

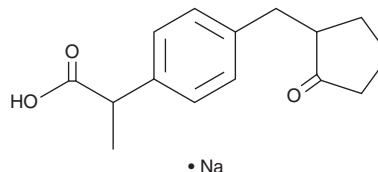
PRODUCT INFORMATION



Loxoprofen (sodium salt)

Item No. 35088

CAS Registry No.: 80382-23-6
Formal Name: α -methyl-4-[(2-oxocyclopentyl)methyl]-benzeneacetic acid, monosodium salt
MF: $C_{15}H_{18}O_3 \cdot Na$
FW: 269.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 223 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

Loxoprofen (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the loxoprofen (sodium salt) in the solvent of choice, which should be purged with an inert gas. Loxoprofen (sodium salt) is soluble in DMSO.

Description

Loxoprofen is a non-steroidal anti-inflammatory drug (NSAID) and prodrug form of loxoprofen-SRS.¹ It inhibits COX-1 and COX-2 in isolated human whole blood ($IC_{50}s = 6.5$ and $13.5 \mu M$, respectively). Loxoprofen reduces acetic acid-induced writhing in mice and carrageenan-induced paw edema and LPS-induced fever in rats ($ED_{30}s = 20.1, 0.7$, and 2.79 mg/kg, respectively).² Formulations containing loxoprofen have been used in the treatment of pain associated with osteoarthritis, myalgia, and post-traumatic swelling.

References

1. Riendeau, D., Salem, M., Styhler, A., *et al.* Evaluation of loxoprofen and its alcohol metabolites for potency and selectivity of inhibition of cyclooxygenase-2. *Bioorg. Med. Chem. Lett.* **14**(5), 1201-1203 (2004).
2. Futaki, N., Yoshikawa, K., Hamasaka, Y., *et al.* NS-398, a novel non-steroidal anti-inflammatory drug with potent analgesic and antipyretic effects, which causes minimal stomach lesions. *Gen. Pharmacol.* **24**(1), 105-110 (1993).

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, 11/15/2022

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