

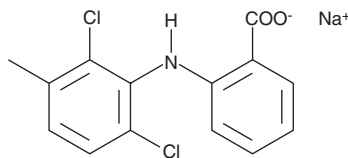
PRODUCT INFORMATION



Meclofenamate (sodium salt)

Item No. 70550

CAS Registry No.: 6385-02-0
Formal Name: 2-[(2,6-dichloro-3-methylphenyl)amino]-
benzoic acid, monosodium salt
Synonym: Meclofenamic Acid
MF: C₁₄H₁₀Cl₂NO₂ • Na
FW: 318.1
Purity: ≥98%
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

Meclofenamate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the meclofenamate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Meclofenamate (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of meclofenamate (sodium salt) in these solvents is approximately 100, 75, and 39 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of meclofenamate (sodium salt) can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of meclofenamate (sodium salt) in PBS (pH 7.2) and PBS (pH 9.0) is approximately 0.05 and 50 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Meclofenamate is a time-dependent, non-selective competitive inhibitor of COX-1 and -2. The IC₅₀ values for inhibition of human recombinant COX-1 and -2 are 1.5 and 9.7 μM, respectively for instantaneous inhibition. However, the IC₅₀ is much lower if pre-incubated with the enzyme.¹

Reference

1. Laneuville, O., Breuer, D.K., DeWitt, D.L., *et al.* Differential inhibition of human prostaglandin endoperoxide H synthases-1 and -2 by nonsteroidal anti-inflammatory drugs. *J. Pharmacol. Exp. Ther.* **271(2)**, 927-934 (1994).

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, 11/10/2022

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