

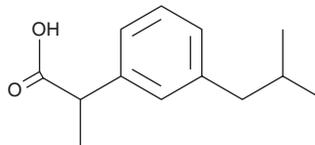
PRODUCT INFORMATION



m-isobutyl Ibuprofen

Item No. 40127

CAS Registry No.: 66622-47-7
Formal Name: α -methyl-3-(2-methylpropyl)-benzeneacetic acid
Synonyms: Ibuprofen Impurity A, *meta*-isobutyl Ibuprofen
MF: C₁₃H₁₈O₂
FW: 206.3
Purity: \geq 98%
Supplied as: An oil
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

m-isobutyl Ibuprofen is supplied as an oil. A stock solution may be made by dissolving the *m*-isobutyl ibuprofen in the solvent of choice, which should be purged with an inert gas. *m*-isobutyl Ibuprofen is soluble in acetonitrile and DMSO.

Description

m-isobutyl Ibuprofen is a potential impurity in commercial preparations of the non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor ibuprofen.¹

Reference

1. Vijay Kumar, R., Rao, V.U., Anil Kumar, N., *et al.* A novel, rapid, and validated stability-indicating UPLC method for the estimation of drotaverine hydrochloride and ibuprofen impurities in oral solid dosage form. *Sci. Pharm.* **83**(4), 567-581 (2015).

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, 02/14/2024

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