

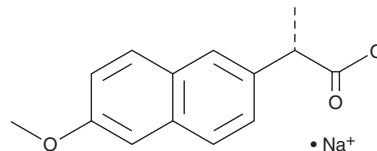
PRODUCT INFORMATION



(+)-Naproxen (sodium salt)

Item No. 37554

CAS Registry No.: 26159-34-2
Formal Name: (αS)-6-methoxy-α-methyl-2-naphthaleneacetic acid, monosodium salt
Synonyms: CG 3117, (S)-Naproxen
MF: C₁₄H₁₃O₃ • Na
FW: 252.2
Purity: ≥95%
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

(+)-Naproxen (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the (+)-naproxen (sodium salt) in the solvent of choice, which should be purged with an inert gas. (+)-Naproxen (sodium salt) is soluble in DMSO and methanol. (+)-Naproxen (sodium salt) is slightly soluble in ethanol.

Description

(+)-Naproxen is a non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor.¹ It inhibits COX-1 and COX-2 in COS-1 cells expressing the human enzymes (IC₅₀s = 4.8 and 28.4 μM, respectively). (+)-Naproxen inhibits thromboxane B₂ (TXB₂; Item No. 19030) and prostaglandin E₂ (PGE₂; Item No. 14010) production *ex vivo* in isolated rat blood (IC₅₀s = 5 and 13 μM, respectively).² Formulations containing (+)-naproxen have been used in the treatment of arthritis, osteoarthritis, gout, and ankylosing spondylitis.

References

1. Laneuville, O., Breuer, D.K., DeWitt, D.L., *et al.* Differential inhibition of human prostaglandin endoperoxide H synthases-1 and -2 by nonsteroidal anti-inflammatory drugs. *J. Pharmacol. Exp. Ther.* **271**(2), 927-934 (1994).
2. Krekels, E.H.J., Angesjö, M., Sjögren, I., *et al.* Pharmacokinetic-pharmacodynamic modeling of the inhibitory effects of naproxen on the time-courses of inflammatory pain, fever, and the *ex vivo* synthesis of TXB₂ and PGE₂ in rats. *Pharm. Res.* **28**(7), 1561-1576 (2011).

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, 10/27/2023

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