

PRODUCT INFORMATION

BSA- α -Linolenate ω -3 Polyunsaturated Fatty Acid Complex (5 mM)

Item No. 44864

Overview and Properties

Synonyms:	Bovine Serum Albumin-ALA, Bovine Serum Albumin- α -Linolenate, BSA-ALA
Supplied as:	5 mM α -Linolenate:0.8 mM BSA (6:1 α -Linolenate:BSA) in 150 mM sodium chloride, pH 7.4
Storage:	-20°C (as supplied)
Stability:	\geq 2 years
Item Origin:	Animal/Bovine

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

Description

BSA- α -Linolenate ω -3 polyunsaturated fatty acid complex is composed of α -linolenic acid (Item No. 90210) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of α -linolenate: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.¹⁻³ 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 μ m 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. Zhang, J., Kris-Etherton, P.M., Thompson, J.T., *et al.* Alpha-linolenic acid increases cholesterol efflux in macrophage-derived foam cells by decreasing stearoyl CoA desaturase 1 expression: Evidence for a farnesoid-X-receptor mechanism of action. *J. Nutr. Biochem.* **23**(4), 400-409 (2012).
3. Hauge, M., Vestmar, M.A., Husted, A.S., *et al.* GPR40 (FFAR1) - Combined Gs and Gq signaling *in vitro* is associated with robust incretin secretagogue action *ex vivo* and *in vivo*. *Mol. Metab.* **4**(1), 3-14 (2015).

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