

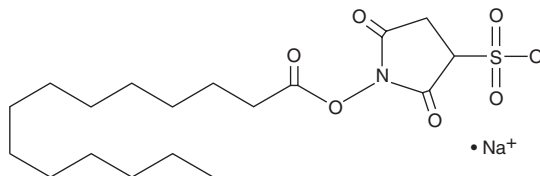
PRODUCT INFORMATION



Sulfosuccinimidyl Myristate (sodium salt)

Item No. 11212

CAS Registry No.: 162823-01-0
Formal Name: tetradecanoic acid, 2,5-dioxo-3-sulfo-1-pyrrolidinyl ester, monosodium salt
Synonym: SSM
MF: $C_{18}H_{30}NO_7S \cdot Na$
FW: 427.5
Purity: $\geq 95\%$
Supplied as: A crystalline 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

Sulfosuccinimidyl myristate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the sulfosuccinimidyl myristate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Sulfosuccinimidyl myristate (sodium salt) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of sulfosuccinimidyl myristate (sodium salt) in these solvents is approximately 2 mg/ml.

Sulfosuccinimidyl myristate (sodium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sulfosuccinimidyl myristate (sodium salt) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Sulfosuccinimidyl myristate (sodium salt) has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SSM is an inhibitor of fatty acid transport.¹ It has been used as an affinity label for certain membrane proteins that bind to fatty acids.¹

Reference

1. Harmon, C.M., Luce, P., Beth, A.H., *et al.* Labeling of adipocyte membranes by sulfo-N-succinimidyl derivatives of long-chain fatty acids: Inhibition of fatty acid transport. *J. Membrane Biol.* **121**(3), 261-268 (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, 02/22/2024

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