-[methyl(2-phenyl-4-thiazolyl)amino]-benzoic acid, ethyl ester
MF: C₁₉H₁₈N₂O₂S
FW: 338.4
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

Neuropathiazol is supplied as a solid. A stock solution may be made by dissolving the neuropathiazol in the solvent of choice, which should be purged with an inert gas. Neuropathiazol is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Neuropathiazol is an inducer of neuronal differentiation.¹ It induces differentiation of primary rat hippocampal neural progenitor cells when used at a concentration of 10 μM. Neuropathiazol (20 mg/kg per day), alone and in combination with cyclophosphamide (Item No. 13849), decreases tumor volume in an SK-N-AS neuroblastoma mouse xenograft model.² It increases intratumoral levels of neuronatin, microtubule-associated protein 2 (MAP2), and neurofilament light polypeptide (NF-L), markers of differentiation, in the same model.

References

1. Warashina, M., Min, K.H., Kuwabara, T., *et al.* A synthetic small molecule that induces neuronal differentiation of adult hippocampal neural progenitor cells. *Angew. Chem. Int. Ed. Engl.* **45(4)**, 591-593 (2006).
2. Xu, H., Zhang, F., Xu, Y., *et al.* Neuropathiazol induces neuronal-like differentiation in neuroblastoma cells via upregulation of PEG5. *Pediatr. Res.* (2025).

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, 09/03/2025

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