

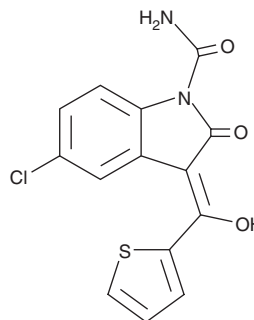
PRODUCT INFORMATION



Tenidap

Item No. 17413

CAS Registry No.: 120210-48-2
Formal Name: (3Z)-5-chloro-2,3-dihydro-3-(hydroxy-2-thienylmethylene)-2-oxo-1H-indole-1-carboxamide
Synonym: CP 66,248
MF: C₁₄H₉ClN₂O₃S
FW: 320.8
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 273, 370 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

Tenidap is supplied as a crystalline solid. A stock solution may be made by dissolving the tenidap in the solvent of choice. Tenidap is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of tenidap in these solvents is approximately 20 and 10 mg/ml, respectively. It is also slightly soluble in ethanol.

Description

Tenidap is a COX-1 selective non-steroidal anti-inflammatory drug (IC₅₀s = <0.03, 1.2, and >30 μM for COX-1, COX-2, and 5-lipoxygenase (5-LO), respectively).¹ It has anti-inflammatory and antirheumatic properties.^{1,2} *In vitro*, it inhibits prostaglandin D₂ (PGD₂), leukotriene B₄ (LTB₄), and prostaglandin E₂ (PGE₂) synthesis (IC₅₀s = 0.02, 18, and 32 μM, respectively).^{3,4} Tenidap also reversibly and dose-dependently activates hK_{ir}2.3 channels in CHO cells (EC₅₀ = 402 nM) and inhibits fatty acid amide hydrolase (FAAH) activity.^{5,6} A formulation containing tenidap was not approved for rheumatoid and osteoarthritis by the FDA due to adverse effects, including bone mineralization loss, as well as liver and kidney toxicity.

References

1. Kirchner, T., Argentieri, D.C., Barbone, A.G., *et al.* Evaluation of the antiinflammatory activity of a dual cyclooxygenase-2 selective/5-lipoxygenase inhibitor, RWJ 63556, in a canine model of inflammation. *J. Pharmacol. Exp. Ther.* **282**(2), 1094-1101 (1997).
2. Blackburn, W.D., Jr., Prupas, H.M., Silverfield, J.C., *et al.* Tenidap in rheumatoid arthritis. A 24-week double-blind comparison with hydroxychloroquine-plus-piroxicam, and piroxicam alone. *Arthritis Rheum.* **38**(10), 1447-1456 (1995).
3. Moilanen, E., Alanko, J., Asmawi, M.Z., *et al.* CP-66,248, a new anti-inflammatory agent, is a potent inhibitor of leukotriene B₄ and prostanoid synthesis in human polymorphonuclear leucocytes *in vitro*. *Eicosanoids* **1**(1), 35-39 (1988).
4. Moore, P.F., Larson, D.L., Otterness, I.G., *et al.* Tenidap, a structurally novel drug for the treatment of arthritis: Antiinflammatory and analgesic properties. *Inflamm. Res.* **45**(2), 54-61 (1996).
5. Liu, Y., Liu, D., Printzenhoff, D., *et al.* Tenidap, a novel anti-inflammatory agent, is an opener of the inwardly rectifying K⁺ channel hK_{ir}2.3. *Eur. J. Pharmacol.* **435**(2-3), 153-160 (2002).
6. Bertolacci, L., Romeo, E., Veronesi, M., *et al.* A binding site for nonsteroidal anti-inflammatory drugs in fatty acid amide hydrolase. *Journal of the American Chemical Society* **135**(1), 22-25 (2013).

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, 12/02/2022

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