

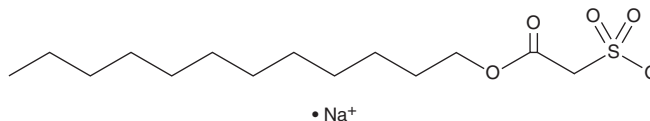
PRODUCT INFORMATION



Sodium Lauryl Sulfoacetate

Item No. 29906

CAS Registry No.: 1847-58-1
Formal Name: 2-sulfo-acetic acid, dodecyl ester, monosodium salt
MF: C₁₄H₂₇O₅S • Na
FW: 330.4
Purity: ≥95%
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.

Description

Sodium lauryl sulfoacetate is an anionic surfactant.¹ It increases proliferation of isolated chicken peripheral blood mononuclear cells (PBMCs) when used at concentrations ranging from 62.5 to 500 µg/ml.² Sodium lauryl sulfoacetate (1, 2, and 4 mg/kg) increases the blood CD4⁺ to CD8⁺ T cell ratio induced by a Newcastle disease virus vaccine in chickens. Formulations containing sodium lauryl sulfoacetate have been used as surfactants in the manufacture of cosmetics.

References

1. Gad, E.A.M., El-Sukkary, M.M.A., and Azzam, E.M.S. Surface and thermodynamic properties of octyl, dodecyl, and cetyl sulfoacetates. *Monatshefte für Chemie* **128**, 1085-1092 (1997).
2. Cheng, D., Zhu, S., and Sun, H. The immune enhancement of sodium lauryl sulfoacetate in chickens. *Vet. Med. Int.* 485060 (2010).

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, 12/08/2022

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