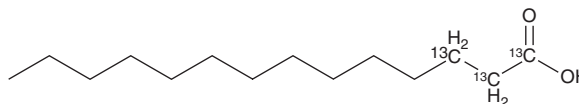


# PRODUCT INFORMATION



## Myristic Acid-<sup>13</sup>C<sub>3</sub> Item No. 45766

**CAS Registry No.:** 202114-49-6  
**Formal Name:** tetradecanoic-1,2,3-<sup>13</sup>C<sub>3</sub> acid  
**Synonyms:** C14:0-<sup>13</sup>C<sub>3</sub>, FA 14:0-<sup>13</sup>C<sub>3</sub>,  
Tetradecanoic Acid-<sup>13</sup>C<sub>3</sub>  
**MF:** C<sub>11</sub>[<sup>13</sup>C]<sub>3</sub>H<sub>28</sub>O<sub>2</sub>  
**FW:** 231.4  
**Purity:** ≥98%  
**Supplied as:** A 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-<sup>13</sup>C<sub>3</sub> is intended for use as an internal standard for the quantification of myristic acid (Item Nos. 13351 | 45445) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Myristic acid-<sup>13</sup>C<sub>3</sub> is supplied as a solid. A stock solution may be made by dissolving the myristic acid-<sup>13</sup>C<sub>3</sub> in the solvent of choice, which should be purged with an inert gas. Myristic acid-<sup>13</sup>C<sub>3</sub> is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### 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.<sup>1,2</sup>

### 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, 05/20/2026

#### 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