

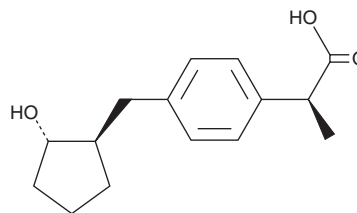
PRODUCT INFORMATION



Loxoprofen-SRS

Item No. 38122

CAS Registry No.: 83648-76-4
Formal Name: 4-[[[(1R,2S)-2-hydroxycyclopentyl]methyl]- α S-methyl-benzeneacetic acid
MF: C₁₅H₂₀O₃
FW: 248.3
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Loxoprofen-SRS is supplied as a solid. A stock solution may be made by dissolving the loxoprofen-SRS in the solvent of choice, which should be purged with an inert gas. Loxoprofen-SRS is soluble in methanol and acetonitrile.

Description

Loxoprofen-SRS is a non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor (IC_{50} s = 0.5 and 0.39 for recombinant human COX-1 and COX-2, respectively) and an active metabolite of the prodrug loxoprofen (Item No. 35088).¹ Intravenous administration of loxoprofen-SRS inhibits acetic acid-induced writhing, reduces carrageenan-induced mechanical hyperalgesia, and decreases carrageenan-induced paw edema in rats (ED_{50} s = 0.43, 1.28, and 8.01 mg/kg, respectively).²

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. Ye, H., Lv, T., Min, T., *et al.* HR1405-01, a safe intravenous NSAID with superior anti-inflammatory and analgesic activities in preclinical trials. *Eur. J. Med. Chem.* **235**:114258, (2022).

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, 03/15/2024

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