

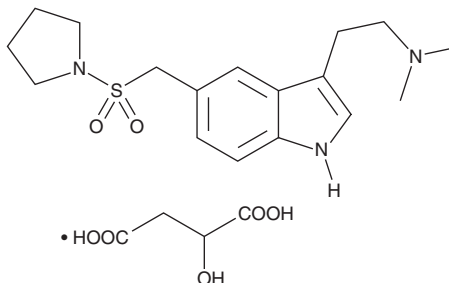
PRODUCT INFORMATION

Almotriptan (malate)

Item No. 25646

CAS Registry No.: 181183-52-8
Formal Name: 2-hydroxy-butanedioic acid, compd. with N,N-dimethyl-5-[(1-pyrrolidinylsulfonyl)methyl]-1H-indole-3-ethanamine

Synonym: PNU 180638E
MF: C₁₇H₂₅N₃O₂S • C₄H₆O₅
FW: 469.6
Purity: ≥98%
UV/Vis.: λ_{max}: 229 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

Almotriptan (malate) is supplied as a crystalline solid. A stock solution may be made by dissolving the almotriptan (malate) in the solvent of choice. Almotriptan (malate) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of almotriptan (malate) in these solvents is approximately 0.1 and 0.2 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of almotriptan (malate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of almotriptan (malate) in PBS, pH 7.2, is approximately 20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Almotriptan is an agonist of the serotonin (5-HT) receptor subtypes 5-HT_{1B} and 5-HT_{1D} (IC₅₀s = 12 and 13 nM, respectively, in a radioligand binding assay).¹ It is selective for human 5-HT_{1B} and 5-HT_{1D} receptors over rat 5-HT_{1A} and human 5-HT_{2A} and 5-HT₄ receptors (IC₅₀s = 0.85, 25.1, and 140 μM, respectively). Almotriptan induces contractions in isolated canine saphenous veins (EC₅₀ = 394 nM) but not isolated rabbit renal or mesenteric arteries. It increases carotid vascular resistance in anesthetized cats (ED₁₀₀ = 11 μg/kg, i.v.) without increasing blood pressure or heart rate.² Formulations containing almotriptan have been used in the treatment of migraine headaches.

References

1. Bou, J., Domènech, T., Puig, J., *et al.* Pharmacological characterization of almotriptan: An indolic 5-HT receptor agonist for the treatment of migraine. *Eur. J. Pharmacol.* **410(1)**, 33-41 (2000).
2. Gras, J., Bou, J., Llenas, J., *et al.* Functional profile of almotriptan in animal models predictive of antimigraine activity. *Eur. J. Pharmacol.* **410(1)**, 43-51 (2000).

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