

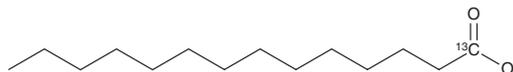
PRODUCT INFORMATION



Myristic Acid-¹³C

Item No. 29463

CAS Registry No.: 57677-52-8
Formal Name: tetradecanoic-1-¹³C acid
Synonyms: C14:0-¹³C, FA14:0-¹³C, Tetradecanoic Acid-¹³C
MF: C₁₃[¹³C]H₂₈O₂
FW: 229.4
Purity: ≥95%
Supplied as: A crystalline 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

Myristic acid-¹³C is supplied as a crystalline solid. A stock solution may be made by dissolving the myristic acid-¹³C in the solvent of choice, which should be purged with an inert gas. Myristic acid-¹³C is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of myristic acid-¹³C in ethanol and DMF is approximately 15 mg/ml and approximately 12 mg/ml in DMSO.

Description

Myristic acid-¹³C is intended for use as an internal standard for the quantification of myristic acid (Item No. 13351) by GC- or LC-MS. 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, 03/16/2026

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