

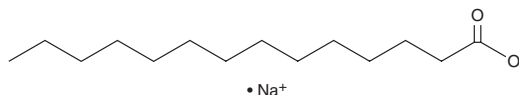
PRODUCT INFORMATION



Myristic Acid (sodium salt)

Item No. 45445

CAS Registry No.: 822-12-8
Formal Name: tetradecanoic acid, monosodium salt
Synonyms: C14:0, FA 14:0, Tetradecanoic Acid
MF: $C_{14}H_{27}O_2 \cdot Na$
FW: 250.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Myristic Acid (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the myristic Acid (sodium salt) in the solvent of choice, which should be purged with an inert gas. Myristic Acid (sodium salt) is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Myristic acid is a 14-carbon saturated fatty acid. It is incorporated into myristoyl coenzyme A (myristoyl-CoA) and transferred by N-myristoyltransferase to the N-terminal glycine of certain proteins either during translation to modify protein activity or post-translationally in apoptotic cells.^{1,2}

References

1. Bhatnagar, R.S., Fütterer, K., Waksman, G., *et al.* The structure of myristoyl-CoA: Protein N-myristoyltransferase. *Biochim. Biophys. Acta* **1441**(2-3), 162-172 (1999).
2. Martin, D.D.O., Beauchamp, E., and Berthiaume, L.G. Post-translational myristoylation: Fat matters in cellular life and death. *Biochimie* **93**(1), 18-31 (2011).

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, 03/13/2026

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