

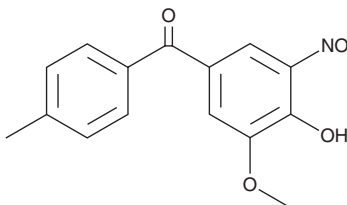
PRODUCT INFORMATION



3-O-methyl Tolcapone

Item No. 45833

CAS Registry No.: 134612-80-9
Formal Name: (4-hydroxy-3-methoxy-5-nitrophenyl)
(4-methylphenyl)-methanone
Synonyms: 3-OMT, Ro 40-7591
MF: C₁₅H₁₃NO₅
FW: 287.3
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

3-O-methyl Tolcapone is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-O-methyl tolcapone in the solvent of choice, which should be purged with an inert gas. 3-O-methyl Tolcapone is sparingly soluble (1-10 mg/ml) in DMSO.

Description

3-O-methyl Tolcapone is a stabilizer of transthyretin (TTR) and a metabolite of tolcapone (Item No. 20768).¹ Unlike tolcapone, 3-O-methyl tolcapone does not inhibit catechol-O-methyltransferase (COMT) and is not cytotoxic to SH-SY5Y or HepG2 cells (EC₅₀s = 226.3 and 262.3 μM, respectively).^{1,2} 3-O-methyl Tolcapone (1 and 10 μM) stabilizes TTR in isolated human plasma and binds to TTR in a cell-free assay (K_d = 33 nM).²

References

1. Heizmann, P., Schmitt, M., Leube, J., *et al.* Determination of the catechol-O-methyltransferase inhibitor tolcapone and three of its metabolites in animal and human plasma and urine by reversed-phase high-performance liquid chromatography with ultraviolet detection. *J. Chromatogr. B Biomed. Sci. Appl.* **730(2)**, 153-160 (1999).
2. Poonsiri, T., Dell'Accantera, D., Loconte, V., *et al.* 3-O-methyltolcapone and its lipophilic analogues are potent inhibitors of transthyretin amyloidogenesis with high permeability and low toxicity. *Int. J. Mol. Sci.* **25(1)**, 479 (2023).

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, 06/16/2026

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