

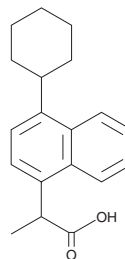
PRODUCT INFORMATION



Vedaprofen

Item No. 15960

CAS Registry No.: 71109-09-6
Formal Name: 4-cyclohexyl- α -methyl-1-naphthaleneacetic acid
Synonym: Quadrisol
MF: C₁₉H₂₂O₂
FW: 282.4
Purity: \geq 95%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 228, 288 nm



Laboratory Procedures

For long term storage, we suggest that vedaprofen be stored as supplied at -20°C. It should be stable for at least two years.

Vedaprofen is supplied as a crystalline solid. A stock solution may be made by dissolving the vedaprofen in the solvent of choice. Vedaprofen is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of vedaprofen in ethanol and DMSO is approximately 10 mg/ml and 25 mg/ml in DMF.

Vedaprofen is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Vedaprofen is a non-steroidal anti-inflammatory drug (NSAID) commonly used in veterinary medicine to combat pain, inflammation, and fever associated with acute and chronic musculoskeletal disorders in horses and dogs.¹ In addition to inhibiting COX and reducing prostaglandin H₂ synthesis, vedaprofen is reported to block the activity of the *E. coli* DNA polymerase III β subunit, preventing DNA replication and repair.²

References

1. Fantoni, D.T., Ida, K.K., de Almeida, T.I., *et al.* A comparison of pre and post-operative vedaprofen with ketoprofen for pain control in dogs. *BMC Vet. Res.* **11:24**, (2015).
2. Yin, Z., Wang, Y., Whittell, L.R., *et al.* DNA replication is the target for the antibacterial effects of nonsteroidal anti-inflammatory drugs. *Chem. Biol.* **21(4)**, 481-487 (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, 01/28/2016

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