

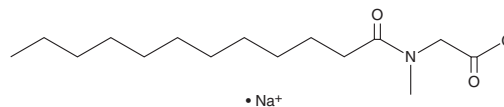
# PRODUCT INFORMATION



## Lauroylsarcosine (sodium salt)

Item No. 33718

**CAS Registry No.:** 137-16-6  
**Formal Name:** N-methyl-N-(1-oxododecyl)-glycine, monosodium salt  
**Synonyms:** Sarkosyl, SDDS, Sodium Lauroyl Sarcosinate, Sodium N-dodecanoyl Sarcosinate  
**MF:** C<sub>15</sub>H<sub>28</sub>NO<sub>3</sub> • Na  
**FW:** 293.4  
**Purity:** ≥90%  
**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

Lauroylsarcosine (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the lauroylsarcosine (sodium salt) in the solvent of choice, which should be purged with an inert gas. Lauroylsarcosine (sodium salt) is slightly soluble in ethanol and methanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of lauroylsarcosine (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. Lauroylsarcosine (sodium salt) is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Lauroylsarcosine is a fatty acid amide and detergent.<sup>1,2</sup> It has been used to solubilize recombinant and membrane proteins.

### References

1. Filip, C., Fletcher, G., Wulff, J.L., *et al.* Solubilization of the cytoplasmic membrane of *Escherichia coli* by the ionic detergent sodium-lauryl sarcosinate. *J. Bacteriol.* **115**(3), 717-722 (1973).
2. Frankel, S., Sohn, R., and Leinwand, L. The use of sarkosyl in generating soluble protein after bacterial expression. *Proc. Natl. Acad. Sci. USA* **88**(4), 1192-1196 (1991).

#### 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, 11/17/2022

#### 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