

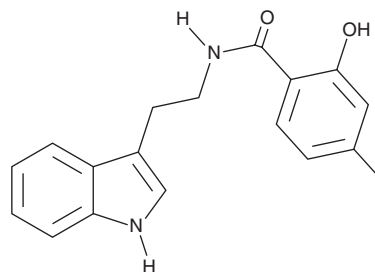
PRODUCT INFORMATION



p-methyl-N-salicyloyl Tryptamine

Item No. 38958

CAS Registry No.: 1371564-45-2
Formal Name: 2-hydroxy-N-[2-(1H-indol-3-yl)ethyl]-4-methyl-benzamide
Synonyms: LZWL02003, *para*-methyl-N-salicyloyl Tryptamine
MF: C₁₈H₁₈N₂O₂
FW: 294.4
Purity: ≥98%
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

p-methyl-N-salicyloyl Tryptamine is supplied as a solid. A stock solution may be made by dissolving the *p*-methyl-N-salicyloyl tryptamine in the solvent of choice, which should be purged with an inert gas. *p*-methyl-N-salicyloyl Tryptamine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of *p*-methyl-N-salicyloyl tryptamine in these solvents is approximately 10 mg/ml.

Description

p-methyl-N-salicyloyl Tryptamine is a neuroprotective agent and derivative of N-salicyloyl tryptamine.¹ It reduces LPS-induced nitric oxide (NO) production in C6 rat glioma cells and BV-2 mouse microglia (IC₅₀s = 7.48 and 4.94 μM, respectively). It also reduces LPS-induced production of TNF-α and prostaglandin E₂ (PGE₂; Item No. 14010) and increases production of IL-10 in BV-2 microglia when used at a concentration of 40 μM. *In vivo*, *p*-methyl-N-salicyloyl tryptamine (25-100 mg/kg) decreases cortical and hippocampal amyloid-β (Aβ) plaque formation and neuronal pyroptosis, as well as improves spatial learning and memory, in a mouse model of Aβ-induced Alzheimer's disease.²

References

1. Fan, X., Li, J., Deng, X., *et al.* Design, synthesis and bioactivity study of N-salicyloyl tryptamine derivatives as multifunctional agents for the treatment of neuroinflammation. *Eur. J. Med. Chem.* **193**, 112217 (2020).
2. Bai, Y., Liu, D., Zhang, H., *et al.* N-salicyloyl tryptamine derivatives as potential therapeutic agents for Alzheimer's disease with neuroprotective effects. *Bioorg. Chem.* **115**, 105255 (2021).

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, 11/26/2025

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