

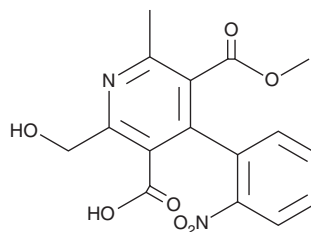
PRODUCT INFORMATION



Hydroxydehydro Nifedipine Carboxylate

Item No. 34406

CAS Registry No.: 34783-31-8
Formal Name: 2-(hydroxymethyl)-6-methyl-4-(2-nitrophenyl)-3,5-pyridinedicarboxylic acid, 5-methyl ester
MF: $C_{16}H_{14}N_2O_7$
FW: 346.3
Purity: $\geq 95\%$
Supplied as: A 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

Hydroxydehydro nifedipine carboxylate is supplied as a solid. A stock solution may be made by dissolving the hydroxydehydro nifedipine carboxylate in the solvent of choice, which should be purged with an inert gas. Hydroxydehydro nifedipine carboxylate is slightly soluble in DMSO.

Description

Hydroxydehydro nifedipine carboxylate is a metabolite of nifedipine (Item No. 11106).¹

Reference

1. Roosemalen, M.C.M., Soons, P.A., Funaki, T., *et al.* High-performance liquid chromatographic determination of the polar metabolites of nifedipine in plasma, blood and urine. *J. Chromatogr.* **565**(1-2), 516-522 (1991).

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, 11/22/2022

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