

PRODUCT INFORMATION



Dehydro Amlodipine (fumarate)

Item No. 26622

CAS Registry No.: 2138811-33-1
Formal Name: 2-[(2-aminoethoxy)methyl]-4-(2-chlorophenyl)-6-methyl-3,5-pyridinedicarboxylic acid, 3-ethyl 5-methyl ester, 2-butenedioate

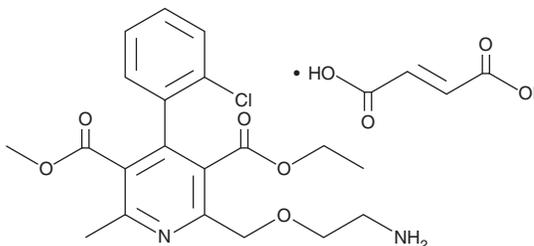
MF: C₂₀H₂₃ClN₂O₅ • C₄H₄O₄
FW: 522.9

Purity: ≥98%

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

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

Dehydro amlodipine (fumarate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dehydro amlodipine (fumarate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Dehydro amlodipine (fumarate) has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Dehydro amlodipine is a potential impurity found in commercial preparations of amlodipine.^{1,2} It is a degradation product formed during storage or when amlodipine (Item No. 14838) is subjected to oxidative or acidic conditions or undergoes photo or electrochemical degradation.

References

1. Kawabe, Y., Nakamura, H., Hino, E., *et al.* Photochemical stabilities of some dihydropyridine calcium-channel blockers in powdered pharmaceutical tablets. *J. Pharm. Biomed. Anal.* **47**(3), 618-624 (2008).
2. Stojilković, Ž.Z., Jadranin, M.B., Durić, S.L.J., *et al.* Investigation of forced and total degradation products of amlodipine besylate by liquid chromatography and liquid chromatography-mass spectrometry. *Chem. Ind. Chem. Eng. Q.* **20**(2), 295-304 (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, 11/29/2022

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