

# PRODUCT INFORMATION



## Myristyl Nicotinate

Item No. 37714

CAS Registry No.: 273203-62-6

Formal Name: 3-pyridinecarboxylic acid,  
tetradecyl ester

Synonyms: Tetradecyl Nicotinate,  
Tetradecyl Pyridine-3-carboxylate

MF:  $C_{20}H_{33}NO_2$

FW: 319.5

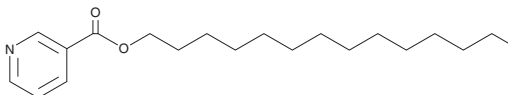
Purity:  $\geq 95\%$

UV/Vis.:  $\lambda_{max}$ : 218 nm

Supplied as: A solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Myristyl nicotinate is supplied as a solid. A stock solution may be made by dissolving the myristyl nicotinate in the solvent of choice, which should be purged with an inert gas. Myristyl nicotinate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of myristyl nicotinate in these solvents is approximately 25, 0.2, and 10 mg/ml, respectively.

### Description

Myristyl nicotinate is a prodrug form of the hydroxycarboxylic acid receptor 2 (HCA2) antagonist nicotinic acid (Item No. 37416).<sup>1</sup> Formulations containing myristyl nicotinate have been used in lotions for the treatment of epithelial photodamage.

### Reference

1. Altaani, B., Haddad, R., Al-Tarakji, A., *et al.* Hydrolysis kinetics of the prodrug myristyl nicotinate. *Pharm. Dev. Technol.* **27(10)**, 1083-1092 (2022).

#### 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, 05/31/2023

#### 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