

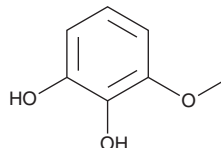
PRODUCT INFORMATION



3-Methoxycatechol

Item No. 36535

CAS Registry No.: 934-00-9
Formal Name: 3-methoxy-1,2-benzenediol
Synonyms: 1,2-Dihydroxy-3-methoxybenzene,
3-Methoxypyrocatechol, NSC 66525,
Pyrogallol 1-monomethyl ether
MF: C₇H₈O₃
FW: 140.1
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

3-Methoxycatechol is supplied as a solid. A stock solution may be made by dissolving the 3-methoxycatechol in the solvent of choice, which should be purged with an inert gas. 3-Methoxycatechol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-methoxycatechol in DMF is approximately 3 mg/ml and approximately 2 mg/ml in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 3-methoxycatechol can be prepared by directly dissolving the 3-methoxycatechol in aqueous buffers. The solubility of 3-methoxycatechol in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

3-Methoxycatechol is a lignan-derived phenol.¹ It is an agonist of G protein-coupled receptor 35 (GPR35) with an EC₅₀ value of 147 μM in a dynamic mass redistribution (DMR) assay using HT-29 cells.² Dietary administration of 3-methoxycatechol (2%) alone, or in a model of multiorgan carcinogenesis induced by nitrosamines, promotes esophageal carcinogenesis in rats.³ 3-Methoxycatechol has been used as a precursor in the enzymatic synthesis of the phenols pyrogallol (Item No. 20347) and purpurogallin (Item No. 29689) and in the electro-organic synthesis of coumestan derivatives.^{1,4}

References

1. Zhang, S., Xiaofeng, W., and Xiao, Y. Conversion of lignin-derived 3-methoxycatechol to the natural product purpurogallin using bacterial P450 GcoAB and laccase CueO. *Appl. Microbiol. Biotechnol.* **106**(2), 593-603 (2022).
2. Deng, H. and Fang, Y. The three catechol benserazide, catechol and pyrogallol are GPR35 agonists. *Pharmaceuticals (Basel)* **6**(4), 500-509 (2013).
3. Hirose, M., Tnaka, H., Takahashi, S., et al. Effects of sodium nitrite and catechol, 3-methoxycatechol, or butylated hydroxyanisole in combination in a rat multiorgan carcinogenesis model. *Cancer Res.* **53**(1), 32-37 (1993).
4. Golabi, S.M. and Nematollahi, D. Electrochemical study of catechol and some 3-substituted catechols in the presence of 4-hydroxy coumarin: Application to the electro-organic synthesis of new coumestan derivatives. *J. Electroanal. Chem.* **420**(1-2), 127-134 (1997).

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, 12/20/2022

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