

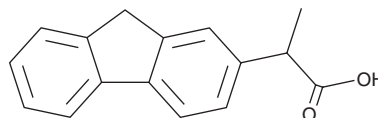
PRODUCT INFORMATION



Cicloprofen

Item No. 33571

CAS Registry No.: 36950-96-6
Formal Name: α -methyl-9H-fluorene-2-acetic acid
Synonyms: (\pm)-Cicloprofen, α -Methylfluorene-2-acetic Acid, NSC 293916, SQ 20,824
MF: C₁₆H₁₄O₂
FW: 238.3
Purity: \geq 95%
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

Cicloprofen is supplied as a solid. A stock solution may be made by dissolving the cicloprofen in the solvent of choice, which should be purged with an inert gas. Cicloprofen is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of cicloprofen in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Cicloprofen is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cicloprofen should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cicloprofen has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cicloprofen is a non-steroidal anti-inflammatory drug (NSAID).^{1,2} It inhibits histamine release induced by ragweed, A23187 (Item No. 11016), or concanavalin A (Item No. 14951) in isolated human leukocytes when used at a concentration of 0.5 mM.¹ Cicloprofen inhibits carrageenan-induced paw edema in rats (ED₅₀ = 67 mg/kg).²

References

- Hearn, T. and Wojnar, R.J. Characterization of augmentation of allergic and non-allergic histamine release by non-steroidal anti-inflammatory/analgesic agents. *Agents Actions* **11(4)**, 352-360 (1981).
- Stiller, E.T., Diassi, P.A., Gerschutz, D., *et al.* Synthesis and antiinflammatory activities of α -methylfluorene-2-acetic acid and related compounds. *J. Med. Chem.* **15(10)**, 1029-1032 (1972).

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, 11/15/2022

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