

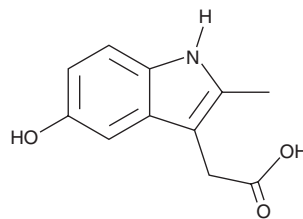
PRODUCT INFORMATION



O-Desmethyl-N-deschlorobenzoyl Indomethacin

Item No. 34581

CAS Registry No.: 50995-53-4
Formal Name: 5-hydroxy-2-methyl-1H-indole-3-acetic acid
MF: $C_{11}H_{11}NO_3$
FW: 205.2
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 231, 230, 279 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

O-Desmethyl-N-deschlorobenzoyl indomethacin is supplied as a solid. A stock solution may be made by dissolving the O-desmethyl-N-deschlorobenzoyl indomethacin in the solvent of choice, which should be purged with an inert gas. O-Desmethyl-N-deschlorobenzoyl indomethacin is soluble in ethanol.

Description

O-Desmethyl-N-deschlorobenzoyl indomethacin is a metabolite of the non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor indomethacin (Item No. 70270).¹ It is formed from indomethacin in isolated rabbit hepatocytes. O-Desmethyl-N-deschlorobenzoyl indomethacin (600 μM) decreases the viability of HL-60 leukemia cells when cultured with glucose oxidase.² It has also been used in the synthesis of prostaglandin D_2 (PGD₂; Item No. 12010) receptor antagonists.³

References

1. Evans, M.A., Papazafiratou, C., Bhat, R., *et al.* Indomethacin metabolism in isolated neonatal and fetal rabbit hepatocytes. *Pediatr. Res.* **15**(11), 1406-1410 (1981).
2. Morgan, A.G.M., Babu, D., Michail, K., *et al.* An evaluation of myeloperoxidase-mediated bio-activation of NSAIDs in promyelocytic leukemia (HL-60) cells for potential cytotoxic selectivity. *Toxicol. Lett.* **280**, 48-56 (2017).
3. Torisu, K., Kobayashi, K., Iwahashi, M., *et al.* Discovery of new chemical leads for prostaglandin D_2 receptor antagonists. *Bioorg. Med. Chem. Lett.* **14**(17), 4557-4562 (2004).

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, 11/23/2022

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