

PRODUCT INFORMATION



BSA-Cholate Bile Acid Complex (5 mM)

Item No. 44505

Synonyms: Bovine Serum Albumin-CA Bile Acid, Bovine Serum Albumin-Cholate Bile Acid, BSA-CA Bile Acid
Supplied as: 5 mM Cholate:0.8 mM BSA (6:1 cholate:BSA) in 150 mM sodium chloride, pH 7.4
Storage: -20°C (as supplied)
Stability: ≥2 years
Item Origin: Animal/Bovine

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

Description

BSA-Cholate bile acid complex (5 mM) is composed of cholic acid (Item No. 20250) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of cholate:BSA. It can be used for efficient bile acid delivery to cells in culture. BSA/BSA-Bile acids are suitable for use in short-term cell culture applications (acute treatment to 18 hours); however, for long-term applications (25+ hours) the product should be filter-sterilized using a 0.2 µm filter and sterile receptacle, which will not affect its performance. For best results, it is recommended that this product be used in conjunction with BSA control for BSA-fatty acid complexes (5 mM) (Item No. 29556).

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, 01/23/2026

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