

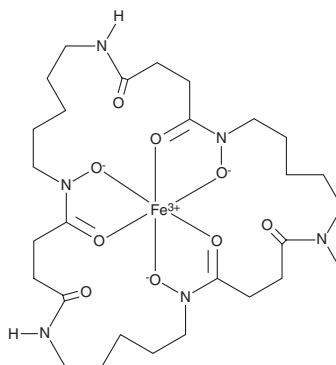
PRODUCT INFORMATION



Ferrioxamine E

Item No. 27275

CAS Registry No.: 20008-20-2
Formal Name: (OC-6-22)-[1,12,23-tri(hydroxy- κ O)-1,6,12,17,23,28-hexaazacyclotritriacontane-2,5,13,16,24,27-hexonato(3-)- κ O², κ O¹³, κ O²⁴]-iron
MF: C₂₇H₄₅FeN₆O₉
FW: 653.5
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ferrioxamine E is supplied as a solid. A stock solution may be made by dissolving the ferrioxamine E in the solvent of choice, which should be purged with an inert gas. Ferrioxamine E is soluble in organic solvents such as DMSO and methanol.

Description

Ferrioxamine E is a bacterial siderophore produced by a variety of bacteria including *S. glaucescens*, *M. luteus*, and *P. mendocina*.¹ Ferrioxamines, including ferrioxamine E, are the sole iron source for *Salmonella*, and have been used as growth enriching agents in the detection of *Salmonella* in food and industrial applications.^{2,3}

References

- Ahmed, E. and Holmstrom, S.J.M. Siderophore production by microorganisms isolated from a podzol soil profile. *Geomicrobiol. J.* **32(5)**, 397-411 (2014).
- Kingsley, R.A., Reissbrodt, R., Rabsch, W., et al. Ferrioxamine-mediated Iron(III) utilization by *Salmonella enterica*. *Appl. Environ. Microbiol.* **65(4)**, 1610-1618 (1999).
- Thammasuvimol, G., Seo, K.H., Song, K.Y., et al. Optimization of ferrioxamine E concentration as effective supplementation for selective isolation of *Salmonella enteritidis* in egg white. *J. Food Prot.* **69(3)**, 634-638 (2006).

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, 12/07/2022

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