

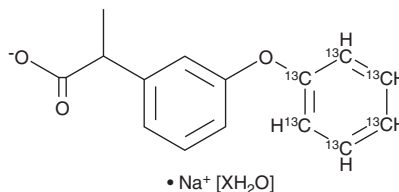
PRODUCT INFORMATION



Fenopropfen-¹³C₆ (sodium salt hydrate)

Item No. 41683

CAS Registry No.: 2687960-29-6
Formal Name: α-methyl-3-(phenoxy-1,2,3,4,5,6-¹³C₆)-benzeneacetic acid, monosodium salt, hydrate
MF: C₉[¹³C]₆H₁₃O₃ • Na [XH₂O]
FW: 270.2
Purity: ≥98%
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.

Description

Fenopropfen-¹³C₆ is intended for use as an internal standard for the quantification of fenopropfen (Item No. 23835) by GC- or LC-MS. Fenopropfen is a non-steroidal anti-inflammatory drug (NSAID).¹ It inhibits spontaneous and estradiol-stimulated prostaglandin F_{2α} (PGF_{2α}; Item Nos. 16010 | 16020) release by 84% in isolated rat uterine horns when used at a concentration of 33 μM. Fenopropfen (10 mg/kg, i.p.) reduces serum PGF_{2α} levels in rats. It reduces acetic acid-induced writhing and stretching and formalin-induced licking behaviors in mice, however, it also induces the formation of gastric lesions in mice.² Formulations containing fenopropfen have been used in the treatment of pain and the signs and symptoms of rheumatoid arthritis and osteoarthritis.

References

1. Patrono, C., Ciabattoni, G., and Grossi-Belloni, D. In vitro and in vivo inhibition of prostaglandin synthesis by fenopropfen, a non steroid anti-inflammatory drug. *Pharmacol. Res. Commun.* **6**(5), 509-518 (1974).
2. Agotegaray, M., Gumilar, F., Boeris, M., et al. Enhanced analgesic properties and reduced ulcerogenic effect of a mononuclear copper(II) complex with fenopropfen in comparison to the parent drug: Promising insights in the treatment of chronic inflammatory diseases. *Biomed. Res. Int.* 505987 (2014).

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, 10/28/2024

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