

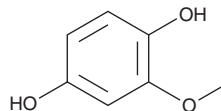
PRODUCT INFORMATION



2-Methoxyhydroquinone

Item No. 36537

CAS Registry No.: 824-46-4
Formal Name: 2-methoxy-1,4-benzenediol
Synonyms: o-Methoxyhydroquinone, MHQ, MOHQ
MF: C₇H₈O₃
FW: 140.1
Purity: ≥90%
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

2-Methoxyhydroquinone is supplied as a solid. A stock solution may be made by dissolving the 2-methoxyhydroquinone in the solvent of choice, which should be purged with an inert gas. 2-Methoxyhydroquinone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2-methoxyhydroquinone in these solvents is approximately 10, 5, and 2 mg/ml, respectively.

Description

2-Methoxyhydroquinone is a phenolic compound. It reduces TNF- α -induced production of chemokine (C-C motif) ligand 2 (CCL2) *in vitro* with an IC₅₀ value of 64.3 μ M.¹ 2-Methoxyhydroquinone has been used as a precursor in the synthesis of the Hsp90 inhibitor geldanamycin (Item No. 13355).²

References

1. Houser, K.R., Johnson, D.K., and Ishmael, F.T. Anti-inflammatory effects of methoxyphenolic compounds on human airway cells. *J. Inflamm. (Lond.)* **9**, 6 (2012).
2. Qin, H.-L. and Panek, J.S. Total synthesis of the Hsp90 inhibitor geldanamycin. *Org. Lett.* **10(12)**, 2477-2479 (2008).

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