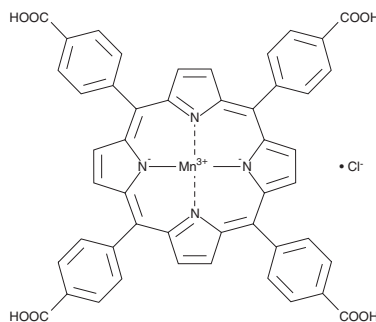


PRODUCT INFORMATION

Mn(III)TBAP (chloride)

Item No. 75850

CAS Registry No.: 55266-18-7
Formal Name: Mn(III)tetrakis(4-benzoic acid) porphyrin, monochloride
MF: $C_{48}H_{28}MnN_4O_8 \cdot Cl$
FW: 879.2
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 467 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Mn(III)TBAP (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the Mn(III)TBAP (chloride) in the solvent of choice, which should be purged with an inert gas. Mn(III)TBAP (chloride) is soluble in organic solvents such as DMSO. The solubility of Mn(III)TBAP (chloride) in this solvent is approximately 0.1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Mn(III)TBAP (chloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. Mn(III)TBAP (chloride) is soluble in neutral or basic aqueous buffers, with or without EDTA. The solubility of Mn(III)TBAP (chloride) in 0.1 M Tris-HCl buffer (1 mM EDTA, pH 9.0) is approximately 10 mg/ml and approximately 0.05 mg/ml in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Mn(III)TBAP is a cell-permeable SOD mimic. At a concentration of 30 μM , it decreases the doubling time of SOD-null *E. coli* by 2-fold.¹

Reference

1. Faulkner, K.M., Liochev, S.I., and Fridovich, I. Stable Mn(III) porphyrins mimic superoxide dismutase *in vitro* and substitute for it *in vivo*. *J. Biol. Chem.* **269**(38), 23471-23476 (1994).

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, 10/05/2022

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