

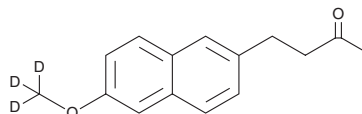
PRODUCT INFORMATION



Nabumetone-d₃

Item No. 30721

CAS Registry No.: 1216770-08-9
Formal Name: 4-(6-(methoxy-d₃)naphthalen-2-yl)butan-2-one
MF: C₁₅H₁₃D₃O₂
FW: 231.3
Chemical Purity: ≥98% (Nabumetone)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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

Nabumetone-d₃ is intended for use as an internal standard for the quantification of nabumetone (Item No. 20251) 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).

Nabumetone-d₃ is supplied as a solid. A stock solution may be made by dissolving the nabumetone-d₃ in the solvent of choice, which should be purged with an inert gas. Nabumetone-d₃ is slightly soluble in chloroform and ethyl acetate.

Description

Nabumetone is a non-steroidal anti-inflammatory drug (NSAID) and a prodrug form of 6-methoxy naphthalene acetic acid (Item No. 70620).¹ Nabumetone (5-45 mg/kg) reduces carrageenan-induced edema in rats, as well as UV-induced erythema in guinea pigs (ED₅₀ = 11 mg/kg).² It reduces phenyl-*p*-quinone-induced writhing in mice (ED₅₀ = 152 mg/kg). Nabumetone (50, 100, and 200 mg/kg) reduces endotoxin-induced pyresis in rabbits.

References

1. Cryer, B. and Feldman, M. Cyclooxygenase-1 and cyclooxygenase-2 selectivity of widely used nonsteroidal anti-inflammatory drugs. *Am. J. Med.* **104**(5), 413-421 (1998).
2. Boyle, E.A., Freeman, P.C., Mangan, F.R., *et al.* Nabumetone (BRL 14777, 4-[6-methoxy-2-naphthyl]-butan-2-one): a new anti-inflammatory agent. *J. Pharm. Pharmacol.* **34**(9), 562-569 (1982).

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, 11/03/2022

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