

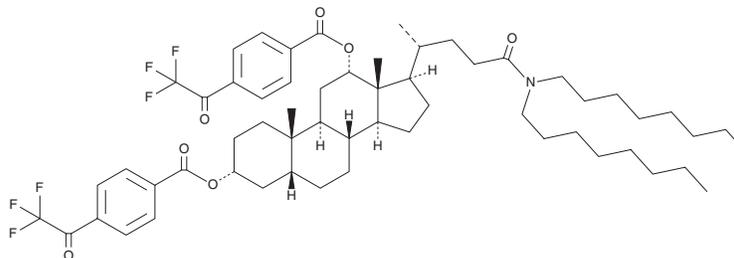
PRODUCT INFORMATION



Carbonate Ionophore VII

Item No. 42078

CAS Registry No.: 222310-82-9
Formal Name: (3 α ,5 β ,12 α)-N,N-dioctyl-3,12-bis[[4-(2,2,2-trifluoroacetyl)benzoyl]oxy]-cholan-24-amide deoxy-3,12-bis(TFAB)CA
Synonym:
MF: C₅₈H₇₉F₆NO₇
FW: 1,016.2
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Carbonate ionophore VII is supplied as a solid. A stock solution may be made by dissolving the carbonate ionophore VII in the solvent of choice, which should be purged with an inert gas. Carbonate ionophore VII is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Carbonate ionophore VII is a carbonate-selective ionophore.¹ It has been used as an ion carrier in carbonate-selective electrodes and a component in the membranes of electric tongues.^{1,2}

References

1. Choi, Y.S., Lvova, L., Shin, J.H., *et al.* Determination of oceanic carbon dioxide using a carbonate-selective electrode. *Anal. Chem.* **74**(10), 2435-2440 (2022).
2. Lenik, J., Wesoły, M., Ciosek, P., *et al.* Evaluation of taste masking effect of diclofenac using sweeteners and cyclodextrin by a potentiometric electronic tongue. *J. Electroanal. Chem.* **780**, 153-159 (2016).

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, 02/11/2025

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