

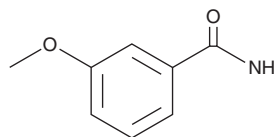
PRODUCT INFORMATION



m-Methoxybenzamide

Item No. 30451

CAS Registry No.: 5813-86-5
Formal Name: 3-methoxy-benzamide
Synonyms: *m*-Anisamide, *meta*-Methoxybenzamide, NSC 28589, NSC 209527
MF: C₈H₉NO₂
FW: 151.2
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

m-Methoxybenzamide is supplied as a crystalline solid. A stock solution may be made by dissolving the *m*-methoxybenzamide in the solvent of choice, which should be purged with an inert gas. *m*-Methoxybenzamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of *m*-methoxybenzamide in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

m-Methoxybenzamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, *m*-methoxybenzamide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. *m*-Methoxybenzamide has a solubility of approximately 0.50 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

m-Methoxybenzamide is an inhibitor of poly(ADP-ribose) polymerase (PARP; K_i = <2 μM).¹ It also inhibits human placental ADP ribosyl transferase activity by 93% when used at a concentration of 100 μM.² *m*-Methoxybenzamide (100, 200, and 300 mg/kg) increases the pain threshold in the paw pinch test in rats.³

References

1. Purnell, M.R. and Whish, W.J. Novel inhibitors of poly(ADP-ribose) synthetase. *Biochem. J.* **185**(3), 775-777 (1980).
2. Burtcher, H.J., Auer, B., Klocker, H., *et al.* Isolation of ADP-ribosyltransferase by affinity chromatography. *Anal. Biochem.* **152**(2), 285-290 (1986).
3. Sanders, R., Dill, R.E., Dorris, R.L., *et al.* Analgesic effects of 3-methoxybenzamide in rats. *J. Pharm. Pharmacol.* **41**(4), 268-269 (1989).

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, 11/02/2022

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