

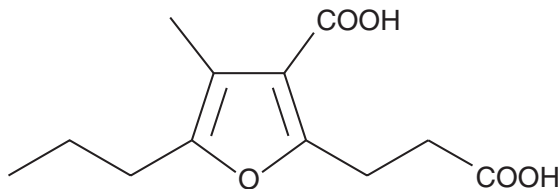
PRODUCT INFORMATION



CMPF

Item No. 10007133

CAS Registry No.: 86879-39-2
Formal Name: 3-carboxy-4-methyl-5-propyl-2-furanpropanoic acid
Synonym: FA 12:4;O3
MF: C₁₂H₁₆O₅
FW: 240.3
Purity: ≥98%
UV/Vis.: λ_{max}: 260 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

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

CMPF is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CMPF should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CMPF has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Furan fatty acids are unique, naturally occurring lipids that are found in significant amounts in dietary phospholipids, such as in salmon roe.¹ CMPF is an endogenous metabolite of furan fatty acids in humans. CMPF is highly albumin-bound and accumulates in the serum of uremic patients to concentrations in excess of 0.2 mM. Its primary effect is to inhibit cellular transport and subsequent deiodination of thyroxine (T₄).^{2,3} CMPF is tightly bound to albumin but only moderately inhibits T₄ binding in a direct manner (10-14% at 0.3 mM). However, CMPF effectively displaces competitive T₄ binding molecules from albumin, such as acidic drugs and free fatty acids.³ Therefore, CMPF may indirectly influence T₄ binding to albumin by increasing the serum concentration of competitive binding molecules, particularly free fatty acids such as oleic acid.³

References

1. Ishii, K., Okajima, H., Okada, Y., *et al.* Studies on furan fatty acids of salmon roe phospholipids. *J. Biochem.* **103**(5), 836-839 (1988).
2. Lim, C.-F., Bernard, B.F., De Jong, M., *et al.* A furan fatty acid and indoxyl sulfate are the putative inhibitors of thyroxine hepatocyte transport in uremia. *J. Clin. Endocrinol. Metab.* **76**(2), 318-324 (1993).
3. Lim, C.-F., Stockigt, J.R., Curtis, A.J., *et al.* A naturally occurring furan fatty acid enhances drug inhibition of thyroxine binding in serum. *Metabolism* **42**(11), 1468-1474 (1993).

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, 03/11/2024

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