

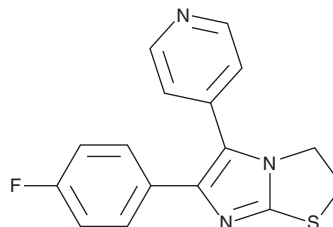
PRODUCT INFORMATION



SKF 86002

Item No. 30202

CAS Registry No.: 72873-74-6
Formal Name: 6-(4-fluorophenyl)-2,3-dihydro-5-(4-pyridinyl)-imidazo[2,1-b]thiazole
MF: C₁₆H₁₂FN₃S
FW: 297.4
Purity: ≥98%
UV/Vis.: λ_{max}: 238 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

SKF 86002 is an anti-inflammatory agent.¹⁻⁴ It inhibits rat seminal vesicle prostaglandin H₂ (PGH₂) synthase (IC₅₀ = 120 μM), as well as prostanoid production by rat basophilic leukemia (RBL-1) cells and human monocytes (IC₅₀s = 70 and 1 μM, respectively).¹ SKF 86002 inhibits leukotriene B₄ (LTB₄) and LTC₄ production induced by A23187 (Item No. 11016) in human neutrophils and monocytes, respectively (IC₅₀ = 20 μM for both). It also inhibits LPS-induced IL-1 production in human monocytes (IC₅₀ = 1.3 μM).² SKF 86002 (10, 30, and 90 mg/kg) reduces hindleg volume in rat models of adjuvant- or collagen-induced arthritis.³ It also decreases serum levels of TNF-α and increases survival in a mouse model of LPS and galactosamine-induced endotoxic shock when administered at a dose of 100 mg/kg.⁴

References

1. Griswold, D.E., Marshall, P.J., Webb, E.F., *et al.* SK&F 86002: A structurally novel anti-inflammatory agent that inhibits lipoxygenase- and cyclooxygenase-mediated metabolism of arachidonic acid. *Biochem. Pharmacol.* **36(20)**, 3463-3470 (1987).
2. Lee, J.C., Griswold, D.E., Votta, B., *et al.* Inhibition of monocyte IL-1 production by the anti-inflammatory compound, SK&F 86002. *Int. J. Immunopharmacol.* **10(7)**, 835-843 (1988).
3. DiMartino, M.J., Griswold, D.E., Berkowitz, B.A., *et al.* Pharmacologic characterization of the antiinflammatory properties of a new dual inhibitor of lipoxygenase and cyclooxygenase. *Agents Actions* **20(1-2)**, 113-123 (1987).
4. Badger, A.M., Olivera, D., Talmadge, J.E., *et al.* Protective effect of SK&F 86002, a novel dual inhibitor of arachidonic acid metabolism, in murine models of endotoxin shock: Inhibition of tumor necrosis factor as a possible mechanism of action. *Circ. Shock* **27(1)**, 51-61 (1989).

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, 05/19/2020

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