

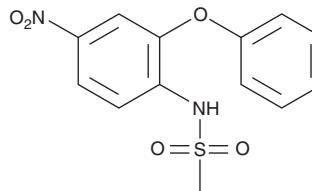
PRODUCT INFORMATION



Nimesulide

Item No. 70640

CAS Registry No.: 51803-78-2
Formal Name: N-(4-nitro-2-phenoxyphenyl)-methanesulfonamide
Synonym: R805
MF: C₁₃H₁₂N₂O₅S
FW: 308.3
Purity: ≥99%
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

Nimesulide is supplied as a crystalline solid. A stock solution may be made by dissolving the nimesulide in the solvent of choice, which should be purged with an inert gas. Nimesulide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of nimesulide in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Nimesulide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nimesulide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nimesulide has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Nimesulide is a non-steroidal anti-inflammatory drug (NSAID) and COX-2 inhibitor (IC₅₀s = 1.27 and 0.03 μM for the human and ovine enzymes, respectively).^{1,2} It is selective for COX-2 over COX-1 (IC₅₀s = 70 and 22 μM for the human and ovine enzymes, respectively). Nimesulide also inhibits sodium-dependent neutral amino acid transporter (B⁰AT1) with an IC₅₀ value of 23 μM for the rat kidney transporter.³ It inhibits infection-induced increases in brain prostaglandin E₂ (PGE₂; Item No. 14010) levels, as well as reduces pyresis (ED₅₀ = 0.3 mg/kg), in yeast-infected rats.⁴ Nimesulide (2.9 mg/kg) inhibits formalin-induced hindpaw thermal hyperalgesia in rats.⁵

References

1. 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**(1), 26-34 (1995).
2. 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**(1), 130-139 (1994).
3. Pochini, L., Seidita, A., Sensi, C., *et al.* Nimesulide binding site in the B⁰AT1 (SLC6A19) amino acid transporter. Mechanism of inhibition revealed by proteoliposome transport assay and molecular modelling. *Biochem. Pharmacol.* **89**(3), 422-430 (2014).
4. Taniguchi, Y., Yokoyama, K., Inui, K., *et al.* Inhibition of brain cyclooxygenase-2 activity and the antipyretic action of nimesulide. *Eur. J. Pharmacol.* **330**, 221-229 (1997).
5. Bianchi, M. and Brogгинi, M. Anti-hyperalgesic effects of nimesulide: Studies in rats and humans. *Int. J. Clin. Pract. Suppl.* **57**(128), 11-19 (2002).

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, 10/04/2022

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