

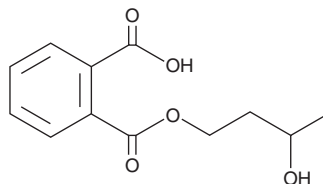
PRODUCT INFORMATION



MHBP

Item No. 46077

CAS Registry No.: 57074-43-8
Formal Name: 1,2-benzenedicarboxylic acid,
1-(3-hydroxybutyl) ester
Synonyms: 3-hydroxy MBP,
Mono-3-hydroxybutyl Phthalate,
Mono-3-hydroxy-*n*-butyl Phthalate
MF: C₁₂H₁₄O₅
FW: 238.2
Purity: ≥95%
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

MHBP is supplied as a solid. A stock solution may be made by dissolving the MHBP in the solvent of choice, which should be purged with an inert gas. MHBP is slightly soluble in methanol and DMSO.

Description

MHBP is a metabolite of the plasticizer and environmental contaminant di-*n*-butyl phthalate (DBP).¹ Unlike other phthalates, MHBP is not associated with negative reproductive health outcomes in patients undergoing assisted reproductive treatment.²

References

1. Silva, M.J., Samandar, E., Reidy, J.A., *et al.* Metabolite profiles of di-*n*-butyl phthalate in humans and rats. *Environ. Sci. Technol.* **41**, 7576-7580 (2007).
2. Xie, G., Zhou, Y., Zheng, Q., *et al.* Association between phthalate exposure and reproductive health in patients undergoing assisted reproductive treatment: A systematic review and meta-analysis. *Reproductive Toxicology* **136**, 108948 (2025).

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, 06/29/2026

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