

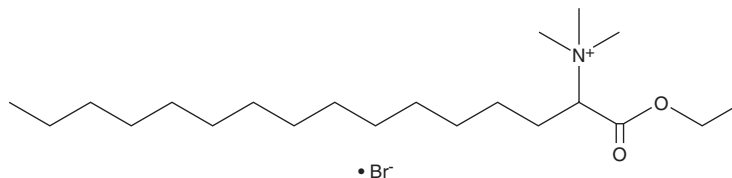
PRODUCT INFORMATION



Carbaethopendecine (bromide)

Item No. 37482

CAS Registry No.: 10567-02-9
Formal Name: 1-ethoxy-N,N,N-trimethyl-1-oxo-2-hexadecanaminium, monobromide
Synonym: Mucoseptonex
MF: C₂₁H₄₄NO₂ • Br
FW: 422.5
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Carbaethopendecine (bromide) is supplied as a crystalline solid. A stock solution may be made by dissolving the carbaethopendecine (bromide) in the solvent of choice, which should be purged with an inert gas. Carbaethopendecine (bromide) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of carbaethopendecine (bromide) in DMSO is approximately 15 mg/ml and approximately 30 mg/ml in ethanol and DMF.

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 carbaethopendecine (bromide) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of carbaethopendecine (bromide) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Carbaethopendecine is a quaternary ammonium-containing cationic surfactant with antibiotic activity.¹ It is active against *S. aureus* and *E. coli* (MIC₅₀s = 0.4 and 2 µg/ml, respectively). Formulations containing carbaethopendecine have been used as antiseptics and disinfectants in veterinary and clinical applications.

Reference

1. Wynne, J.H. and Fulmer, P.A. *Biocides for the Battlefield – Interim Report*. US Naval Research Laboratory (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, 03/28/2023

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