

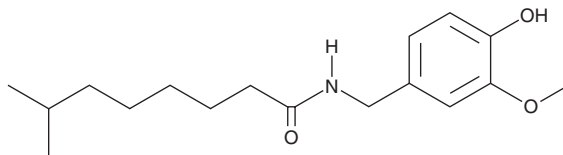
PRODUCT INFORMATION



Nordihydrocapsaicin

Item No. 33407

CAS Registry No.: 28789-35-7
Formal Name: N-[(4-hydroxy-3-methoxyphenyl)methyl]-7-methyl-octanamide
Synonym: NDHC
MF: $C_{17}H_{27}NO_3$
FW: 293.4
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Nordihydrocapsaicin is supplied as a crystalline solid. A stock solution may be made by dissolving the nordihydrocapsaicin in the solvent of choice, which should be purged with an inert gas. Nordihydrocapsaicin is soluble in organic solvents such as chloroform and DMSO.

Description

Nordihydrocapsaicin is a capsaicinoid and a major pungency principle that has been found in *C. annuum*.^{1,2}

References

1. Topuz, A. and Ozdemir, F. Assessment of carotenoids, capsaicinoids and ascorbic acid composition of some selected pepper cultivars (*Capsicum annuum* L.) grown in Turkey. *J. Food Compos. Anal.* **20**(7), 596-602 (2007).
2. Masada, Y., Hashimoto, K., Inoue, T., *et al.* Analysis of the pungent principles of *Capsicum annuum* by combined gas chromatography-mass spectrometry. *J. Food. Sci. Technol.* **36**(6), 858-860 (1971).

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, 12/07/2022

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