

PRODUCT INFORMATION

BSA-*trans*-Vaccenate Monounsaturated Fatty Acid Complex (1 mM) Item No. 44252

Overview and Properties

Synonyms:	Bovine Serum Albumin- <i>trans</i> -Vaccenate, BSA- <i>trans</i> -11-Octadecenoic Acid
Supplied in:	1 mM <i>trans</i> -Vaccenate:0.17 mM BSA (6:1 <i>trans</i> -Vaccenate:BSA) in 150 mM sodium chloride, pH 7.4
Storage:	-20°C
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-*trans*-Vaccenate monounsaturated fatty acid complex is composed of *trans*-vaccenic acid (Item No. 15301) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of *trans*-vaccenate: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 (1 mM) (Item No. 34932).

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. Sarnyai, F., Donkó, M.B., Mátyási, J., *et al.* Cellular toxicity of dietary trans fatty acids and its correlation with ceramide and diglyceride accumulation. *Food Chem. Toxicol.* **124**, 324-335 (2019).

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