

PRODUCT INFORMATION

BSA-Docosahexaenoate-d₅ Polyunsaturated Fatty Acid Complex (1 mM) Item No. 43591

Synonyms:	Bovine Serum Albumin-DHA-d ₅ , Bovine Serum Albumin-Docosahexaenoate-d ₅ , BSA-Cervonate-d ₅ , BSA-DHA -d ₅
Supplied as:	1 mM Docosahexanoate d ₅ :0.17 mM BSA (6:1 docosahexanoate d ₅ :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-Docosahexaenoate-d₅ Polyunsaturated Fatty Acid Complex (1 mM) is composed of docosahexaenoic acid-d₅ (DHA-d₅; Item No. 10005057) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of docosahexaenoate-d₅:BSA. It is intended for the quantification of DHA-d₅ uptake and incorporation by GC- or LC-MS. BSA-Docosahexaenoate-d₅ Polyunsaturated Fatty Acid Complex (1 mM) can be used for efficient fatty acid delivery to cells in culture for the purpose of monitoring lipid metabolism, inflammatory signaling pathways, fatty acid uptake and efflux, and ferroptotic cell death.¹⁻⁵ Cayman's BSA-Docosahexaenoate-d₅ 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 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 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.* Cell culture models of fatty acid overload: Problems and solutions. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1863**(2), 143-151 (2018).
2. Katsnelson, G. and Ceddia, R.B. Docosahexaenoic and eicosapentaenoic fatty acids differentially regulate glucose and fatty acid metabolism in L6 rat skeletal muscle cells. *Am. J. Physiol. Cell Physiol.* **319**(6), C1120-C1129 (2020).
3. Low, Y.L., Pan, Y., Short, J.L., *et al.* Profiling the expression of fatty acid-binding proteins and fatty acid transporters in mouse microglia and assessing their role in docosahexaenoic acid-d₅ uptake. *Prostaglandins Leukot. Essent. Fatty Acids* **171**:102303, (2021).
4. Ochiai, Y., Uchida, Y., Tachikawa, M., *et al.* Amyloid beta₂₅₋₃₅ impairs docosahexaenoic acid efflux by down-regulating fatty acid transport protein 1 (FATP1/SLC27A1) protein expression in human brain capillary endothelial cells. *J. Neurochem.* **150**(4), 385-401 (2019).
5. von Krusenstiern, A.N., Robson, R.N., Qian, N., *et al.* Identification of essential sites of lipid peroxidation in ferroptosis. *Nat. Chem. Biol.* **19**(6), 719-730 (2023).

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, 06/18/2026

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