

PRODUCT INFORMATION



12:0 N-Biotinyl Fatty Acid

Item No. 25714

CAS Registry No.: 135447-73-3

Formal Name: 12-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]-dodecanoic acid

Synonym: N-Biotinyl Dodecanoic Acid

MF: $C_{22}H_{39}N_3O_4S$

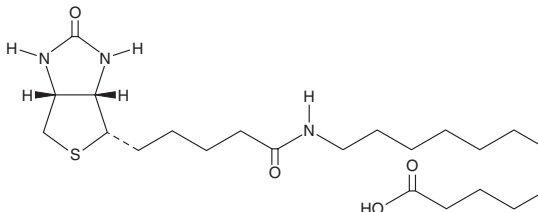
FW: 441.6

Purity: $\geq 98\%$

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

12:0 N-Biotinyl fatty acid is supplied as a solid. A stock solution may be made by dissolving the 12:0 N-biotinyl fatty acid in the solvent of choice, which should be purged with an inert gas. 12:0 N-Biotinyl fatty acid is soluble in the organic solvent DMSO.

Description

12:0 N-Biotinyl fatty acid is a biotinylated form of lauric acid (Item No. 10006626). It is an intermediate in the synthesis of N-biotinylated phenylarsonous acids.¹

Reference

1. Moaddel, R., Sharma, A., Huseni, T., *et al.* Novel biotinylated phenylarsonous acids as bifunctional reagents for spatially close thiols: Studies on reduced antibodies and the agonist binding site of reduced Torpedo nicotinic receptors. *Bioconjug. Chem.* **10**(4), 629-637 (1999).

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, 11/28/2022

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