

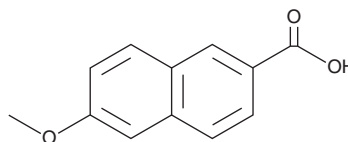
# PRODUCT INFORMATION



## 6-methoxy-2-Naphthoic Acid

Item No. 35374

**CAS Registry No.:** 2471-70-7  
**Formal Name:** 6-methoxy-2-naphthalenecarboxylic acid  
**Synonyms:** 6-methoxy Naphthalene Acetic Acid,  
6-methoxy-2-Naphthylacetic Acid,  
6-MNA, NSC 408786  
**MF:** C<sub>12</sub>H<sub>10</sub>O<sub>3</sub>  
**FW:** 202.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 237, 243 nm  
**Supplied as:** A 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

6-methoxy-2-Naphthoic acid (6-MNA) is supplied as a solid. A stock solution may be made by dissolving the 6-MNA in the solvent of choice, which should be purged with an inert gas. 6-MNA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 6-MNA in these solvents is approximately 3 mg/ml.

### Description

6-MNA is an active metabolite of the non-steroidal anti-inflammatory drug (NSAID) nabumetone (Item No. 20251).<sup>1,2</sup> It inhibits COX-1 and LPS-induced COX-2 activities in isolated human whole blood (IC<sub>50</sub>s = 31.01 and 19.84 μM, respectively).<sup>2</sup> 6-MNA is also a potential impurity in commercial preparations of naproxen.<sup>3</sup>

### References

1. Ray, J.E. and Day, R.O. High-performance liquid chromatographic determination of a new anti-inflammatory agent, nabumetone, and its major metabolite in plasma using fluorimetric detection. *J. Chromatogr.* **336(1)**, 234-238 (1984).
2. Cryer, B. and Feldman, M. Cyclooxygenase-1 and cyclooxygenase-2 selectivity of widely used nonsteroidal anti-inflammatory drugs. *Am. J. Med.* **104(5)**, 413-421 (1998).
3. Hrk, M., Babu, K.S., Lanka, A.R.P., *et al.* Simultaneous estimation of related compounds in esomeprazole and naproxen tablets by using ion pair reverse phase HPLC. *Int. J. Pharm. Pharm. Sci.* **7(2)**, 460-464 (2015).

#### 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, 11/30/2022

#### 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