

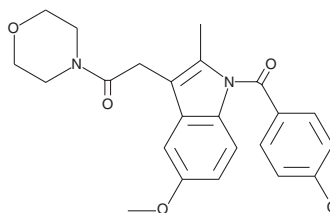
# Product Information



## IMMA

Item No. 70275

**Formal Name:** 1-(4-chlorobenzoyl)-5-methoxy-2-methyl-1H-indole-3-acetic acid, morpholineamide  
**Synonyms:** BML-190, Indomethacin Morpholinylamide  
**MF:** C<sub>23</sub>H<sub>23</sub>ClN<sub>2</sub>O<sub>4</sub>  
**FW:** 426.9  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that IMMA be stored as supplied at -20°C. It should be stable for at least one year.

IMMA is supplied as a crystalline solid. A stock solution may be made by dissolving the IMMA in an organic solvent purged with an inert gas. IMMA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of IMMA in these solvents is approximately 1.25, 12.75, 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 IMMA can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of IMMA in PBS (acidic or basic) is approximately 50 µg/ml. We do not recommend storing the aqueous solution for more than one day.

IMMA is a selective CB<sub>2</sub> receptor agonist. The binding constant for the CB<sub>2</sub> receptor is 435 nM compared to >20,000 nM for the cannabinoid receptor.<sup>1</sup>

### Reference

1. Gallant, M., Dufresne, C., Gareau, Y., *et al.* New class of potent ligands for the human peripheral cannabinoid receptor. *Bioorg. Medicinal Chem. Letters* **6**, 2263-2268 (1996).

### Related Products

1-Arachidonoyl Glycerol - Item No. 62150 • 2-Arachidonoyl Glycerol - Item No. 62160 • 2-Arachidonoyl Glycerol ether - Item No. 62165 • 2-Linoleoyl Glycerol - Item No. 62260 • AM251 - Item No. 71670 • Arachidonoyl Ethanolamide - Item No. 90050 • Arachidonoyl Glycine - Item No. 90051 • Arvanil - Item No. 90052 • Arachidonoyl Dopamine - Item No. 90057 • R-1 Methanandamide - Item No. 90070 • S-1 Methanandamide - Item No. 90072 • R-2 Methanandamide - Item No. 90074 • S-2 Methanandamide - Item No. 90076 • Cannabidiol - Item No. 90080 • Mead Acid Ethanolamide - Item No. 90195 • Palmitoyl Ethanolamide - Item No. 90350 • Arachidonoyl Cyclopropylamide - Item No. 91053

**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, 08/24/2010

### 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