

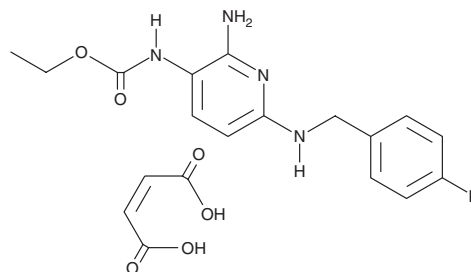
PRODUCT INFORMATION



Flupirtine (maleate)

Item No. 16674

CAS Registry No.: 75507-68-5
Formal Name: N-[2-amino-6-[[[(4-fluorophenyl)methyl]amino]-3-pyridinyl]-carbamic acid, ethyl ester, 2Z-butenedioate
MF: C₁₅H₁₇FN₄O₂ • C₄H₄O₄
FW: 420.4
Purity: ≥95%
UV/Vis.: λ_{max}: 204, 250, 348 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

Flupirtine (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the flupirtine (maleate) in the solvent of choice, which should be purged with an inert gas. Flupirtine (maleate) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of flupirtine (maleate) in these solvents is approximately 20 and 30 mg/ml, respectively.

Flupirtine (maleate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, flupirtine (maleate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Flupirtine (maleate) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Flupirtine is an activator of voltage-gated potassium channel 7 (K_v7/KCNQ).¹⁻³ It induces relaxation of precontracted pulmonary arteries isolated from wild-type and serotonin transporter-overexpressing (SERT⁺) mice.² Flupirtine (30 mg/kg per day) decreases mean right ventricular pressure and right ventricular hypertrophy in hypoxia-induced and SERT⁺ mouse models of pulmonary arterial hypertension. It increases the paw withdrawal threshold in a rat model of streptozotocin-induced diabetic neuropathy when administered at a dose of 10 mg/kg and increases paw withdrawal latency in a rat model of carrageenan-induced paw inflammation when used in combination with morphine.³ Flupirtine also indirectly antagonizes NMDA receptors via its effects on potassium channels.^{1,4}

References

- Devulder, J. Flupirtine in pain management: Pharmacological properties and clinical use. *CNS Drugs* **25(10)**, 867-881 (2010).
- Morecroft, I., Murray, A., Nilsen, M., et al. Treatment with the K_v7 potassium channel activator flupirtine is beneficial in two independent mouse models of pulmonary hypertension. *Br. J. Pharmacol.* **157(7)**, 1241-1249 (2009).
- Goodchild, C.S., Kolosov, A., Tucker, A.P., et al. Combination therapy with flupirtine and opioid: Studies in rat pain models. *Pain Med.* **9(7)**, 928-938 (2008).
- Kornhuber, J., Bleich, S., Wiltfang, J., et al. Flupirtine shows functional NMDA receptor antagonism by enhancing Mg²⁺ block via activation of voltage independent potassium channels. *J. Neural Transm. (Vienna)* **106(9-10)**, 857-867 (1999).

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, 11/30/2022

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