

PRODUCT INFORMATION



BSA-Laurate Saturated Fatty Acid Complex (5 mM)

Item No. 44990

Synonym:	Bovine Serum Albumin-Laurate, BSA-Dodecanoic Acid
Supplied as:	5 mM Laurate:0.8 mM BSA (6:1 laurate:BSA) in 150 mM sodium chloride, pH 7.4
Storage:	-20°C; avoid freeze/thaw cycles by aliquoting protein
Stability:	≥1 year
Item Origin:	Animal/Bovine

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

Description

BSA-Laurate saturated fatty acid complex is composed of lauric acid (Item No. 10006626) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of laurate:BSA. It has been used for efficient fatty acid delivery to cells in culture for the purpose of monitoring fatty acid oxidation or similar processes in various cellular metabolic studies.^{1,2} BSA/BSA-Fatty 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).

References

1. Alsabeeh, N., Chausse, B., Kakimoto, P.A., *et al.* Cell culture models of fatty acid overload: Problems and solutions. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1863**(2), 143-151 (2018).
2. Tsenkova, M., Brauer, M., Pozdeev, V.I., *et al.* Ketogenic diet suppresses colorectal cancer through the gut microbiome long chain fatty acid stearate. *Nat. Commun.* **16**(1), 1792 (2025).

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, 01/22/2026

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