

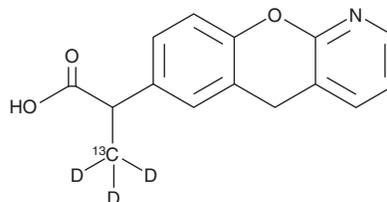
PRODUCT INFORMATION



Pranoprofen-¹³C-d₃

Item No. 34009

Formal Name:	α-methyl- ¹³ C-d ₃ -5H-[1]benzopyrano[2,3-b]pyridine-7-acetic acid
Synonyms:	Pranoprofen- ¹³ C-d ₃ , Y 8004- ¹³ C-d ₃
MF:	C ₁₄ [¹³ C]H ₁₀ D ₃ NO ₃
FW:	259.3
Chemical Purity:	≥95% (Pranoprofen)
Deuterium Incorporation:	≥99% deuterated forms (d ₁ -d ₃); ≤1% d ₀
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

Pranoprofen-¹³C-d₃ is intended for use as an internal standard for the quantification of Pranoprofen (Item No. 17875) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Description

Pranoprofen is a non-steroidal anti-inflammatory drug (NSAID) and COX-1 inhibitor (IC₅₀ = 3.1 μM).^{1,2} It decreases *B. pertussis*-induced production of prostaglandin E₂ (PGE₂; Item No. 14010) in isolated rat peritoneal leukocytes (IC₅₀ = 0.39 μM).² Ocular application of pranoprofen reduces corneal epithelial damage in a mouse model of corneal alkali burns.¹

References

1. Chen, M., Gureeye, A., Cissé, Y., *et al.* The therapeutic effects and possible mechanism of pranoprofen in mouse model of corneal alkali burns. *J. Ophthalmol.* 7485912 (2020).
2. Naka, Y., Yamada, I., Ochi, H., *et al.* Structural optimization of alkyl-substituted 2-(7-5H-[1]benzopyrano[2,3-b]pyridyl)propionic acid as anti-inflammatory agents. *Bioorg. Med. Chem. Lett.* 5(9), 959-962 (1995).

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, 10/19/2022

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