

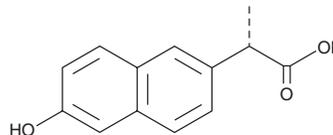
PRODUCT INFORMATION



(S)-O-Desmethyl Naproxen

Item No. 34199

CAS Registry No.: 52079-10-4
Formal Name: (αS)-6-hydroxy-α-methyl-2-naphthaleneacetic acid
Synonym: (S)-6-Desmethyl Naproxen
MF: C₁₃H₁₂O₃
FW: 216.2
Purity: ≥98%
UV/Vis.: λ_{max}: 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

(S)-O-Desmethyl naproxen is supplied as a solid. A stock solution may be made by dissolving the (S)-O-desmethyl naproxen in the solvent of choice, which should be purged with an inert gas. (S)-O-Desmethyl naproxen is soluble in DMSO.

Description

(S)-O-Desmethyl naproxen is a metabolite of the non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor (+)-naproxen ((S)-naproxen; Item No. 70290).¹ It is formed from (S)-naproxen by the cytochrome P450 (CYP) isoforms CYP1A2 and CYP2C9.

Reference

1. Miners, J.O., Soultter, S., Tukey, R.H., *et al.* Cytochromes P450, 1A2, and 2C9 are responsible for the human hepatic O-demethylation of R- and S-naproxen. *Biochem. Pharmacol.* **51(8)**, 1003-1008 (1996).

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, 10/03/2022

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