

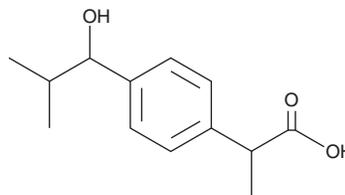
PRODUCT INFORMATION



1-hydroxy Ibuprofen

Item No. 36295

CAS Registry No.: 53949-53-4
Formal Name: 4-(1-hydroxy-2-methylpropyl)- α -methylbenzeneacetic acid
MF: C₁₃H₁₈O₃
FW: 222.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 219 nm
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

1-hydroxy Ibuprofen is supplied as a solid. A stock solution may be made by dissolving the 1-hydroxy ibuprofen in the solvent of choice, which should be purged with an inert gas. 1-hydroxy Ibuprofen is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1-hydroxy ibuprofen in these solvents is approximately 10 mg/ml.

Description

1-hydroxy Ibuprofen is a metabolite of the non-steroidal anti-inflammatory drug (NSAID) and non-selective COX inhibitor ibuprofen (Item Nos. 70280 | 16793 | 16794).¹

Reference

1. Neunzig, I., Göhring, A., Drăgan, C.-A., *et al.* Production and NMR analysis of the human ibuprofen metabolite 3-hydroxyibuprofen. *J. Biotechnol.* **157**(3), 417-420 (2012).

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, 11/14/2022

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