

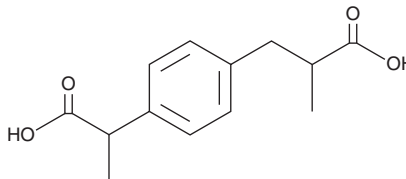
PRODUCT INFORMATION



Ibuprofen Carboxylic Acid

Item No. 26594

CAS Registry No.: 15935-54-3
Formal Name: 4-(2-carboxypropyl)- α -methylbenzeneacetic acid
Synonym: Carboxyibuprofen
MF: C₁₃H₁₆O₄
FW: 236.3
Purity: \geq 98% (mixture of diastereomers)
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ibuprofen carboxylic acid is supplied as a solid. A stock solution may be made by dissolving the ibuprofen carboxylic acid in the solvent of choice. Ibuprofen carboxylic acid is soluble in organic solvents such as chloroform and ethyl acetate, which should be purged with an inert gas.

Description

Ibuprofen carboxylic acid is a major metabolite of ibuprofen (Item Nos. 70280 | 16793 | 16794).¹ It prevents fructose-, cyanate-, and prednisolone-induced inactivation of catalase and fructose- and cyanate-induced inactivation of fumarase when used at a concentration of 2 mM and penetrates into the lens of isolated bovine eyes.² Ibuprofen carboxylic acid has been found in bank filtrate and surface water, as well as influent and effluent wastewater, and is considered a micropollutant.³

References

1. Mills, R.F., Adams, S.S., Cliffe, E.E., *et al.* The metabolism of ibuprofen. *Xenobiotica* **3(9)**, 589-598 (1973).
2. Shyadehi, A.Z. and Harding, J.J. Protective effects of ibuprofen and its major metabolites against in vitro inactivation of catalase and fumarase: Relevance to cataract. *Biochim. Biophys. Acta.* **1587(1)**, 31-35 (2002).
3. Hermes, N., Jewell, K.S., Wick, A., *et al.* Quantification of more than 150 micropollutants including transformation products in aqueous samples by liquid chromatography-tandem mass spectrometry using scheduled multiple reaction monitoring. *J. Chromatogr. A.* **1531**, 64-73 (2018).

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, 10/27/2022

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