

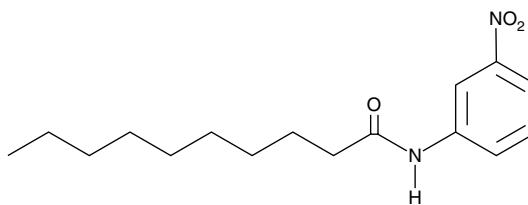
Product Information



Decanoyl *m*-Nitroaniline

Item No. 90349

CAS Registry No.: 72298-61-4
Formal Name: N-(3-nitrophenyl)-decanamide
Synonym: DemNA
MF: C₁₆H₂₄N₂O₃
FW: 292.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 202, 241 nm



Laboratory Procedures

For long term storage, we suggest that decanoyl *m*-nitroaniline (DemNA) be stored as supplied at -20°C. It should be stable for at least two years.

DemNA is supplied as a crystalline solid. A stock solution may be made by dissolving the DemNA in an organic solvent purged with an inert gas. DemNA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of DemNA in these solvents is approximately 20 mg/ml.

DemNA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DemNA should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. DemNA has a solubility of approximately 0.6 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

DemNA is one of several nitroaniline fatty acid amides which can be used to measure fatty acid amide hydrolase (FAAH) activity.¹ FAAH is a relatively unselective enzyme in that it accepts a variety of amide head groups other than the ethanolamine of its endogenous substrate anandamide (AEA). It also will hydrolyze fatty acid amides with fewer carbons and fewer double bonds than arachidonate. Exposure of DemNA to FAAH activity results in the release of the yellow colorimetric dye *m*-nitroaniline (ε = 13,500 at 410 nm). This allows the fast and convenient measurement of FAAH activity using a 96 well plate spectrophotometer.

Reference

1. Patricelli, M.P. and Cravatt, B.F. Characterization and manipulation of the acyl chain selectivity of fatty acid amide hydrolase. *Biochemistry* **40**, 6107-6115 (2001).

Related Products

Arachidonoyl *p*-Nitroaniline - Item No. 10168 • Arachidonoyl Ethanolamide - Item No. 90050 • Arachidonoyl *m*-Nitroaniline - Item No. 90059 • Fatty Acid Amide Hydrolase Polyclonal Antibody - Item No. 101600

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/20/2011