

PRODUCT INFORMATION

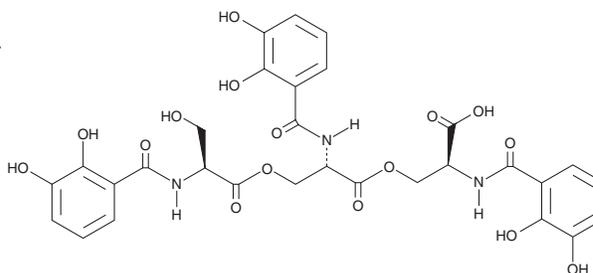


apo-Enterobactin

Item No. 31760

CAS Registry No.: 30414-16-5
Formal Name: N-(2,3-dihydroxybenzoyl)-O-[N-(2,3-dihydroxybenzoyl)-O-[N-(2,3-dihydroxybenzoyl)-L-seryl]-L-seryl]-L-serine

MF: C₃₀H₂₉N₃O₁₆
FW: 687.6
Purity: ≥85%
UV/Vis.: λ_{max}: 251 nm
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

apo-Enterobactin is supplied as a solid. A stock solution may be made by dissolving the apo-enterobactin in the solvent of choice, which should be purged with an inert gas. apo-Enterobactin is soluble in methanol and DMSO.

Description

apo-Enterobactin is a derivative of the siderophore enterobactin.¹

Reference

1. Singh, V., Yeoh, B.S., Xiao, X., *et al.* Interplay between enterobactin, myeloperoxidase and lipocalin 2 regulates *E. coli* survival in the inflamed gut. *Nat. Commun.* **6**, 7113 (2015).

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, 12/02/2022

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