

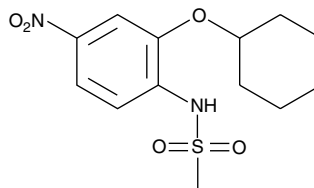
Product Information



NS-398

Item No. 70590

CAS Registry No.: 123653-11-2
Formal Name: N-[2-(cyclohexyloxy)-4-nitrophenyl]-methanesulfonamide
MF: C₁₃H₁₈N₂O₅S
FW: 314.4
Purity: ≥98%
Stability: ≥2 years at room temperature
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 238, 296, 337 nm



Laboratory Procedures

For long term storage, we suggest that NS-398 be stored as supplied at room temperature. It will be stable for at least two years.

NS-398 is supplied as a crystalline solid. A stock solution may be made by dissolving the NS-398 in an organic solvent purged with an inert gas. NS-398 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NS-398 in these solvents is approximately 0.5, 25, and 30 mg/ml, respectively.

NS-398 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NS-398 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NS-398 has a solubility of approximately 400 µg/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.

NS-398 is a selective inhibitor of cyclooxygenase-2 (COX-2). The IC₅₀ values for human recombinant COX-1 and -2 are 75 and 1.77 µM, respectively.¹ The IC₅₀ values for ovine COX-1 and -2 are 220 and 0.15 µM, respectively.²

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. Johnson, J.L., Wimsatt, J., Buckel, S.D., *et al.* Purification and characterization of prostaglandin H synthase-2 from sheep placental cotyledons. *Arch. Biochem. Biophys.* **324**, 26-34 (1995).

Related Products

CAY10404 - Item No. 70210 • Indomethacin heptyl ester - Item No. 70271 • N-(2-phenylethyl)-Indomethacin amide - Item No. 70272 • N-(3-pyridyl)-Indomethacin amide - Item No. 70274 • N-(4-acetamidophenyl)-Indomethacin amide - Item No. 70278 • APHS - Item No. 70330 • Nimesulide - Item No. 70640 • DuP-697 - Item No. 70645 • Niflumic Acid - Item No. 70650 • SC-58125 - Item No. 70655 • COX (ovine) Inhibitor Screening Assay - Item No. 560101 • COX Inhibitor Screening Assay - Item No. 560131 • Chemiluminescent COX (ovine) Inhibitor Screening Assay - Item No. 760101 • Colorimetric COX (ovine) Inhibitor Screening Assay - Item No. 760111

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, 05/02/2011