

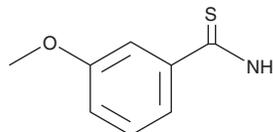
PRODUCT INFORMATION



3-methoxythio Benzamide

Item No. 10553

CAS Registry No.: 64559-06-4
Formal Name: 3-methoxy-benzenecarbothioamide
MF: C₈H₉NOS
FW: 167.2
Purity: ≥98%
UV/Vis.: λ_{max}: 253, 299 nm
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-methoxythio Benzamide is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-methoxythio benzamide in the solvent of choice, which should be purged with an inert gas. 3-methoxythio Benzamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-methoxythio benzamide in ethanol is approximately 3 mg/ml and approximately 30 mg/ml in DMSO and DMF.

3-methoxythio Benzamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3-methoxythio benzamide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 3-methoxythio Benzamide has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

3-methoxythio Benzamide is a synthetic intermediate useful for pharmaceutical synthesis. Thiobenzamides, including 3-methoxythio benzamide, are potent hepatotoxins *in vivo*.¹

Reference

1. Ikehata, K., Duzhak, T.G., Galeva, N.A., *et al.* Protein targets of reactive metabolites of thiobenzamide in rat liver *in vivo*. *Chem. Res. Toxicol.* **21**(7), 1432-1442 (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, 11/07/2022

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