

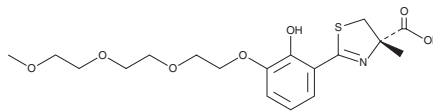
PRODUCT INFORMATION



Deferitazole

Item No. 33329

CAS Registry No.: 945635-15-4
Formal Name: 4S,5-dihydro-2-[2-hydroxy-3-[2-[2-(2-methoxyethoxy)ethoxy]ethoxy]phenyl]-4-methyl-4-thiazolecarboxylic acid
Synonyms: FBS0701, SPD-602
MF: C₁₈H₂₅NO₇S
FW: 399.5
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥5 years



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

Laboratory Procedures

Deferitazole is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as acetonitrile purged with an inert gas can be used. Deferitazole is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Deferitazole is an iron chelator.¹ It decreases the levels of iron in iron-citrate complexes when used at a concentration of 50 μM. Deferitazole (25 μM) also reduces the number of *P. falciparum* oocytes, an effect that can be prevented by the addition of ferric chloride, and the percent of flagellated *P. falciparum* male gametocytes *in vitro*.² It decreases the urine and fecal levels of iron excreted from iron-loaded *C. apella* monkeys when administered at a dose of 150 μmol/kg.³

References

1. Hider, R.C., Kong, X., Abbate, V., *et al.* Deferitazole, a new orally active iron chelator. *Dalton Trans.* **44(11)**, 5197-5204 (2015).
2. Ferrer, P., Vega-Rodriguez, J., Tripathi, A.K., *et al.* Antimalarial iron chelator FBS0701 blocks transmission by *Plasmodium falciparum* gametocyte activation inhibition. *Antimicrob. Agents Chemother.* **59(3)**, 1418-1426 (2015).
3. Bergeron, R.J., Wiegand, J., McManis, J.S., *et al.* Design, synthesis, and testing of non-nephrotoxic desazadesferrithiocin polyether analogues. *J. Med. Chem.* **51(13)**, 3913-3923 (2008).

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, 11/18/2024

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