

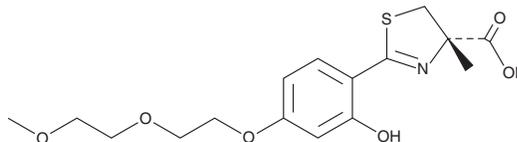
PRODUCT INFORMATION



SP-420

Item No. 28498

CAS Registry No.: 911714-45-9
Formal Name: (4S)-4,5-dihydro-2-[2-hydroxy-4-[2-(2-methoxyethoxy)ethoxy]phenyl]-4-methyl-4-thiazolecarboxylic acid
Synonym: (S)-4'-(HO)-DADFT-norPE
MF: C₁₆H₂₁NO₆S
FW: 355.4
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 273, 309 nm
Supplied as: A crystalline 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

SP-420 is supplied as a crystalline solid. A stock solution may be made by dissolving the SP-420 in the solvent of choice, which should be purged with an inert gas. SP-420 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SP-420 in ethanol and DMSO is approximately 15 mg/ml and approximately 10 mg/ml in DMF.

SP-420 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SP-420 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SP-420 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SP-420 is an iron chelator and a derivative of desferriothiocin.¹ It induces iron excretion in rat bile and urine with an iron-clearing efficiency (ICE) value of 26.7% when administered at a dose of 300 μmol/kg and decreases iron levels in rat liver and heart at 384 μmol/kg per day. SP-420 (75 μmol/kg) also induces iron excretion in iron-overloaded *C. apella* monkeys (ICE = 26.3%).

Reference

- Bergeron, R.J., Bharti, N., Wiegand, J., *et al.* The impact of polyether chain length on the iron clearing efficiency and physiochemical properties of desferrithiocin analogues. *J. Med. Chem.* **53(7)**, 2843-2853 (2010).

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, 06/01/2023

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