

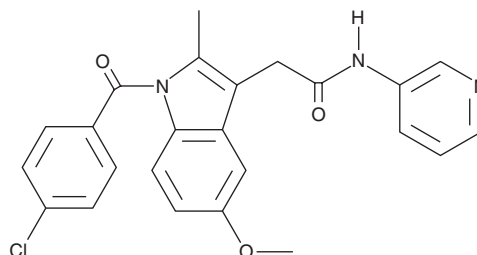
# PRODUCT INFORMATION



## N-(3-pyridyl)-Indomethacin amide

Item No. 70274

**CAS Registry No.:** 261766-29-4  
**Formal Name:** N-(3-pyridyl)-1-(4-chlorobenzoyl)-5-methoxy-2-methyl-1H-indole-3-acetamide  
**Synonym:** N-3PyIA  
**MF:** C<sub>24</sub>H<sub>20</sub>ClN<sub>3</sub>O<sub>3</sub>  
**FW:** 433.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 240 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

N-(3-pyridyl)-Indomethacin amide is supplied as a crystalline solid. A stock solution may be made by dissolving the N-(3-pyridyl)-indomethacin amide in the solvent of choice, which should be purged with an inert gas. N-(3-pyridyl)-Indomethacin amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-(3-pyridyl)-indomethacin amide in these solvents is approximately 3, 5.4, and 8.8 mg/ml, respectively.

N-(3-pyridyl)-Indomethacin amide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-(3-pyridyl)-indomethacin amide should first be dissolved in DMF and then diluted with the aqueous buffer of choice. N-(3-pyridyl)-Indomethacin amide has a solubility of approximately 5.2 mg/ml in a 5:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

N-(3-pyridyl)-indomethacin amide (N-3PyIA) is one of several aromatic amides of indomethacin reported to be potent and selective reversible inhibitors of COX-2.<sup>1</sup> N-3PyIA selectively inhibits human recombinant COX-2 with an IC<sub>50</sub> of 0.052 μM. It is about 1,300 times less potent as an inhibitor of ovine recombinant COX-1.<sup>1</sup> However, N-3PyIA is not selective as an inhibitor of ovine seminal vesicular<sup>2</sup>

### References

1. Kalgutkar, A.S., Marnett, A.B., Crews, B.C., *et al.* Ester and amide derivatives of the nonsteroidal antiinflammatory drug, indomethacin, as selective cyclooxygenase-2 inhibitors. *J. Med. Chem.* **43(15)**, 2860-2870 (2000).
2. Johnson, J.L. [Unpublished] (1999).

#### 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, 04/02/2024

#### 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