

PRODUCT INFORMATION



Oxtriphylline

Item No. 24006

CAS Registry No.: 4499-40-5

Formal Name: 2-hydroxy-N,N,N-trimethyl-ethanaminium salt with 3,9-dihydro-1,3-dimethyl-1H-purine-2,6-dione

Synonyms: Choline Theophylline, Cholinophylline, Zy 15061

MF: $C_7H_7N_4O_2 \cdot C_5H_{14}NO$

FW: 283.3

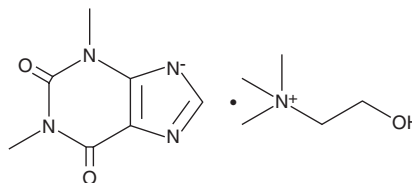
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 272 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

Oxtriphylline is supplied as a crystalline solid. Aqueous solutions of oxtriphylline can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of oxtriphylline in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Oxtriphylline is a choline salt form of theophylline (Item No. 23760). It dose-dependently decreases histamine-induced contractions of isolated guinea pig trachea and increases cAMP levels in guinea pig tracheal smooth muscle cells *in vitro*.¹ Formulations containing oxtriphylline were previously used in the treatment of asthma.

Reference

1. Bilčíková, L., Bauer, V., and Kolena, J. The effects of methylxanthines, ethymizol, ephedrine and papaverine on guinea pig and dog trachea. *Gen. Physiol. Biophys.* **6(2)**, 137-148 (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, 11/16/2022

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