

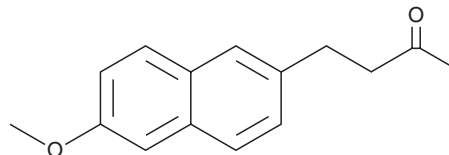
PRODUCT INFORMATION



Nabumetone

Item No. 20251

CAS Registry No.: 42924-53-8
Formal Name: 4-(6-methoxy-2-naphthalenyl)-2-butanone
Synonym: BRL 14777
MF: $C_{15}H_{16}O_2$
FW: 228.3
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 230 nm
Supplied as: A crystalline 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 is supplied as a crystalline solid. A stock solution may be made by dissolving the nabumetone in the solvent of choice. Nabumetone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of nabumetone in these solvents is approximately 20 mg/ml in ethanol and 30 mg/ml in DMSO and DMF.

Nabumetone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nabumetone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nabumetone has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

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 ($ED_{50} = 11$ mg/kg).² It reduces phenyl-*p*-quinone-induced writhing in mice ($ED_{50} = 152$ 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, 12/19/2022

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