

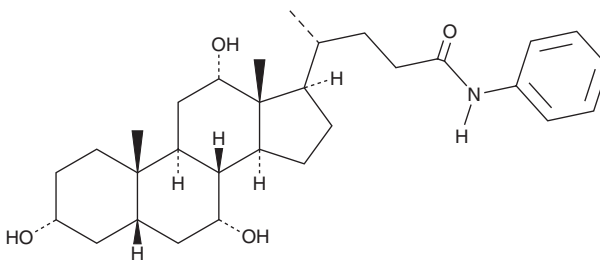
PRODUCT INFORMATION



Cholic Acid anilide

Item No. 26144

CAS Registry No.: 6912-96-5
Formal Name: (5 β)-3 α ,7 α ,12 α -trihydroxy-N-phenyl-cholan-24-amide
Synonym: CaPA
MF: C₃₀H₄₅NO₄
FW: 483.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 243 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

Cholic acid anilide is supplied as a crystalline solid. A stock solution may be made by dissolving the cholic acid anilide in the solvent of choice, which should be purged with an inert gas. Cholic acid anilide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of cholic acid anilide in ethanol is approximately 1 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Cholic acid anilide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cholic acid anilide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cholic acid anilide has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cholic acid anilide is a synthetic bile acid and derivative of cholic acid (Item No. 20250) that inhibits the germination of *C. difficile* strain R20291 spores *in vitro* (IC₅₀ = 1.8 μ M).¹

Reference

1. Sharma, S.K., Yip, C., Esposito, E.X., *et al.* The design, synthesis, and characterizations of spore germination inhibitors effective against an epidemic strain of *Clostridium difficile*. *J. Med. Chem.* **61(15)**, 6759-6778 (2018).

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, 08/24/2023

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