

PRODUCT INFORMATION

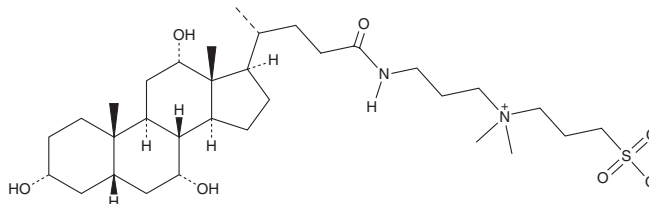


CHAPS

Item No. 44746

CAS Registry No.: 75621-03-3
Formal Name: N,N-dimethyl-N-(3-sulfopropyl)-3-[[[(3 α ,5 β ,7 α ,12 α)-3,7,12-trihydroxy-24-oxocholan-24-yl]amino]-1-propanaminium, inner salt

MF: C₃₂H₅₈N₂O₇S
FW: 614.9
Purity: \geq 98%
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

CHAPS is supplied as a solid. Aqueous solutions of CHAPS can be prepared by directly dissolving the solid in aqueous buffers. CHAPS is soluble (\geq 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

CHAPS is a zwitterionic detergent that has been used in membrane protein solubilization.^{1,2}

References

1. Rodi, P.M., Bocco Gianello, M.D., Corregido, M.C., *et al.* Comparative study of the interaction of CHAPS and Triton X-100 with the erythrocyte membrane. *Biochim Biophys Acta* **1838**(3), 859-866 (2014).
2. Kline, T., Park, H., and Meyerson, L.R. CHAPS solubilization of a G-protein sensitive 5-HT_{1A} receptor from bovine hippocampus. *Life Sci.* **45**(21), 1997-2005 (1989).

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, 12/17/2025

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