

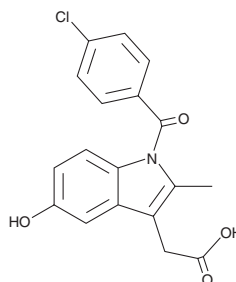
# PRODUCT INFORMATION



## 5-hydroxy Indomethacin

Item No. 34578

**CAS Registry No.:** 2504-32-7  
**Formal Name:** 1-(4-chlorobenzoyl)-5-hydroxy-2-methyl-1H-indole-3-acetic acid  
**Synonym:** O-Desmethyl Indomethacin  
**MF:** C<sub>18</sub>H<sub>14</sub>ClNO<sub>4</sub>  
**FW:** 343.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 231 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

5-hydroxy Indomethacin is supplied as a solid. A stock solution may be made by dissolving the 5-hydroxy indomethacin in the solvent of choice, which should be purged with an inert gas. 5-hydroxy Indomethacin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 5-hydroxy indomethacin in these solvents is 3, 15, and 30 mg/ml, respectively.

### Description

5-hydroxy Indomethacin is a metabolite of indomethacin (Item No. 70270).<sup>1</sup> It is formed from indomethacin in rabbit hepatic microsomes.

### Reference

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).

#### 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, 11/09/2022

#### 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](http://WWW.CAYMANCHEM.COM)