

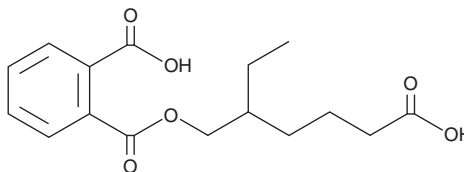
PRODUCT INFORMATION



MECPP

Item No. 46076

CAS Registry No.: 40809-41-4
Formal Name: 1,2-benzenedicarboxylic acid,
1-(5-carboxy-2-ethylpentyl) ester
Synonyms: 5-cx MEHP, 5-carboxy MEPP,
Mono-(2-ethyl-5-carboxypentyl) Phthalate
MF: C₁₆H₂₀O₆
FW: 308.3
Purity: ≥95%
Supplied as: A semi-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

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

Description

MECPP is an inactive urinary metabolite of the plasticizer and environmental contaminant di-(2-ethylhexyl) phthalate (DEHP) and *rac*-mono(ethylhexyl) phthalate (MEHP; Item No. 35605).^{1,2} It is associated with lower systolic blood pressure in female adolescents and lower serum LDL-cholesterol (LDL-C) in male adolescents.³ MECPP is also associated with lower mental development indices in female infants.⁴

References

1. Chiang, H.-c., Kuo, Y.-T., Shen, C.-C., *et al.* Mono(2-ethylhexyl)phthalate accumulation disturbs energy metabolism of fat cells. *Arch. Toxicol.* **90**(3), 589-601 (2016).
2. Kambia, N., Séverin, I., Farce, A., *et al.* Comparative effects of di-(2-ethylhexyl)phthalate and di-(2-ethylhexyl)terephthalate metabolites on thyroid receptors: In vitro and in silico studies. *Metabolites* **11**(2), 94 (2021).
3. Berghuis, S.A., Bocca, G., Bos, A.F., *et al.* Adolescent urinary concentrations of phthalate metabolites and indices of overweight and cardiovascular risk in Dutch adolescents. *Environ. Int.* **194**, 109167 (2024).
4. Téllez-Rojo, M.M., Cantoral, A., Cantonwine, D.E., *et al.* Prenatal urinary phthalate metabolites levels and neurodevelopment in children at two and three years of age. *Sci. Total Environ.* **461-462**, 386-390 (2013).

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, 06/30/2026

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