

PRODUCT INFORMATION

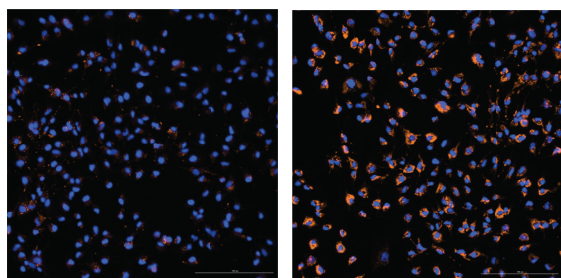
BSA- α -Linolenate ω -3 Polyunsaturated Fatty Acid Complex (1 mM)

Item No. 39063

Supplied as: 1 mM α -linolenate:0.17 mM BSA (6:1 α -linolenate:BSA) in 150 mM sodium chloride, pH 7.4
Synonyms: Bovine Serum Albumin-ALA, Bovine Serum Albumin- α -Linolenate, BSA-ALA, BSA-9(Z),12(Z),15(Z)-Octadecatrienoic Acid
Storage: -20°C
Stability: $\geq$ 2 years
Item Origin: Animal/Bovine

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

Image



BSA Control

BSA- α -Linolenate

Huh7 hepatocytes were plated and treated with 200 μ M BSA- α -Linolenate ω -3 Polyunsaturated Fatty Acid Complex (Item No. 39063; right) or BSA Control (Item No. 34932; left) for 24 hours. On day two, cells were incubated with Hoechst (blue) and Nile Red (red) for 15 minutes before imaging, showing lipid droplet accumulation.

Description

BSA- α -Linolenate ω -3 Polyunsaturated Fatty Acid Complex (1 mM) 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 can be used for efficient fatty acid delivery to cells in culture for the purpose of monitoring lipid metabolism and inflammatory signaling pathways and studying the role of α -linolenic acid in neuroprotection.¹⁻⁴ Cayman's BSA- α -Linolenate ω -3 Polyunsaturated Fatty Acid Complex (1 mM) is 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 Cayman's BSA Control for BSA-Fatty Acid Complexes (1 mM) (Item No. 34932), prepared with fatty acid-free BSA.

References

1. Alsabeeh, N., Chausse, B., Kakimoto, P.A., et al. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1863**(2), 143-151 (2018).
2. Zhang, J., Kris-Etherton, P.M., Thompson, J.T., et al. *J. Nutr. Biochem.* **23**(4), 400-409 (2012).
3. Hauge, M., Vestmar, M.A., Husted, A.S., et al. *Mol. Metab.* **4**(1), 3-14 (2015).
4. Blondeau, N., Lipsky, R.H., Bourourou, M., et al. *Biomed Res. Int.* 519830 (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, 07/13/2026

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