

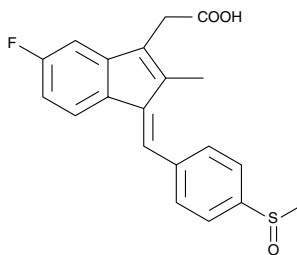
Product Information



Sulindac

Item No. 10004386

CAS Registry No.: 38194-50-2
Formal Name: 5-fluoro-2-methyl-1Z-[[4-(methylsulfinyl)phenyl]methylene]-1H-indene-3-acetic acid
MF: C₂₀H₁₇FO₃S
FW: 356.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 226, 258, 286, 329 nm



Laboratory Procedures

For long term storage, we suggest that sulindac be stored as supplied at -20°C. It should be stable for at least two years.

Sulindac is supplied as a crystalline solid. A stock solution may be made by dissolving the sulindac in an organic solvent purged with an inert gas. Sulindac is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of sulindac in these solvents is approximately 2 mg/ml in ethanol and 30 mg/ml in DMSO and DMF.

Sulindac is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sulindac should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Sulindac has a solubility of approximately 0.5 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Many non-steroidal anti-inflammatory drugs (NSAIDs) are potent but non-selective inhibitors of both COX-1 and COX-2 in humans.¹ Sulindac is one of the older NSAIDs, an isostere of indomethacin developed before the inducible form of cyclooxygenase (COX-2) was discovered.² Although a number of NSAIDs have been found to protect against digestive tract cancers, sulindac has an extensive epidemiology documenting reduced human colorectal cancer. In murine models, sulindac was found not only to inhibit the enzymatic activity of polyp-associated COX-2, but also to downregulate the expression of colonic COX-2 protein to control levels.³

References

1. Barnett, J., Chow, J., Ives, D., *et al.* Purification, characterization and selective inhibition of human prostaglandin G/H synthase 1 and 2 expressed in the baculovirus system. *Biochim. Biophys. Acta* **1209**, 130-139 (1994).
2. Shen, T.-Y. and Winter, C.A. Chemical and biological studies on indomethacin, sulindac and their analogs. *Adv. Drug Res.* **12**, 89-98 (1977).
3. Boolbol, S.K., Dannenberg, A.J., Chadburn, A., *et al.* Cyclooxygenase-2 overexpression and tumor formation are blocked by sulindac in a murine model of familial adenomatous polyposis. *Cancer Res.* **56**, 2556-2560 (1996).

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 06/20/2011