

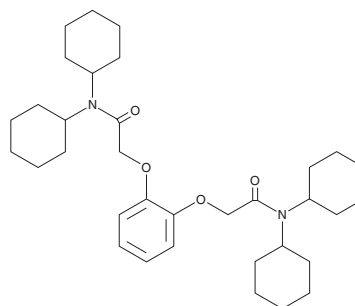
PRODUCT INFORMATION



Sodium Ionophore III

Item No. 22465

CAS Registry No.: 81686-22-8
Formal Name: 2,2'-[1,2-phenylenebis(oxy)]bis[N,N-dicyclohexyl]-acetamide
Synonym: ETH 2120
MF: C₃₄H₅₂N₂O₄
FW: 552.8
Purity: ≥98%
UV/Vis.: λ_{max}: 276 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

Sodium ionophore III is supplied as a crystalline solid. A stock solution may be made by dissolving the sodium ionophore III in the solvent of choice, which should be purged with an inert gas. Sodium ionophore III is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

Sodium ionophore III is a sodium-selective ionophore.^{1,2} It has been used as component of sodium-selective liquid membrane electrodes for the quantification of sodium in aqueous solutions and blood serum.

References

1. Maruizumi, T., Wegmann, D., Suter, G., *et al.* Neutral carrier-based Na⁺-selective electrode for application in blood serum. *Mikrochim. Acta [Wien]* **88(5-6)**, 331-336 (1986).
2. Metzger, E., Dohner, R., and Simon, W. Lithium/sodium ion concentration ratio measurements in blood serum with lithium and sodium ion selective liquid membrane electrodes. *Anal. Chem.* **59(13)**, 1600-1603 (1987).

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, 09/30/2022

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