

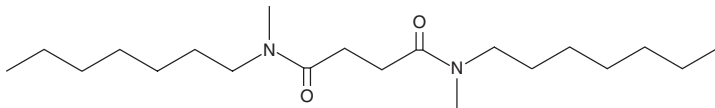
PRODUCT INFORMATION



Magnesium Ionophore I

Item No. 17432

CAS Registry No.: 75513-72-3
Formal Name: N¹,N⁴-diheptyl-N¹,N⁴-dimethyl-butanediamide
Synonyms: ETH 1117
MF: C₂₀H₄₀N₂O₂
FW: 340.5
Purity: ≥98%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Magnesium ionophore I is supplied as a neat oil. A stock solution may be made by dissolving the magnesium ionophore I in the solvent of choice, which should be purged with an inert gas. Magnesium ionophore I is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Magnesium ionophore I is a magnesium sensor that has been used in the development of magnesium-selective microelectrodes to measure magnesium concentrations in biological samples.^{1,2}

References

- Ikeda, K., Morizono, T., Kusakari, J., *et al.* Magnesium ion activity in the mammalian endolymph measured with ion-selective microelectrodes. *Arch. Otorhinolaryngol.* **245**(3), 142-144 (1988).
- Alvarez-Leefmans, F.J., Gamiño, S.M., Giraldez, F., *et al.* Intracellular free magnesium in frog skeletal muscle fibres measured with ion-selective micro-electrodes. *J. Physiol.* **378**, 461-483 (1986).

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, 08/15/2025

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