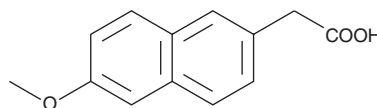


# PRODUCT INFORMATION

## 6-methoxy Naphthalene Acetic Acid

Item No. 70620

**CAS Registry No.:** 23981-47-7  
**Formal Name:** 6-methoxy-2-Naphthaleneacetic acid  
**Synonym:** 6-MNA  
**MF:** C<sub>13</sub>H<sub>12</sub>O<sub>3</sub>  
**FW:** 216.2  
**Purity:** ≥99%  
**UV/Vis.:** λ<sub>max</sub>: 231 nm  
**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

6-methoxy Naphthalene acetic acid (6-MNA) is supplied as a crystalline 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 55, 24, and 25 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 6-MNA can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 6-MNA in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

6-MNA is a competitive, non-selective COX inhibitor.<sup>1-3</sup> The K<sub>i</sub> values for ovine COX-1 and -2 are 21 and 19 μM, respectively.<sup>2</sup> The 50% B/B<sub>0</sub> values are 70 and 20 μM for human recombinant COX-1 and -2, respectively.<sup>1</sup>

### 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).
3. 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**, 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, 11/08/2023

#### 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